

Thermo Scientific

TRACE GC Ultra

Gas Chromatograph

Standard Operating Procedures

PN 317 092 00, Revision May 2010_B



TRACE™ GC Ultra - Standard Operating Procedures

May 2010 Edition

Part Number 31709200

© 2007-2010 Thermo Fisher Scientific Inc. All rights reserved.

Printed in Italy

Published by Thermo Fisher Scientific S.p.A., Strada Rivoltana, 20090 Rodano - Milan - Italy

Tel: +39 02 95059355 Fax: +39 02 95059388

Printing History: First Edition, released June 1998.

Second Edition, released November 1998.

Third Edition, released June 1999.

Fourth Edition, released May 2001

Fifth Edition, released January 2002

Sixth Edition, released, May 2003

Seventh Edition, released April 2004

Eighth Edition, released January 2005

Ninth Edition, released June 2005

Tenth Edition, released December 2005

Eleventh Edition, released June 2006

Twelfth Edition, released January 2007

Thirteenth Edition, released May 2007

Fourteenth Edition, released February 2008

Fifteenth Edition, released April 2009

Sixteenth Edition, released September 2009

Seventeenth Edition, released May 2010_B

Disclaimer

Technical Information contained in this publication is for reference purposes only and is subject to change without notice. Every effort has been made to supply complete and accurate information; however, Thermo Fisher Scientific assumes no responsibility and will not be liable for any errors, omissions, damage, or loss that might result from any use of this manual or the information contained therein (even if this information is properly followed and problems still arise).

This publication is not part of the Agreement of Sale between Thermo Fisher Scientific and the purchaser of a TRACE™ GC Ultra system. In the event of any conflict between the provisions of this document and those contained in Thermo Fisher Scientific's Terms and Conditions, the provisions of the Terms and Conditions shall govern.

Reference to System Configurations and Specifications supercede all previous information and are subject to change without notice.

Trademarks

TRACE™ GC Ultra is a trademark of Thermo Fisher Scientific Inc., and its subsidiaries. Other brand and product names may be trademarks or registered trademarks of their respective companies.

Contents

About This Manual	xiii
Overview	xiii
Conventions Used in This Manual.....	xiv
Abbreviations for Injectors and Detectors	xvi
Instrument Markings and Symbols	xvii
Safety Information	xix
Precaution for Gases	xix
Precaution for the Electron Capture Detector.....	xix
Using the TRACE GC Ultra Document Set.....	xx

SECTION I

SOPs Overview

Chapter 1

General Overview	25
Scope	26
Parts Referenced.....	27
Getting Started	28

SECTION II

SOPs Using FID

Chapter 2

Checkout Using FID with S/SL Injector	37
SOP Number: P0292/07/E - 01 September 2009	38
Scope.....	38
Parts Referenced	38
Analytical Conditions Required for Splitless Injection	39
Recommended Initial Operations.....	40

Chapter 3

Checkout Using FID with OC Injector	47
SOP Number: P0293/08/E - 01 September 2009	48
Scope.....	48
Parts Referenced	48
Analytical Conditions Required for On-Column Injection.....	49
Recommended Initial Operations.....	50

Chapter 4

Checkout Using FID with PKD Injector57
SOP Number: P0307/07/E - 01 September 2009 58
Scope 58
Parts Referenced 58
Analytical Conditions Required for Packed Injector 60
Recommended Initial Operations 61

Chapter 5

Checkout Using FID with PPKD Injector69
SOP Number: P0308/07/E - 01 September 2009 70
Scope 70
Parts Referenced 70
Analytical Conditions Required for Purged Packed Injector 72
Recommended Initial Operations 73

Chapter 6

Checkout Using FID with PTV Injector81
SOP Number: P0309/07/E - 01 September 2009 82
Scope 82
Parts Referenced 82
Analytical Conditions Required for PTV Splitless Injection 83
Recommended Initial Operations 84

SECTION III

SOPs Using ECD

Chapter 7

Checkout Using ECD with S/SL Injector95
SOP Number: P0294/07/E - 01 September 2009 96
Scope 96
Parts Referenced 96
Analytical Conditions Required for Splitless Injection 97
Recommended Initial Operations 98

Chapter 8

Checkout Using ECD with OC Injector107
SOP Number: P0295/08/E - 01 September 2009 108

Scope.....	108
Parts Referenced	108
Analytical Conditions Required for On-Column Injection.....	109
Recommended Initial Operations.....	110
Chapter 9	
Checkout Using ECD with PKD Injector.....	119
SOP Number: P0310/07/E - 01 September 2009	120
Scope.....	120
Parts Referenced	120
Analytical Conditions Required for Packed Injector.....	122
Recommended Initial Operations.....	123
Chapter 10	
Checkout Using ECD with PPKD Injector.....	131
SOP Number: P0311/07/E - 01 September 2009	132
Scope.....	132
Parts Referenced	132
Analytical Conditions Required for Purged Packed Injector.....	134
Recommended Initial Operations.....	135
Chapter 11	
Checkout Using ECD with PTV Injector.....	143
SOP Number: P0312/07/E - 01 September 2009	144
Scope.....	144
Parts Referenced	144
Analytical Conditions Required for PTV Splitless Injection.....	145
Recommended Initial Operations.....	146
SECTION IV	
SOPs Using NPD	
Chapter 12	
Checkout Using NPD with S/SL Injector.....	157
SOP Number: P0296/08/E - 01 September 2009	158
Scope.....	158
Parts Referenced	158
Analytical Conditions Required for Splitless Injection	159

Contents

Recommended Initial Operations	160
Chapter 13	
Checkout Using NPD with OC Injector	169
SOP Number: P0297/08/E - 01 September 2009	170
Scope	170
Parts Referenced	170
Analytical Conditions Required for On-Column Injection	171
Recommended Initial Operations	172
Chapter 14	
Checkout Using NPD with PKD Injector.....	181
SOP Number: P0313/08/E - 01 September 2009	182
Scope	182
Parts Referenced	182
Analytical Conditions Required for Packed Injector.....	184
Recommended Initial Operations	185
Chapter 15	
Checkout Using NPD with PPKD Injector	195
SOP Number: P0314/08/E - 01 September	196
Scope	196
Parts Referenced	196
Analytical Conditions Required for Purged Packed Injector	198
Recommended Initial Operations	199
Chapter 16	
Checkout Using NPD with PTV Injector	209
SOP Number: P0315/08/E - 01 September 2009	210
Scope	210
Parts Referenced	210
Analytical Conditions Required for PTV Splitless Injection	211
Recommended Initial Operations	212

SECTION V SOPs Using FPD

Chapter 17	
Checkout Using FPD with S/SL Injector	225
SOP Number: P0316/08/E - 01 September 2009	226
Scope	226
Parts Referenced	226
Analytical Conditions Required for Splitless Injection	227
Recommended Initial Operations.....	228
Chapter 18	
Checkout Using FPD with OC Injector.....	237
SOP Number: P0317/09/E - 01 September 2009	238
Scope.....	238
Parts Referenced	238
Analytical Conditions Required for On-Column Injection.....	239
Recommended Initial Operations.....	240
Chapter 19	
Checkout Using FPD with PKD Injector	249
SOP Number: P0318/08/E - 01 September 2009	250
Scope.....	250
Parts Referenced	250
Analytical Conditions Required for Packed Injector	252
Recommended Initial Operations.....	253
Chapter 20	
Checkout Using FPD with PPKD Injector	263
SOP Number: P0319/08/E - 01 September 2009	264
Scope.....	264
Parts Referenced	264
Analytical Conditions Required for Purged Packed Injector.....	266
Recommended Initial Operations.....	267
Chapter 21	
Checkout Using FPD with PTV Injector.....	277
SOP Number: P0320/08/E - 01 September 2009	278
Scope.....	278
Parts Referenced	278
Analytical Conditions Required for PTV Splitless Injection.....	279
Recommended Initial Operations.....	280

**SECTION VI
SOPs Using PID**

Chapter 22

Checkout Using PID with S/SL Injector293
 SOP Number: P0321/09/E - 01 September 2009 294
 Scope 294
 Parts Referenced 294
 Analytical Conditions Required for Splitless Injection 296
 Recommended Initial Operations 297

Chapter 23

Checkout Using PID with OC Injector307
 SOP Number: P0322/09/E - 01 September 2009 308
 Scope 308
 Parts Referenced 308
 Analytical Conditions Required for On-Column Injection 309
 Recommended Initial Operations 310

Chapter 24

Checkout Using PID with PTV Injector321
 SOP Number: P0325/08/E - 01 September 2009 322
 Scope 322
 Parts Referenced 322
 Analytical Conditions Required for PTV Splitless Injection 324
 Recommended Initial Operations 325

**SECTION VII
SOPs Using TCD**

Chapter 25

Checkout Using TCD with S/SL Injector339
 SOP Number: P0326/11/E - 01 September 2009 340
 Scope 340
 Parts Referenced 340
 Analytical Conditions Required for Splitless Injection 341
 Recommended Initial Operations 342

Chapter 26

Checkout Using TCD with PKD Injector	349
SOP Number: P0328/11/E - 01 September 2009	350
Scope	350
Parts Referenced	350
Analytical Conditions Required for Packed Injector	352
Recommended Initial Operations.....	353

Chapter 27

Checkout Using TCD with PPKD Injector	361
SOP Number: P0329/11/E - 01 September 2009	362
Scope	362
Parts Referenced	362
Analytical Conditions Required for Purged Packed Injector.....	364
Recommended Initial Operations.....	365

SECTION VIII**SOPs Using PDD****Chapter 28**

Checkout Using PDD with S/SL Injector.....	375
SOP Number: P0381/05/E - 01 September 2009	376
Scope	376
Parts Referenced	376
Analytical Conditions Required for Splitless Injection	377
Recommended Initial Operations.....	378

Chapter 29

Checkout Using PDD with OC Injector	385
SOP Number: P0382/06/E - 01 September 2009	386
Scope	386
Parts Referenced	386
Analytical Conditions Required for On-Column Injection.....	387
Recommended Initial Operations.....	388

SECTION IX**SOPs Using FID-NPD-FPD in Stacked Configuration**

Chapter 30

Checkout Using Tandem FID.....397
SOP Number: P0383/05/E - 01 September 2009 398
Scope 398
Checkout Overview 399
Important Considerations 400
 FID Gas Required..... 400
 Column Installation 400
 FID Detector and Signal Menus 400

Chapter 31

Checkout Using Tandem NPD.....405
SOP Number: P0384/06/E - 01 September 2009 406
Scope 406
Checkout Overview 407
Important Considerations 408
 NPD Gas Required 408
 Column Installation 408
 NPD Detector and Signal Menus 408

Chapter 32

Checkout Using Tandem FPD413
SOP Number: P0385/06/E - 01 September 2009 414
Scope 414
Checkout Overview 415
Important Considerations 416
 FPD Gas Required..... 416
 Column Installation 416
 FPD Detector and Signal Menus 416

SECTION X

SOPs for Large Volume Applications

Chapter 33

Checkout Using FID with LVSL Injector423
SOP Number: P0430/04/E - 01 September 2009 424
Scope 424
Parts Referenced..... 424

Analytical Conditions Required for LVSL Injection.....	426
Recommended Initial Operations.....	427
Chapter 34	
Checkout Using FID with LVOC Injector.....	435
SOP Number: TE P0609/02/E-01 September 2009	436
Scope.....	436
Parts Referenced	436
Analytical Conditions Required for LVOC Injection	438
Analytical Condition for the GC.....	438
Recommended Initial Operations.....	440
Appendix A	
Customer Communication	445
How To Contact Us.....	445
Reader Survey	446
Glossary.....	447
Index	453

Contents

About This Manual

Overview

This manual is organized as follows:

Section **I**, *SOPs Overview*, contains a general description of the Standard Operating Procedures.

Section **II**, *SOPs Using FID*, contains the procedure to test the TRACE GC Ultra with the Flame Ionization Detector (FID) using different injectors.

Section **III**, *SOPs Using ECD*, contains the procedure to test the TRACE GC Ultra with the Electron Capture Detector (ECD) using different injectors.

Section **IV**, *SOPs Using NPD*, contains the procedure to test the TRACE GC Ultra with the Nitrogen Phosphorus Detector (NPD) using different injectors.

Section **V**, *SOPs Using FPD*, contains the procedure to test the TRACE GC Ultra with the Flame Photometric Detector (FPD) using different injectors.

Section **VI**, *SOPs Using PID*, contains the procedure to test the TRACE GC Ultra with the Photoionization Detector (PID) using different injectors.

Section **VII**, *SOPs Using TCD*, contains the procedure to test the TRACE GC Ultra with the Thermal Conductivity Detector (TCD) using different injectors.

Section **VIII**, *SOPs Using PDD*, contains the procedure to test the TRACE GC Ultra with the Pulsed Discharge Detector (PDD) using different injectors.

Section **IX**, *SOPs Using FID-NPD-FPD in Stacked Configuration*, contains the procedures to test the TRACE GC Ultra with the Flame Ionization Detector (FID), Nitrogen-Phosphorus Detector (NPD) or Flame Photometric Detector (FPD) in series (stacked configuration) with the Electron Capture Detector ECD using different injectors.

Section **X**, *SOPs for Large Volume Applications*, contains the procedures to test the TRACE GC Ultra for large volume application by using different injectors.



Appendix A, *Customer Communication*, contains contact information for Thermo Fisher Scientific offices worldwide. Use the *Reader Survey* in this section to give us feedback on this manual and help us improve the quality of our documentation.

The *Glossary* contains definitions of terms used in this guide and the help diskette. This also includes abbreviations, acronyms, metric prefixes, and symbols.

The *Index* contains an alphabetical list of key terms and topics in this guide, including cross references and the corresponding page numbers.

Conventions Used in This Manual

The following symbols and typographical conventions are used throughout this manual.

Bold	Bold text indicates names of windows, menus, dialog boxes, buttons, and fields.
<i>Italic</i>	Italic indicates cross references, first references to important terms defined in the glossary, and special emphasis.
Monospace	Monospace, or Courier, indicates filenames and filepaths, or to indicate text the user should enter with the keyboard.
Monospace Bold	Monospace Bold indicates messages or prompts displayed on the computer screen or on a digital display.
»	This symbol illustrates menu paths to select, such as File»Open...
KEY NAME	Bold, uppercase sans serif font indicates the name of a key on a keyboard or keypad, such as ENTER .
 CAUTION	This symbol alerts you to an action or procedure that, if performed improperly, could damage the instrument.
 NOTE	This symbol alerts you to important information related to the text in the previous paragraph.



WARNING!

This symbol alerts you to an action or procedure that, if performed improperly, could result in damage to the instrument or possible physical harm to the user. This symbol may be followed by icons indicating special precautions that should be taken to avoid injury.



This symbol indicates danger from high temperature surfaces or substances.



This symbol indicates an explosion hazard.



This symbol indicates the presence of radioactive material.



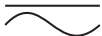







Abbreviations for Injectors and Detectors








	Abbreviation
Split/splitless Injector	S/SL
Large Volume Splitless	LVSL
Cold On-Column Injector	OCI
Packed Column Injector	PKD
Packed Column with Septum Purge	PPKD
Programmable Temperature Vaporizing Injector	PTV

Detector	Abbreviation
Flame Ionization Detector	FID
Electron Capture Detector	ECD
Nitrogen-Phosphorus Detector	NPD
Flame Photometric Detector	FPD
Photoionization Detector	PID
Thermal Conductivity Detector	TCD
Pulsed Discharge Detector	PDD

Instrument Markings and Symbols

The following table explains the symbols used on Thermo Fisher Scientific instruments. Only a few of them are used on the TRACE GC Ultra gas chromatograph.

Symbol	Description
	Direct Current
	Alternating Current
	Both direct and alternating current
	Three-phase alternating current
	Earth (ground) terminal
	Protective conductor terminal
	Frame or chassis terminal
	Equipotentiality
	On (Supply)
	Off (Supply)

Symbol	Description
	Equipment protected throughout by DOUBLE INSULATION or REINFORCED INSULATION (Equivalent to Class II of IEC 536)
	Caution, risk of electric shock
	Caution, hot surface
	Caution (refer to accompanying documents)
	In-position of a bistable push control
	Out-position of a bistable push control
	Symbol in compliance to the Directive 2002/96/EC on Waste Electrical and Electronic Equipment (WEEE) placed on the european market after August, 13, 2005.

Safety Information

Precaution for Gases

Before using gases, carefully read the indications of hazard and the warning reported in the Safety sheet supplied by the manufacturer with reference to the relevant CAS number (Chemical Abstract Service).



WARNING! Hydrogen is a harmful gas that mixed with air may give rise to an explosion hazard. The use of hydrogen requires the operator's extreme caution and the recourse to special precautions due to the risk involved.



Precaution for the Electron Capture Detector



WARNING! The Electron Capture Detector (ECD) contains a ^{63}Ni beta-emitting radioactive source of 370 MBq (10 mCi). For no reason should the detector be opened or handled by the operator. Any maintenance or service operations involving even partial disassembling of the instrument must be performed **ONLY** by qualified personnel at the laboratory expressly authorized by Thermo Fisher Scientific and specially licensed to handle radioactive material.



Using the TRACE GC Ultra Document Set

The TRACE GC Ultra Document Set (CD-Rom PN 317 095 00) includes all manuals in electronic format, and serves as your library for information about the TRACE GC Ultra hardware and software.

The TRACE GC Ultra Document Set (PN 317 093 00) as paper copy is also available. Furthermore, Thermo Fisher Scientific part numbers (PN) for the paper copy manuals are provided for each book title.

Site Preparation and Installation Manual (PN 317 091 90)

This manual and diskette describes how to set up a workspace for the TRACE GC Ultra and how to connect the TRACE GC Ultra to the gas supplies and peripheral devices.

Acceptance Package (PN 317 092 20)

This folder contains required shipping documents and quality report forms.

Getting Started (PN 317 092 30)

This guide contains sequences for checking configuration, installing detectors, and making a first analysis with the TRACE GC Ultra.

Operating Manual (PN 317 091 70)

This manual provides descriptions of the TRACE GC Ultra hardware and software and instructions for their use.

UFM Ultra Fast Module Device (PN 317 093 98)

This manual provides descriptions of the TRACE GC Ultra equipped with the UFM device, and instructions for its use. The relevant *Standard Operating Procedure* is provided in a separated document PN 317 094 09.

Quick Reference Card (PN 317 092 40)

This reference card contains guidelines for carrier gas use and injection sequences.

K-Factor Quick Reference (P/N 317 092 41)

This card indicates the theoretical K-Factors related to the carrier gas and the column in use.

Preventive Maintenance Schedule (PN 317 092 80)

This document provides a list of recommended scheduled maintenance and a year-long log book to record maintenance, observations, supply lists, and service records.

Maintenance and Troubleshooting Manual (PN 317 091 80)

This manual contains instructions for diagnosing and resolving operational problems.

Standard Operating Procedures (PN 317 092 00)

This manual contains instructions, operating sequences, and test criteria for final testing of the TRACE GC Ultra.

Spare Parts Catalog (PN 317 092 10)

This catalog contains a list of spare parts for the TRACE GC Ultra.

SECTION

I

SOPs Overview

The *SOPs Overview* section contains a general description of the Standard Operating Procedures.

Chapter 1 *General Overview*, contains a guideline to apply correctly the SOPs.



General Overview

Chapter at a Glance...

Scope	26
Parts Referenced	27
Getting Started	28

Operating Procedures

Test Column Conditioning	31
--------------------------------	----

Scope

The Standard Operating Procedures (SOP) described in this book are a series of instructions, operations and test criteria derived from our quality policy procedures used for final testing of the TRACE GC Ultra. The SOPs have been developed to test and verify instrument complete analytical performance after the installation has been completed. This will help you as a guideline, to check if your TRACE GC Ultra continues to perform according to the original checkout testing specifications carried out in the factory premises. However, these tests alone cannot define if the instrument is not performing according to the original specifications. The checkout is carried out injecting a standard solution into a test column under analytical conditions set according to the injector(s) and detector(s) hardware provided with the GC. Before starting the test checkout, refer to the Parts Referenced and the Analytical Condition required.



NOTE

Each SOP has a proper Registration and Revision Number (e.g. P0292/01/E - 12 June 1998), according to our Quality Management policy.

If your GC is equipped with the Ultra Fast Module, please refer to the relevant SOP (PN 317 094 09).

For specific operating or maintenance questions, please refer to the following manuals:

- TRACE GC Ultra Operating Manual PN 317 091 70
- TRACE GC Ultra Maintenance and Troubleshooting Manual PN 317 091 80
- TRACE GC Ultra Site Preparation and Installation Manual PN 317 091 90
- TRACE GC Ultra Getting Started Manual PN 317 092 30

Parts Referenced

The SOPs require the following parts:

Table 1-1. SOPs Parts Referenced

	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long; 0.32 mm ID; 0.25 µm film thickness.	260 800 01
	Graphite ferrule for 0.32 mm ID column	290 134 87
Syringes	10 µl size; 70 mm needle length for S/SL injections	365 001 03
	10 µl size; 75 mm needle length for OC injections	365 020 07
	10 µl size; 50 mm needle length for PTV, PKD and PPKD injections	365 005 25
Test Mixtures	Test Mixture for FID checkout	338 190 20
	Test Mixture for ECD checkout	338 190 11
	Test Mixture for NPD checkout	338 190 06
	Test Mixture for FPD checkout	338 190 06
	Test Mixture for PID checkout	338 190 06
	Test Mixture for TCD checkout	338 190 16
	Test Mixture for PDD checkout	338 190 32
Gases	Gas Chromatographic-grade purity <i>Carrier Gas</i> = Helium <i>Fuel Gases</i> = Hydrogen - Air <i>Make-up Gas</i> = Nitrogen <i>Discharge gas for PDD</i> = Helium ultrapure (At least 99.999% of purity)	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Getting Started

Before starting checkout, perform the following preliminary operations sequentially:

1. **Gas Supply Connections**

Connect the gas supplies following the instructions reported in Chapter 4 of the TRACE GC Ultra Site Preparation and Installation Manual

2. **Data Handling Connections**

Connect your data handling following the instructions reported in Chapter 5 of the TRACE GC Ultra Site Preparation and Installation Manual

3. **Column Installation**

Install the test column according to the injector installed in your GC following the **Operating Sequences** reported in Chapter 14 “*Columns*” of the TRACE GC Ultra Operating Manual.

- *Preparing a Capillary Column*
- *Connecting a Capillary Column to a Split/Splitless Injector*
- *Connecting a Capillary Column to an Cold On-Column Injector*
- *Connecting a Capillary Column to an Packed Column Injector*
- *Connecting a Capillary Column to an Purged Packed Injector*
- *Connect a Capillary Column to an Programmable Temperature Vaporizing Injector*

4. **Glass Liner and Septum Installation**

Install the glass liner following the **Operating Sequences** “*Install a Liner and Septum*” reported in the TRACE GC Ultra Operating Manual.

- Chapter 5: *Split/Splitless Injector*
- Chapter 9: *Packed Column Injector*

- Chapter 10: *Purged Packed Injector*
- Chapter 11: *Programmable Temperature Vaporizing Injector*

5. Column Leak Test

Perform the column leak check following the **Operating Sequence** “*Performing a Leak Check*” reported in Chapter 14 “*Columns*” of the TRACE GC Ultra Operating Manual.

6. Column Evaluation

Set column length, nominal ID and film thickness to calculate the column K factor. It is also possible to manually set the carrier gas flow measured at the end of the column to obtain the effective K factor.

Perform column evaluation following the **Operating Sequence** “*Performing a Column Evaluation*” reported in Chapter 14 “*Columns*” of the TRACE GC Ultra Operating Manual.

7. Column Conditioning



CAUTION

When performing column conditioning, the column should be connected only to the injector leaving the column outlet disconnected to avoid the possibility of contamination of the detector base body.

Column conditioning consists of passing a flow of carrier gas through the column and heating it to a temperature of 20-50 °C above the maximum temperature that will be used for running the analysis. If the Cold On-Column Injector is used, ensure that the injection valve is closed. Refer to Table 1-2 for the parameter setting and to the **Operating Procedure** “*Test Column Conditioning*” on page 31, to perform the operation.

8. Detector Connections

This operation should be carried out at the end of the column conditioning. Connect the test column to the detector following the **Operating Sequences** reported in Chapter 14 “*Columns*” of the TRACE GC Ultra Operating Manual.

- *Connecting a Capillary Column to an FID, FPD or NPD Detector*
- *Connecting a Capillary Column to an ECD Detector*

- *Connect a Capillary Column to a PID Detector*
- *Connecting a Capillary Column to a TCD Detector*
- *Connecting a Capillary Column to a PDD Detector*

OPERATING PROCEDURE

Test Column Conditioning

Refer to Table 1-2 for the parameters setting:

Table 1-2. Column Conditioning Parameters

Gases	Carrier Gas: Helium = 30 kPa Constant Pressure
Oven Program	Initial Temperature = 50 °C Initial Time = 1 min. Ramp 1 = 20 °C/min. Final Temperature = 250 °C Final Time = 30 min.
Injector	Temperature = according to the injector in use

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	S/SL, OCI, PKD, PPKD or PTV
Left carrier or Right carrier	He (helium)

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to `Pressure mode`, press **MODE/TYPE** to access the selection menu, then select `Constant pressure`. Scroll to `Pressure` and set the pressure value to have the required carrier gas flow rate `Col.flow`.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col.flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

- Using **OVEN**, enter the Column Oven Control Table and set the oven temperature and the Oven Program required.

OVEN		
Temp	50.0	50.0
Initial Time		1.00
Ramp 1		20.0
Final temp		250
Final time 1		30.0<
Ramp 2		Off

- Using **LEFT INLET** or **RIGHT INLET**, enter the appropriate Injector Control Table and set the required temperature setpoint Temp. In the case of Cold On-Column Injector, this step is not required.

LEFT INLET (S/SL) ¹		
Temp	230	230<
Pressure	30.0	30.0
Mode:		Splitless

1. These settings could also be for a right inlet.

or

LEFT INLET (OCI) ¹		
Pressure	30.0	30.0<
Sec. Cool time		Off

1. These settings could also be for a right inlet.

or

LEFT INLET (PKD) ¹		
Temp	200	200
Pressure	30.0	30.0

1. These settings could also be for a right inlet.

or

LEFT INLET (PPKD) ¹		
Temp	200	200
Pressure	30.0	30.0
Mode:	Wide bore<	
Constant sept purge?		Y<

1. These settings could also be for a right inlet.

or

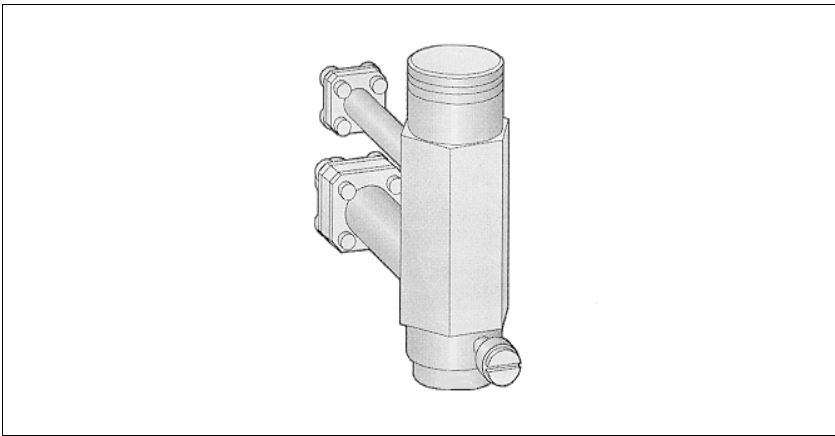
RIGHT INLET (PTV)		
Temp	70	70
Pressure	30.0	30.0
Mode:	PTV Splitless	

Press **PREP RUN** then **START** on the GC to begin the column conditioning.

SECTION

II

SOPs Using FID



The *SOPs Using FID* section, contains the procedures to test the TRACE GC Ultra with the Flame Ionization Detector (FID) using different injectors.

Chapter 2, *Checkout Using FID with S/SL Injector*.

Chapter 3, *Checkout Using FID with OC Injector*.

Chapter 4, *Checkout Using FID with PKD Injector*.

Chapter 5, *Checkout Using FID with PPKD Injector*.

Chapter 6, *Checkout Using FID with PTV Injector*.

Checkout Using FID with S/SL Injector

Chapter at a Glance...

SOP Number: P0292/07/E - 01 September 2009	38
Scope	38
Parts Referenced	38
Analytical Conditions Required for Splitless Injection	39
Recommended Initial Operations	40

Operating Procedures

FID-S/SL Checkout in Splitless Mode	41
---	----

SOP Number: P0292/07/E - 01 September 2009

Scope

Use the following procedure to verify proper FID operation with the Split/Splitless Injector.

Parts Referenced

Table 2-1. FID-S/SL Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 mt long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	3 mm ID for splitless injection	453 200 32
Liner Seal	Graphite seal for glass liner	290 334 06
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
Septum	Standard septum for S/SL injector	313 032 11
Syringe	10 µl size; 70 mm needle length	365 001 03
Test Mixture	Three components in n-Hexane: <i>Component</i> <i>Concentration</i> Dodecane 20 µg/ml Tetradecane 20 µg/ml Hexadecane 20 µg/ml	338 190 20
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for Splitless Injection

Table 2-2. FID-S/SL Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 35 ml/min Air = 350 ml/min Make-up Gas: Nitrogen = 30 ml/min
Oven Program	Initial Temperature = 50 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 200 °C Final Time = 1 minute
Injector	Operating Mode = Splitless Temperature = 230 °C Splitless Time = 0.8 minutes Split Flow = 60 ml/min Constant Septum Purge = Yes
Detector	Base Temperature = 250 °C Detector Signal Range = 10 ⁰
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Replace the glass liner.
The glass liner currently installed in your injector should be carefully removed and replaced with the 3 mm ID glass liner for splitless application, as required for the checkout, with the appropriate liner seal.
2. Replace the septum
A new septum should be installed properly in your injector.
3. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
4. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
5. Perform Column Evaluation and Leak Test.
6. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

FID-S/SL Checkout in Splitless Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	S/SL
Left carrier or Right carrier	He (helium)
Left detector or Right detector	FID

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to Pressure and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	50.0	50.0
Initial Time		1.00
Ramp 1		20.0
Final temp		200
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate Split/Splitless Injector Control Table. Set the required temperature *Temp* setpoint. Verify to

operate in **Splitless** mode. If not, scroll to **Mode**, press **MODE/TYPE** to access the selection menu, then select **Splitless**. Scroll to **Splitless time** to set the required setpoint.

LEFT INLET (S/SL) ¹		
Temp	230	230
Pressure	30.0	30.0
Mode:	Splitless	
Total flow	(63.0)	
Split Flow	60.0	60.0
Splitless time	0.80	0.80
Constant sept purge?		Y<

1. These settings could also be for a right inlet.

4. Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate FID Detector Control Table. Set the required temperature **Base Temp** and the detector gases **H2**, **Air** and **Mkup** required setpoints.

LEFT DETECTOR (FID) ¹		
Flame		Off
Base temp	250	250
Signal pA		(5.5)
Ign.thresh		2.0
Flameout retry		Off
H2	35	35
Air	350	350
Mkup N2	30	30<

1. These settings could also be for a right detector.

5. Ignite the FID flame scrolling to **Flame** and pressing **ON**.
6. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate FID Detector Signal Control Table. Observe the FID flame signal at the display. This is the flame-on background offset. Scroll to **Range** and set the electrometer amplifier input range required.

LEFT SIGNAL (FID) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline Comp	Off

1. These settings could also be for a right signal.

7. Activate your Data System and set the parameters required for the checkout.
8. In the FID Detector Signal Control Table, scroll to `Auto zero?` and turn it **YES**.
9. Perform a blank analysis injecting pure hexane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 2-3 according to the data handling in use.

10. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
11. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
12. Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 2.1*.

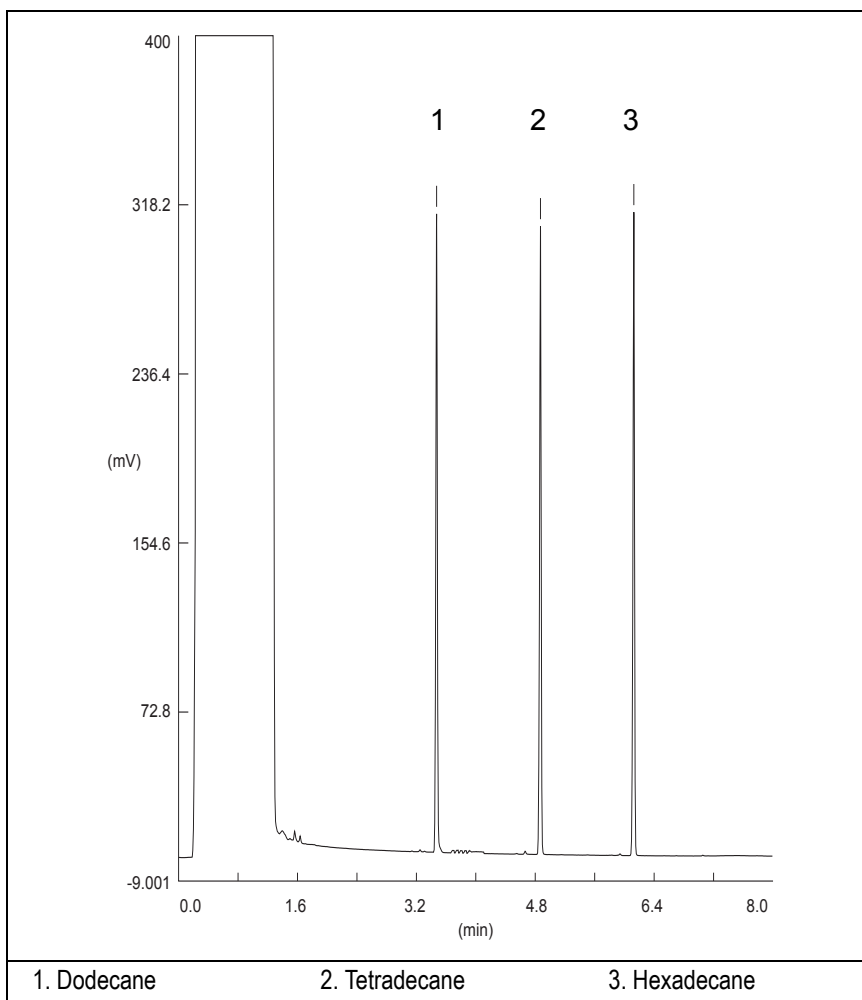





Figure 2-1. FID-Splitless Injection


13. The following acceptance criteria indicate successful completion of FID-S/SL checkout according to the data handling in use.
14. If these criteria are not met, repeat the test.


Table 2-3. FID-S/SL Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 30	< 300
	Wander (μV)	< 50	< 500
	Drift ($\mu\text{V}/\text{h}$)	< 100	< 1 000
	Analytical Results	Analog (1V Full Scale) Area Counts (0.1 μVs)	Digital (10V Full Scale) Area Counts (0.1 μVs)
	Components	> 4 000 000 for each component	> 40 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1	1 ± 0.1
			

Computing-integrator (e.g. ChromJet)	
	

CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 30
	Wander (μV)	< 50
	Drift ($\mu\text{V}/\text{h}$)	< 100
	Analytical Results (1V Full Scale) - Area Counts (0.01 μVs)	
	Components	> 40 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 300
Wander (μV)	< 500
Drift ($\mu\text{V}/\text{h}$)	< 1 000
Analytical Results (10V Full Scale) - Area Counts (μVs)	
Components	> 4 000 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 3 000
Wander Counts)	< 5 000
Drift (Counts/h)	< 10 000
Analytical Results Area Counts (Cts*s)	
Components	> 40 000 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

Analytical Acceptance Comments	
1	When the make-up gas is not used, the acceptance values will result to be 2.5 times lower than the values reported in Table 2-3.
2	When helium is used as make-up gas, the acceptance values will result to be 10 times lower than the values reported in Table 2-3.

Checkout Using FID with OC Injector

Chapter at a Glance...

SOP Number: P0293/08/E - 01 September 2009	48
Scope	48
Parts Referenced	48
Analytical Conditions Required for On-Column Injection	49
Recommended Initial Operations	50

Operating Procedures

FID-OCI Checkout in On-Column Mode.....	51
---	----

SOP Number: P0293/08/E - 01 September 2009

Scope

Use the following procedure to verify proper FID operation with the Cold On-Column Injector.

Parts Referenced

Table 3-1. FID-OCI Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film.thickness.	260 800 01
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Vespel Ferrule	Vespel Ferrule for 0.32 mm ID Column	290 134 60
Retaining Nut	M4 capillary column retaining nut	350 324 23
Syringe	10 µl size; 75 mm needle length	365 020 07
Test Mixture	Three components in n-Hexane: <i>Component</i> <i>Concentration</i> Dodecane 20 µg/ml Tetradecane 20 µg/ml Hexadecane 20 µg/ml	338 190 20
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	
Syringe	10 µl size; 80 mm needle length	365 020 19
Pre-column	2 m long; 0.53 mm ID	260 603 75
Press-fit set	Set of Press-fir connectors for TRACE OC	350 038 45

Analytical Conditions Required for On-Column Injection

Table 3-2. FID-OCI Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 35 ml/min Air = 350 ml/min Make-up Gas: Nitrogen = 30 ml/min
Oven Program	Initial Temperature = 70 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 200 °C Final Time = 1 minute
Injector	Secondary Cooling = 0.2 minutes
Detector	Base Temperature = 250 °C Detector Signal Range = 10 ⁰
Injected Volume	1 µl of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
2. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
In case of automatic On-column for TriPlus sampler, install the pre-column and connect it to the test column by press-fit connector.
3. Install and connect the TriPlus sampler and its components.
4. Perform Column Evaluation and Leak Test.
5. Connect your data handling.
Verify that your data handling is properly connected to your GC system.
6. Verify the opening/closing of the OC injector actuator by using the proper commands.
7. Verify the alignment of the syringe on the OC injector.

OPERATING PROCEDURE

FID-OCI Checkout in On-Column Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	OCI
Left carrier or Right carrier	He (helium)
Left detector or Right detector	FID

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to Pressure and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature *Temp* and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		200
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate Cold On-Column Injector Control Table. Scroll to *Sec. cool time* and set the required secondary cooling time.

LEFT INLET (OCI) ¹		
Pressure	30.0	30.0
Sec. Cool Time		0.2<

1. These settings could also be for a right inlet.

4. Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate FID Detector Control Table. Set the required temperature *Base Temp* and the detector gases *H2*, *Air* and *Mkup* (if available) required setpoints.

LEFT DETECTOR (FID) ¹		
Flame		Off
Base temp	250	250
Signal pA		(5.5)
Ign. thresh		2.0
Flameout retry		Off
H2	35	35
Air	350	350
Mkup N2	30	30<

1. These settings could also be for a right detector.

5. Ignite the FID flame scrolling to *Flame* and pressing **ON**.
6. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate FID Detector Signal Control Table. Observe the FID flame signal at the display. This is the flame-on background offset. Scroll to *Range* and set the electrometer amplifier input range required.

LEFT SIGNAL (FID) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline comp	Off

1. These settings could also be for a right signal.

7. Activate your Data System and set the parameters required for the checkout.
8. In the FID Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
9. Perform a blank analysis injecting pure hexane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 3-3 according to the data handling in use.

10. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
11. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
12. Perform the analysis

Manual injection

- Inject the test mixture and press **START** on the GC to begin the checkout .

Automatic injection with TriPlus sampler

- Fill a vial with the standard mix and place that vial in the sample tray.
- Load the method for OC and perform the sampling.

The resulting chromatogram should look like the one shown in *Figure 3.1*.

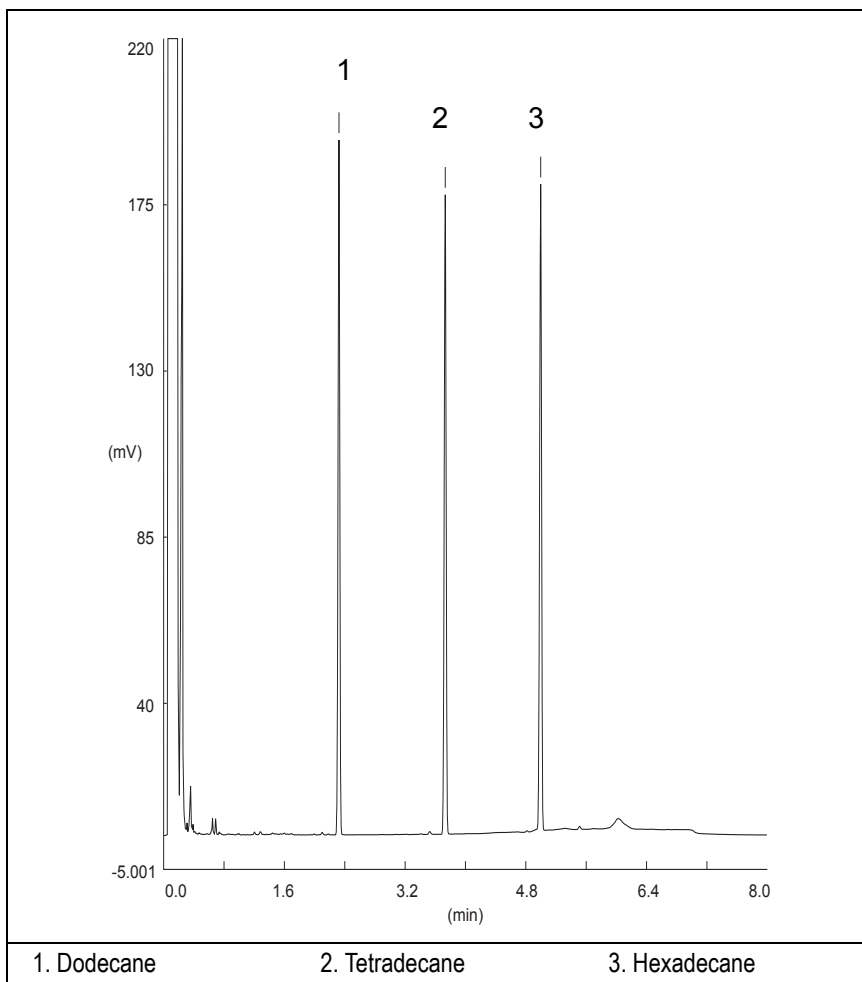





Figure 3-1. FID-On-Column Injection


13. The following criteria indicate successful completion of FID-OCI checkout.
14. If these criteria are not met, repeat the test.


Table 3-3. FID-OCI Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 30	< 300
	Wander (μV)	< 50	< 500
	Drift ($\mu\text{V}/\text{h}$)	< 100	< 1 000
	Analytical Results	Analog (1V Full Scale) Area Counts (0.1 μVs)	Digital (10V Full Scale) Area Counts (0.1 μVs)
	Components	> 2 500 000 for each component	> 25 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1	1 ± 0.1
			

Computing-integrator (e.g. ChromJet)	
	

CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 30
	Wander (μV)	< 50
	Drift ($\mu\text{V}/\text{h}$)	< 100
	Analytical Results (1V Full Scale) - Area Counts (0.01 μVs)	
	Components	> 25 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 300
Wander (μV)	< 500
Drift ($\mu\text{V}/\text{h}$)	< 1 000
Analytical Results (10V Full Scale) - Area Counts (μVs)	
Components	> 2 500 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 3 000
Wander Counts)	< 5 000
Drift (Counts/h)	< 10 000
Analytical Results Area Counts (Cts*s)	
Components	> 25 000 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

Analytical Acceptance Comments	
1	When the make-up gas is not used, the acceptance values will result to be 2.5 times lower than the values reported in Table 3-3.
2	When helium is used as make-up gas, the acceptance values will result to be 10 times lower than the values reported in Table 3-3.

4

Checkout Using FID with PKD Injector

Chapter at a Glance...

SOP Number: P0307/07/E - 01 September 2009	58
Scope	58
Parts Referenced	58
Analytical Conditions Required for Packed Injector.....	60
Recommended Initial Operations	61

Operating Procedures

FID-PKD Checkout	62
------------------------	----

SOP Number: P0307/07/E - 01 September 2009

Scope

Use the following procedure to verify proper FID operation with Packed Injector.

Parts Referenced

Table 4-1. FID-PKD Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	2 mm ID tapered	453 220 50
Liner Seal	Graphite seal for glass liner	290 334 05
Pre-column	Fused Silica Capillary Column 0.5 mt long 0.53 mm ID	260 603 75
Retaining Nut	M4 capillary column retaining nut	350 324 23
Press-fit connections	For columns 0.53/0.32 mm ID	350 438 16
Injection side adapter	For Wide bore column	347 003 03
Retaining Nut	For Injection side adapter	350 024 04
Detector side adapter	For Wide bore column	347 103 04
Ferrule	6 mm ID double brass ferrule	290 341 37
Nut	1/4" G-6 mm ID nut	350 201 18
Graphite Ferrule	Graphite ferrule for 0.53 mm ID Column Graphite ferrule for 0.32 mm ID Column	290 134 86 290 134 87
Septum	Standard septum for Packed Injector	313 032 26
Syringe	10 µl size; 50 mm needle length	365 005 25

Table 4-1. FID-PKD Parts Referenced (Continued)

Part	Description	Part Number	
Test Mixture	Three components in n-Hexane:	338 190 20	
	<i>Component</i> <i>Concentration</i>		
	Dodecane 20 µg/ml		
	Tetradecane 20 µg/ml		
Hexadecane 20 µg/ml			
Gases	Chromatographic-grade purity		
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator		

Analytical Conditions Required for Packed Injector

Table 4-2. FID-PKD Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 35 ml/min Air = 350 ml/min Make-up Gas: Nitrogen = 30 ml/min
Oven Program	Iso Temperature = 50 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 200 °C Final Time = 1 minute
Injector	Operating Mode = Packed Temperature = 200 °C
Detector	Base Temperature = 250 °C Detector Signal Range = 10 ⁰
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Insert the glass liner.
Remove the column and insert the 2 mm ID tapered glass liner, as required for the checkout, from the bottom of the injector with the appropriate liner seal. Fix the liner by using the appropriate adapter for capillary column. Refer to the TRACE GC Ultra Maintenance and Troubleshooting Manual.
2. Mount the adapter for capillary column on the detector base body.
3. Replace the septum
A new septum should be installed properly in your injector.
4. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
5. Install the precolumn.
Connect the precolumn to the injector.
6. Install the test column
Connect the test column to the precolumn by using the press fit connections provided.
7. Perform the manual leak check following the Operating Procedure “*Performing a Leak Check*” in Chapter 14 of the TRACE GC Ultra Operating Manual.
8. Connect the other end of the test column to the detector base body.
9. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

FID-PKD Checkout

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	PKD
Left carrier or Right carrier	He (helium)
Left detector or Right detector	FID

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to Pressure and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	50.0	50.0
Initial Time		1.00
Ramp 1		20.0
Final temp		200
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate PKD Injector Control Table. Set the required temperature setpoint **Temp**. Verify to operate in

Packed mode. If not, scroll to Mode, press **MODE/TYPE** to access the selection menu, then select Packed.

LEFT INLET (PKD) ¹		
Temp	200	200
Pressure	30.0	30.0
Mode:		Packed

1. These settings could also be for a right inlet.

- Use **LEFT DETECTOR** or **RIGHT DETECTOR**, entry the appropriate FID Detector Control Table. Set the required temperature Base Temp and the detector gases H₂, Air and Mkup required setpoints.

LEFT DETECTOR (FID) ¹		
Flame		Off
Base temp	250	250
Signal pA		(5.5)
Ign.thresh		2.0
Flameout retry		Off
H ₂	35	35
Air	350	350
Mkup N ₂	30	30<

1. These settings could also be for a right detector.

- Ignite the FID flame scrolling to Flame and pressing **ON**.
- Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate FID Detector Signal Control Table. Observe the FID flame signal at the display. This is the flame-on background offset. Scroll to Range and set the electrometer amplifier input range required.
- Activate your Data System and set the parameters required for the checkout.
- In the FID Detector Signal Control Table, scroll to Auto zero? and turn it **YES**.

LEFT SIGNAL (FID) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline comp	Off

1. These settings could also be for a right signal.

9. Perform a blank analysis injecting pure hexane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 4-3 according to the data handling in use.

10. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
11. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
12. Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 4.1*

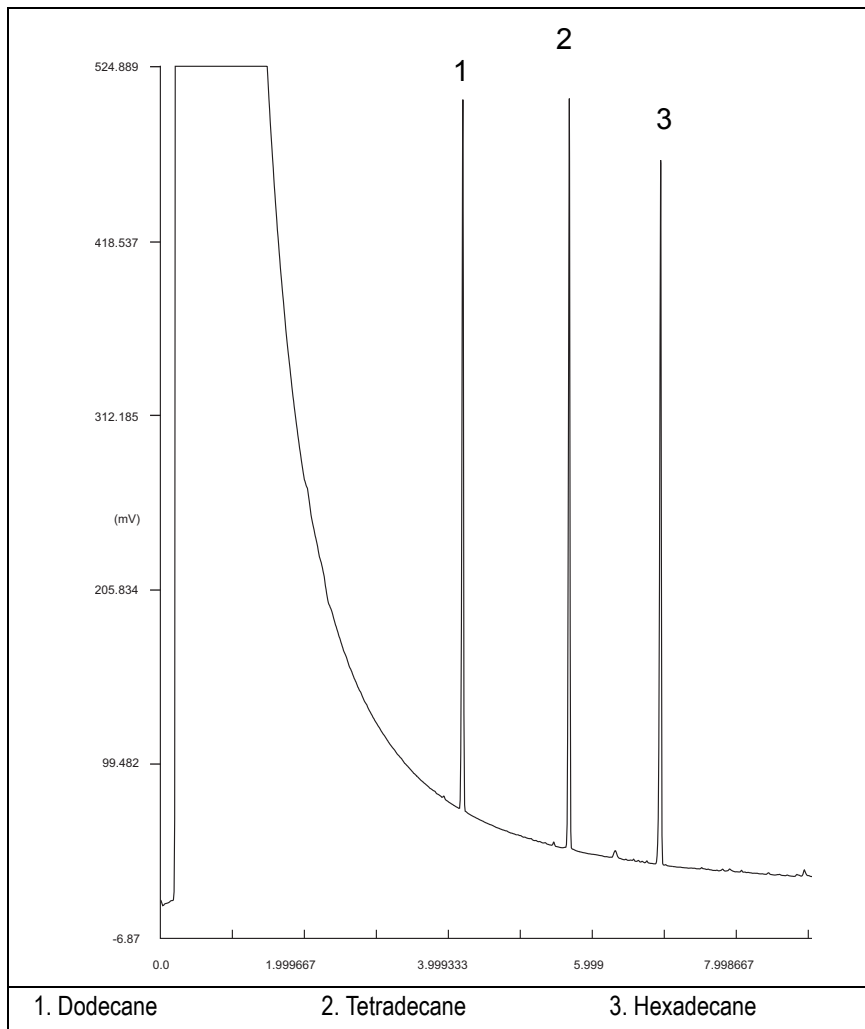





Figure 4-1. FID-Packed Injection


- 13. The following criteria indicate successful completion of FID-PKD checkout.
- 14. If these criteria are not met, repeat the test.


Table 4-3. FID-PKD Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 30	< 300
	Wander (μV)	< 50	< 500
	Drift (μV/h)	< 100	< 1 000
	Analytical Results	Analog (1V Full Scale) Area Counts (0.1 μVs)	Digital (10V Full Scale) Area Counts (0.1 μVs)
	Components	> 3 600 000 for each component	> 36 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1	1 ± 0.1
			

Computing-integrator (e.g. ChromJet)	
	

CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 30
	Wander (μV)	< 50
	Drift (μV/h)	< 100
	Analytical Results (1V Full Scale) - Area Counts (0.01 μVs)	
	Components	> 36 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 300
Wander (μV)	< 500
Drift ($\mu\text{V}/\text{h}$)	< 1 000
Analytical Results (10V Full Scale) - Area Counts (μVs)	
Components	> 3 600 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 3 000
Wander Counts)	< 5 000
Drift (Counts/h)	< 10 000
Analytical Results Area Counts (Cts*s)	
Components	> 36 000 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

Analytical Acceptance Comments	
1	When the make-up gas is not used, the acceptance values will result to be 2.5 times lower than the values reported in Table 4-3.
2	When helium is used as make-up gas, the acceptance values will result to be 10 times lower than the values reported in Table 4-3.

5

Checkout Using FID with PPKD Injector

Chapter at a Glance...

SOP Number: P0308/07/E - 01 September 2009	70
Scope	70
Parts Referenced.....	70
Analytical Conditions Required for Purged Packed Injector	72
Recommended Initial Operations	73

Operating Procedures

FID-PPKD Checkout.....	74
------------------------	----

SOP Number: P0308/07/E - 01 September 2009

Scope

Use the following procedure to verify proper FID operation with Purged Packed Injector.

Parts Referenced

Table 5-1. FID-PPKD Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	2 mm ID tapered	453 220 50
Liner Seal	Graphite seal for glass liner	290 334 05
Pre-column	Fused Silica Capillary Column 0.5 mt long 0.53 mm ID	260 603 75
Retaining Nut	M4 capillary column retaining nut	350 324 23
Press-fit connections	For columns 0.53/0.32 mm ID	350 438 16
Injection side adapter	For Wide bore column	347 003 03
Retaining Nut	For Injection side adapter	350 024 04
Detector side adapter	For Wide bore column	347 103 04
Ferrule	6 mm ID double brass ferrule	290 341 37
Nut	1/4" G-6 mm ID nut	350 201 18
Graphite Ferrule	Graphite ferrule for 0.53 mm ID Column Graphite ferrule for 0.32 mm ID Column	290 134 86 290 134 87
Septum	Standard septum for Purged Packed Injector	313 032 26
Syringe	10 µl size; 50 mm needle length	365 005 25

Table 5-1. FID-PPKD Parts Referenced (Continued)

Part	Description	Part Number	
Test Mixture	Three components in n-Hexane:	338 190 20	
	<i>Component</i> <i>Concentration</i>		
	Dodecane 20 µg/ml		
	Tetradecane 20 µg/ml		
Hexadecane 20 µg/ml			
Gases	Chromatographic-grade purity		
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator		

Analytical Conditions Required for Purged Packed Injector

Table 5-2. FID-PPKD Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 35 ml/min Air = 350 ml/min Make-up Gas: Nitrogen = 30 ml/min
Oven Program	Initial Temperature = 50 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 200 °C Final Time = 1 minute
Injector	Operating Mode = Wide bore Temperature = 200 °C
Detector	Base Temperature = 250 °C Detector Signal Range = 10 ⁰
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Insert the glass liner.
Remove the column and insert the 2 mm ID tapered glass liner, as required for the checkout, from the bottom of the injector with the appropriate liner seal. Fix the liner by using the appropriate adapter for capillary column. Refer to the TRACE GC Ultra Maintenance and Troubleshooting Manual.
2. Mount the adapter for capillary column on the detector base body.
3. Replace the septum
A new septum should be installed properly in your injector.
4. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
5. Install the precolumn.
Connect the precolumn to the injector.
6. Install the test column
Connect the test column to the precolumn by using the press fit connections provided.
7. Perform the manual leak check following the Operating Procedure “*Performing a Leak Check*” in Chapter 14 of the TRACE GC Ultra Operating Manual.
8. Connect the other end of the test column to the detector base body.
9. Perform Column Evaluation.
10. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

FID-PPKD Checkout

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	PPKD
Left carrier or Right carrier	He (helium)
Left detector or Right detector	FID

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select con pres. Scrool to Pressure and set the pressure value to have the required carrier gas flow rate Col. flow.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	50.0	50.0
Initial Time		1.00
Ramp 1		20.0
Final temp		200
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate PPKD Injector Control Table. Set the required temperature setpoint Temp. Verify to operate in

Wide bore mode. If not, scroll to *Mode*, press **MODE/TYPE** to access the selection menu, then select *Widebore*.

LEFT INLET (PPKD) ¹		
Temp	200	200
Pressure	30.0	30.0
Mode:	Wide bore<	
Constant sept purge?		Y<

1. These settings could also be for a right inlet.

- Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate FID Detector Control Table. Set the required temperature *Base Temp*, and the detector gases *H2*, *Air* and *Mkup* required setpoints.

LEFT DETECTOR (FID) ¹		
Flame		Off
Base temp	250	250
Signal pA		(5.5)
Ign.thresh		2.0
Flameout retry		Off
H2	35	35
Air	350	350
Mkup N2	30	30<

1. These settings could also be for a right detector.

- Ignite the FID flame scrolling to *Flame* and pressing **ON**.
- Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate FID Detector Signal Control Table. Observe the FID flame signal at the display. This is the flame-on background offset. Scroll to *Range* and set the electrometer amplifier input range required.

LEFT SIGNAL (FID) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline comp	Off

1. These settings could also be for a right signal.

7. Activate your Data System and set the parameters required for the checkout.
8. In the FID Detector Signal Control Table, scroll to `Auto zero?` and turn it **YES**.
9. Perform a blank analysis injecting pure hexane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 5-3 according to the data handling in use.

10. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
11. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
12. Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 5.1*

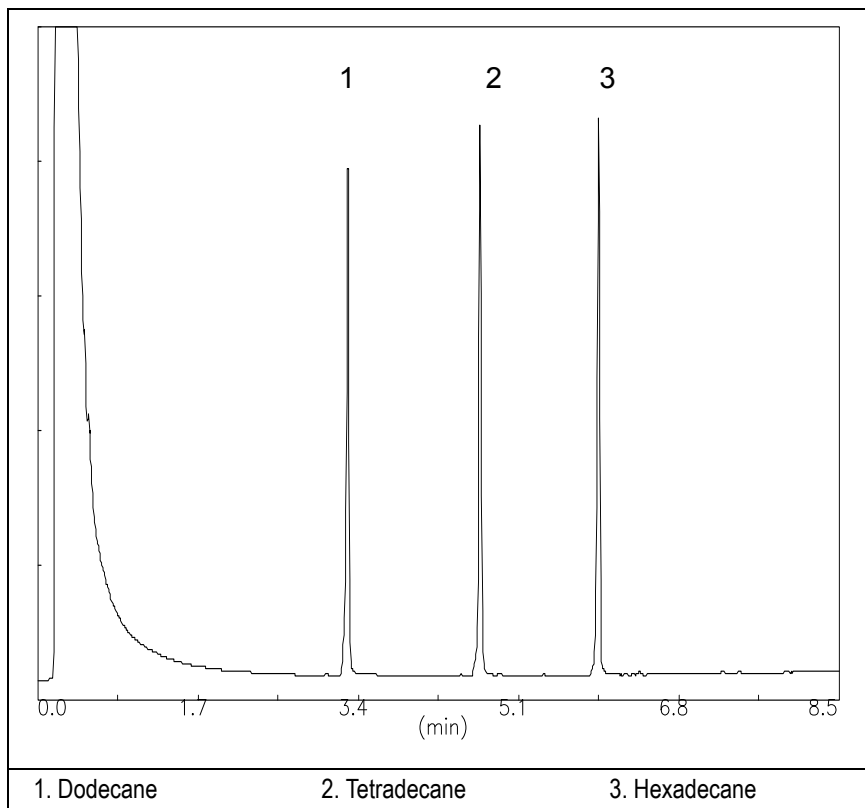





Figure 5-1. FID-PPKD Injection


13. The following criteria indicate successful completion of FID-PPKD checkout.
14. If these criteria are not met, repeat the test.


Table 5-3. FID-PPKD Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 30	< 300
	Wander (μV)	< 50	< 500
	Drift (μV/h)	< 100	< 1 000
	Analytical Results	Analog (1V Full Scale) Area Counts (0.1 μVs)	Digital (10V Full Scale) Area Counts (0.1 μVs)
	Components	> 3 600 000 for each component	> 36 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1	1 ± 0.1
			

Computing-integrator (e.g. ChromJet)	
	

CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 30
	Wander (μV)	< 50
	Drift (μV/h)	< 100
	Analytical Results (1V Full Scale) - Area Counts (0.01 μVs)	
	Components	> 36 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 300
Wander (μV)	< 500
Drift ($\mu\text{V/h}$)	< 1 000
Analytical Results (10V Full Scale) - Area Counts (μVs)	
Components	> 3 600 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 3 000
Wander Counts)	< 5 000
Drift (Counts/h)	< 10 000
Analytical Results Area Counts (Cts*s)	
Components	> 36 000 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

Analytical Acceptance Comments	
1	When the make-up gas is not used, the acceptance values will result to be 2.5 times lower than the values reported in Table 5-3.
2	When helium is used as make-up gas, the acceptance values will result to be 10 times lower than the values reported in Table 5-3.

6

Checkout Using FID with PTV Injector

Chapter at a Glance...

SOP Number: P0309/07/E - 01 September 2009	82
Scope	82
Parts Referenced	82
Analytical Conditions Required for PTV Splitless Injection	83
Recommended Initial Operations	84

Operating Procedures

FID-PTV Checkout in PTV Splitless Mode	85
--	----

SOP Number: P0309/07/E - 01 September 2009

Scope

Use the following procedure to verify proper FID operation with the Programmable Temperature Vaporizing Injector.

Parts Referenced

Table 6-1. FID-PTV Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Liner	Silcosteel 2 mm ID	453 220 44
Liner Seal	Graphite seal for liner	290 034 17
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
Septum	Standard septum for PTV injector (set of 10)	313 132 25
Syringe	10 µl size; 50 mm needle length	365 005 25
Test Mixture	Three components in n-Hexane: <i>Component</i> <i>Concentration</i> Dodecane 20 µg/ml Tetradecane 20 µg/ml Hexadecane 20 µg/ml	338 190 20
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for PTV Splitless Injection

Table 6-2. FID-PTV Analytical Conditions

Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 35 ml/min Air = 350 ml/min Make-up Gas: Nitrogen = 30 ml/min
Oven Program	Initial Temperature = 50 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 200 °C Final Time = 1 minute
Injector	Operating Mode = PTV splitless Splitless Time = 0.8 minutes Split Flow = 50 ml/min Constant Septum Purge = Yes Inject Temp = 50 °C Inject Time = 0.1 minute Transfer ramp = 10 °C/sec Transfer Temperature = 260 °C Transfer time = 1 minute
Detector	Base Temperature = 250 °C Detector Signal Range = 10 ⁰
Injected Volume	1 µl of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Replace the liner.
The liner currently installed in your injector should be carefully removed and replaced with the 2 mm ID Silcosteel liner, as required for the checkout, with the appropriate liner seal.
2. Replace the septum
A new septum should be installed properly in your injector.
3. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
4. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
5. Perform Column Evaluation and Leak Test.
6. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

FID-PTV Checkout in PTV Splitless Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Right inlet	PTV
Right carrier	He (helium)
Right detector	FID

1. Use **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to **Flow** mode, press **MODE/TYPE** to access the selection menu, then select **con pres**. Scroll to **Pressure** and set the pressure value to have the required carrier gas flow rate **Col.flow**.

RIGHT CARRIER ¹		
Pressure	30.0	30.0
Col.flow	3.00	
Lin. veloc.		(60.9) <

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	50.0	50.0
Initial Time		1.00
Ramp 1		20.0
Final temp		200
Final time 1		1.00 <
Ramp 2		Off

- Use **RIGHT INLET** to display the appropriate Programmable Temperature Vaporizing Injector Control Table. Set the required temperature setpoint Temp. Verify to operate in **PTV splitless** mode. If not, scroll to Mode, press **MODE/TYPE** to access the selection menu, then select PTV splitless. Scroll to Splitless time to set the required setpoint.

RIGHT INLET (PTV)		
Temp	50	50
Pressure	30.0	30.0
Mode:	PTV Splitless	
Total flow	(53.0)	
Split Flow	50.0	50.0
Splitless time	0.80	0.80
Constant sept purge?		Y
Inject phase menu:		Y<

- Scroll to Inject phase menu. Press **MODE/TYPE** to enter the **PTV Phase Menu**.

PTV PHASE MENU	
Ramped pressure?	N
Inject temp	50
Inject time	0.1
Transfer ramp	10
Transfer temp	260
Transfer time	1.00<

- Select Ramped pressure? **NO**. Set the required Inject temp and *Inject time* setpoints as required. Then, set the Transfer ramp, the Transfer temp and the Transfer time required setpoints.
- Use **RIGHT DETECTOR** to display the appropriate FID Detector Control Table. Set the required temperature Base Temp. and the detector gases H2, Air and Mkup required setpoints.

7. Ignite the FID flame scrolling to Flame and pressing **ON**.

RIGHT DETECTOR (FID)		
Flame		Off
Base temp	250	250
Signal pA		(5.5)
Ign.thresh		2.0
Flameout retry		Off
H2	35	35
Air	350	350
Mkup N2	30	30<

8. Use **RIGHT SIGNAL** to display the appropriate FID Detector Signal Control Table. Observe the FID flame signal at the display. This is the flame-on background offset. Scroll to Range and set the electrometer amplifier input range required.

RIGHT SIGNAL (FID)	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline comp	Off

9. Activate your Data System and set the parameters required for the checkout.
10. In the FID Detector Signal Control Table, scroll to Auto zero? and turn it **ON**.
11. Activate your Data System and set the parameters required for the checkout.
12. In the FID Detector Signal Control Table, scroll to Auto zero? and turn it **YES**.

13. Perform a blank analysis injecting pure hexane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



Refer to the Acceptance Values reported in the Table 6-3 according to the data handling in use.

14. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
15. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
16. Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 6.1*

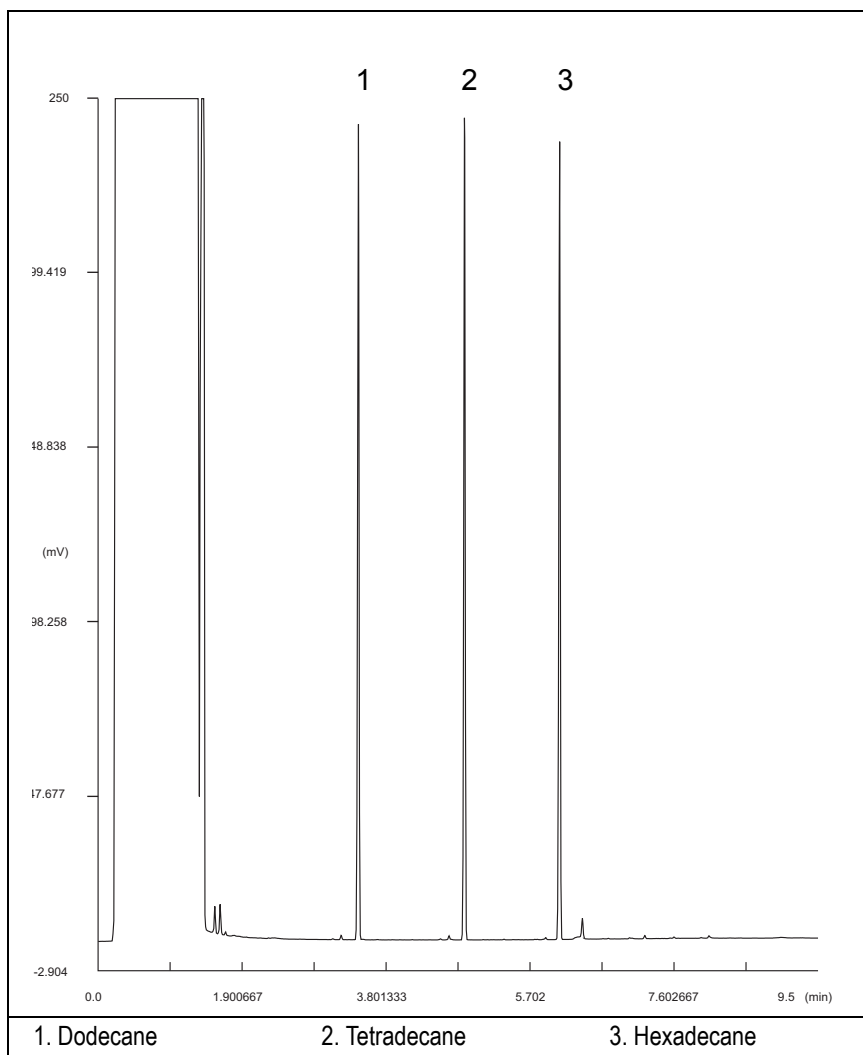





Figure 6-1. FID-PTV Injection


17. The following criteria indicate successful completion of FID-PTV checkout.
18. If these criteria are not met, repeat the test.


Table 6-3. FID-PPKD Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 30	< 300
	Wander (μV)	< 50	< 500
	Drift (μV/h)	< 100	< 1 000
	Analytical Results	Analog (1V Full Scale) Area Counts (0.1 μVs)	Digital (10V Full Scale) Area Counts (0.1 μVs)
	Components	> 2 500 000 for each component	> 25 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1	1 ± 0.1
			

Computing-integrator (e.g. ChromJet)	
	

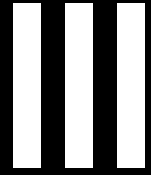
CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 30
	Wander (μV)	< 50
	Drift (μV/h)	< 100
	Analytical Results (1V Full Scale) - Area Counts (0.01 μVs)	
	Components	> 25 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 300
Wander (μV)	< 500
Drift ($\mu\text{V}/\text{h}$)	< 1 000
Analytical Results (10V Full Scale) - Area Counts (μVs)	
Components	> 2 500 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

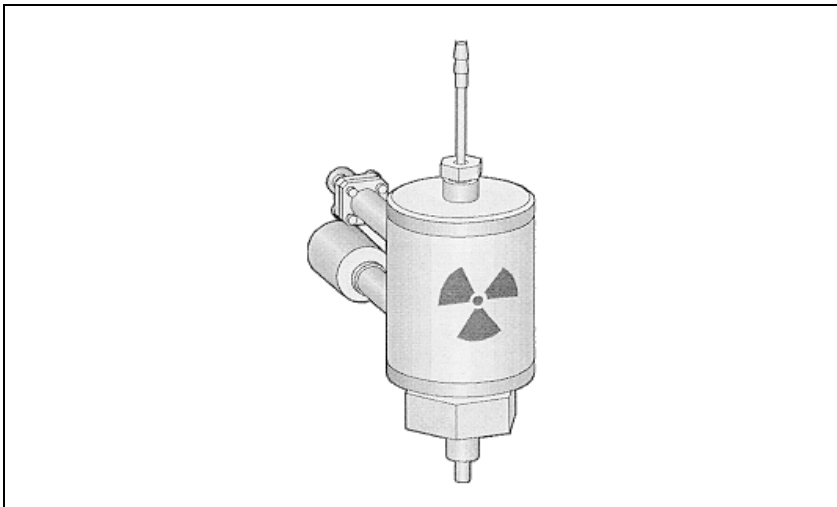
XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 3 000
Wander Counts)	< 5 000
Drift (Counts/h)	< 10 000
Analytical Results Area Counts (Cts*s)	
Components	> 25 000 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

Analytical Acceptance Comments	
1	When the make-up gas is not used, the acceptance values will result to be 2.5 times lower than the values reported in Table 6-3.
2	When helium is used as make-up gas, the acceptance values will result to be 10 times lower than the values reported in Table 6-3.

SECTION



SOPs Using ECD



The *SOPs Using ECD* section, contains the procedures to test the TRACE GC Ultra with the Electron Capture Detector (ECD) using different injectors.

Chapter 7 *Checkout Using ECD with S/SL Injector.*

Chapter 8 *Checkout Using ECD with OC Injector.*

Chapter 9 *Checkout Using ECD with PKD Injector*

Chapter 10 *Checkout Using ECD with PPKD Injector.*

Chapter 11 *Checkout Using ECD with PTV Injector.*



Checkout Using ECD with S/SL Injector

Chapter at a Glance...

SOP Number: P0294/07/E - 01 September 2009	96
Scope	96
Parts Referenced	96
Analytical Conditions Required for Splitless Injection	97
Recommended Initial Operations	98

Operating Procedures

ECD-S/SL Checkout in Splitless Mode	99
---	----

SOP Number: P0294/07/E - 01 September 2009

Scope

Use the following procedure to verify proper ECD operation with the Split/Splitless Injector.

Parts Referenced

Table 7-1. ECD-S/SL Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long; 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	3 mm ID for splitless injections	453 200 32
Liner Seal	Graphite seal glass liner	290 334 06
Graphite Ferrule	Graphite Ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
Septum	Standard septum for S/SL injector	313 032 11
Syringe	10 µl size; 70 mm needle length	365 001 03
Test Mixture	Two components in Iso-octane <i>Component Concentration</i> Lindane 0.030 µg/ml Aldrin 0.030 µg/ml	338 190 11
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for Splitless Injection

Table 7-2. ECD-S/SL Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Make-up: Nitrogen = 30 ml/min
Oven Program	Initial Temperature = 70 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 220 °C Final Time = 1 minute
Injector	Operating Mode = Splitless Temperature = 230 °C Splitless Time = 0.8 minutes Split Flow = 60 ml/min Constant Septum Purge = Yes
Detector	Base Temperature = 250 °C ECD temperature = 300 °C Reference Current = 1 nA Pulse Amplitude = 50 V Pulse Width = 1µs
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Replace the glass liner.
The glass liner currently installed in your injector should be carefully removed and replaced with the 3 mm ID glass liner for splitless application, as required for the checkout, with the appropriate liner seal.
2. Replace the septum
A new septum should be installed properly in your injector.
3. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
4. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
5. Perform Column Evaluation and Leak Test.
6. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

ECD-S/SL Checkout in Splitless Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	S/SL
Left carrier or Right carrier	He (helium)
Left detector or Right detector	ECD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to Pressure and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		220
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate Split/Splitless Injector Control Table. Set the required temperature setpoint *Temp*. Verify to

operate in **Splitless** mode. If not, scroll to **Mode**, press **MODE/TYPE** to access the selection menu, then select **Splitless**. Scroll to **Splitless time** to set the required setpoint.

LEFT INLET (S/SL) ¹		
Temp	230	230
Pressure	30.0	30.0
Mode:	Splitless	
Total flow	(63.0)	
Split Flow	60.0	60.0
Splitless time	0.80	0.80
Constant sept purge?		Y<

1. These settings could also be for a right inlet.

4. Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate ECD Detector Control Table. Set the required temperature **Base Temp.** and the **Mkup** gas required setpoints.

LEFT DETECTOR (ECD) ¹		
Base temp	250	250
ECD Temp	300	300
Ref current nA	1.0	
Freq kHz	(2.20)	
Pulse amp V	50	
Pulse width μ s	1.0	
Mkup (N2)	30	30<

1. These settings could also be for a right detector.

5. Set the **Reference Current** to 1.0 nA.
6. Set the **Pulse Amplitude** to 50 V.
7. Scroll to **Pulse Width** and press **ENTER** to open the menu selection. Select the pulse width to 1 μ s then press **ENTER**.

8. Observe the ECD frequency value at the display. A base frequency value between 1 kHz and 3 kHz should be displayed.

If the ECD frequency value is less than 1kHz, reduce gradually the pulse amplitude till the base frequency value is about 1 kHz (the pulse amplitude value should not be less than 20V). Then, if necessary modify the reference current in order to have a base frequency of about 1-1,5 kHz. Let the detector signal to stabilize.

9. Activate your Data System and set the parameters required for the checkout.
10. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the ECD Detector Signal Control Table. Scroll to `Auto zero?` and turn it **YES**.

LEFT SIGNAL (ECD) ¹	
Output	(1000)
Offset	100<
Auto zero?	Y/N
Baseline comp?	Off

1. These settings could also be for a right signal.

11. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 7-3 according to the data handling in use.

12. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
13. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
14. When the GC is ready, inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 7.1*.

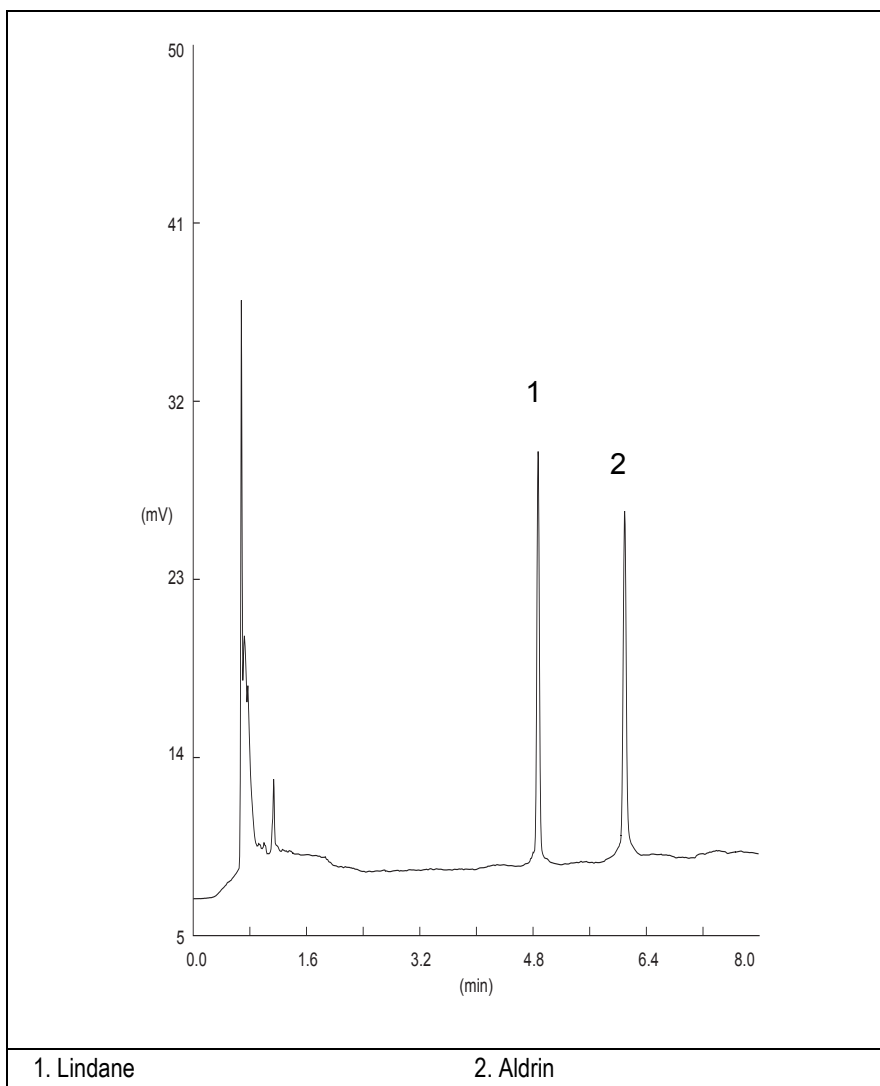


Figure 7-1. ECD-Splitless Injection

15. Establish the integration parameters and the peak table identifying the test mix components.
16. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Lindane component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Lindane.
- Repeat the procedure to calculate the signal-to-noise ratio also for Aldrin.
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Aldrin.



NOTE

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows:

Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:

- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.




CAUTION


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


17. The following criteria indicate successful completion of ECD-S/SL checkout.


18. If these criteria are not met, repeat the test.

Table 7-3. ECD-S/SL Acceptance Criteria

CHROM-CARD				
Acceptance Values	Baseline Parameters		Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)		< 40	< 400
	Wander (μV)		< 80	< 800
	Drift ($\mu\text{V/h}$)		< 200	< 2 000
	Analytical Results			
	Lindane Signal-to-Noise Ratio			> 4 000
	Aldrin Signal-to-Noise Ratio			> 4 000
				

CHROMQUEST			
Acceptance Values	Baseline Parameters (1V Full Scale)		
	Noise (μV)		< 40
	Wander (μV)		< 80
	Drift ($\mu\text{V/h}$)		< 200
	Analytical Results		
	Lindane Signal-to-Noise Ratio		> 4 000
	Aldrin Signal-to-Noise Ratio		> 4 000
			

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 400
Wander (μV)	< 800
Drift ($\mu\text{V/h}$)	< 2 000
Analytical Results	
Lindane Signal-to-Noise Ratio	> 4 000
Aldrin Signal-to-Noise Ratio	> 4 000
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 4 000
Wander (Counts)	< 8 000
Drift (Counts/h)	< 20 000
Analytical Results	
Lindane Signal-to-Noise Ratio	> 4 000
Aldrin Signal-to-Noise Ratio	> 4 000
	

Analytical Acceptance Comments	
1	Using Chrom-Card, set the signal-to-noise report parameters as described in the current procedure.
2	Using ChromQuest, Atlas, Xcalibur or a Computing integrator (e.g. ChromJet), calculate the S/N ratio as <i>Peak Height (counts)/noise (counts)</i> .

Checkout Using ECD with OC Injector

Chapter at a Glance...

SOP Number: P0295/08/E - 01 September 2009	108
Scope	108
Parts Referenced	108
Analytical Conditions Required for On-Column Injection	109
Recommended Initial Operations	110

Operating Procedures

ECD-OCI Checkout in On-Column Mode	111
--	-----

SOP Number: P0295/08/E - 01 September 2009

Scope

Use the following procedure to verify proper ECD operation with the On-Column Injector.

Parts Referenced

Table 8-1. ECD-OCI Parts Referenced

	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long; 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
Vespel Ferrule	Vespel ferrule for 0.32 mm ID Column	290 134 60
Syringe	10 µl size; 75 mm needle length	365 020 07
Test Mixture	Two components in Iso-octane <i>Component Concentration</i> Lindane 0.030 µg/ml Aldrin 0.030 µg/ml	338 190 11
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	
Syringe	10 µl size; 80 mm needle length	365 020 19
Pre-column	2 m long; 0.53 mm ID	260 603 75
Press-fit set	Set of Press-fir connectors for TRACE OC	350 038 45

Analytical Conditions Required for On-Column Injection

Table 8-2. ECD-OCI Analytical Conditions

Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Make-up: Nitrogen = 30 ml/min
Oven Program	Initial Temperature = 85 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 220 °C Final Time = 1 minute
Injector	Secondary Cooling = 3 seconds
Detector	Base Temperature = 250 °C ECD Temperature = 300 °C Reference Current = 1 nA Pulse Amplitude = 50 V Pulse Width = 1µs
Injected Volume	1 µl of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Connect the required gas lines.
Verify the required gas supplies are properly connected to your GC.
2. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
In case of automatic On-column for TriPlus sampler, install the pre-column and connect it to the test column by press-fit connector.
3. Install and connect the TriPlus sampler and its components.
4. Perform Column Evaluation and Leak Test.
5. Connect your data handling.
Verify that your data handling is properly connected to your GC system.
6. Verify the opening/closing of the OC injector actuator by using the proper commands.
7. Verify the alignment of the syringe on the OC injector.

OPERATING PROCEDURE

ECD-OCI Checkout in On-Column Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	OCI
Left carrier or Right carrier	He (helium)
Left detector or Right detector	ECD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to **Pressure** and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	85.0	85.0
Initial Time		1.00
Ramp 1		20.0
Final temp		220
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate Cold On-Column Injector Control Table. Scroll to **Sec. cool time** and set the required secondary cooling time.

LEFT INLET (OCI) ¹		
Pressure	30.0	30.0
Sec. Cool Time		3.00<

1. These settings could also be for a right inlet.

4. Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate ECD Detector Control Table. Set the required temperature Base Temp. and the Mkup gas required setpoints.

LEFT DETECTOR (ECD) ¹		
Base temp	250	250
ECD Temp	300	300
Ref current nA		1.0
Freq kHz		(2.20)
Pulse amp V		50
Pulse width μ s		1.0
Mkup (N2)	30	30<

1. These settings could also be for a right detector.

5. Set the Reference Current to 1.0 nA.
6. Set the Pulse Amplitude to 50 V.
7. Scroll to Pulse Width and press **ENTER** to open the menu selection. Select the pulse width to 1 μ s then press **ENTER**.
8. Observe the ECD frequency value at the display. A base frequency value between 1 kHz and 3 kHz should be displayed. If the ECD frequency value is less than 1kHz, reduce gradually the pulse amplitude till the base frequency value is about 1 kHz (the pulse amplitude value should not be less than 20V). Then, if necessary modify the reference current in order to have a base frequency of about 1-1,5 kHz. Let the detector signal to stabilize.
9. Activate your Data System and set the parameters required for the checkout.
10. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate ECD Detector Signal Control Table. Scroll to Auto zero? and turn it **YES**.

LEFT SIGNAL (ECD) ¹	
Output	(1000)
Offset	100<
Auto zero?	Y/N
Baseline comp?	Off

1. These settings could also be for a right signal.

11. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 8-3 according to the data handling in use.

12. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
13. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
14. Perform the analysis

Manual injection

- Inject the test mixture and press **START** on the GC to begin the checkout run.

Automatic injection with TriPlus sampler

- Fill a vial with the standard mix and place that vial in the sample tray.
- Load the method for OC and perform the sampling.

The resulting chromatogram should look like the one shown in *Figure 8.1*.

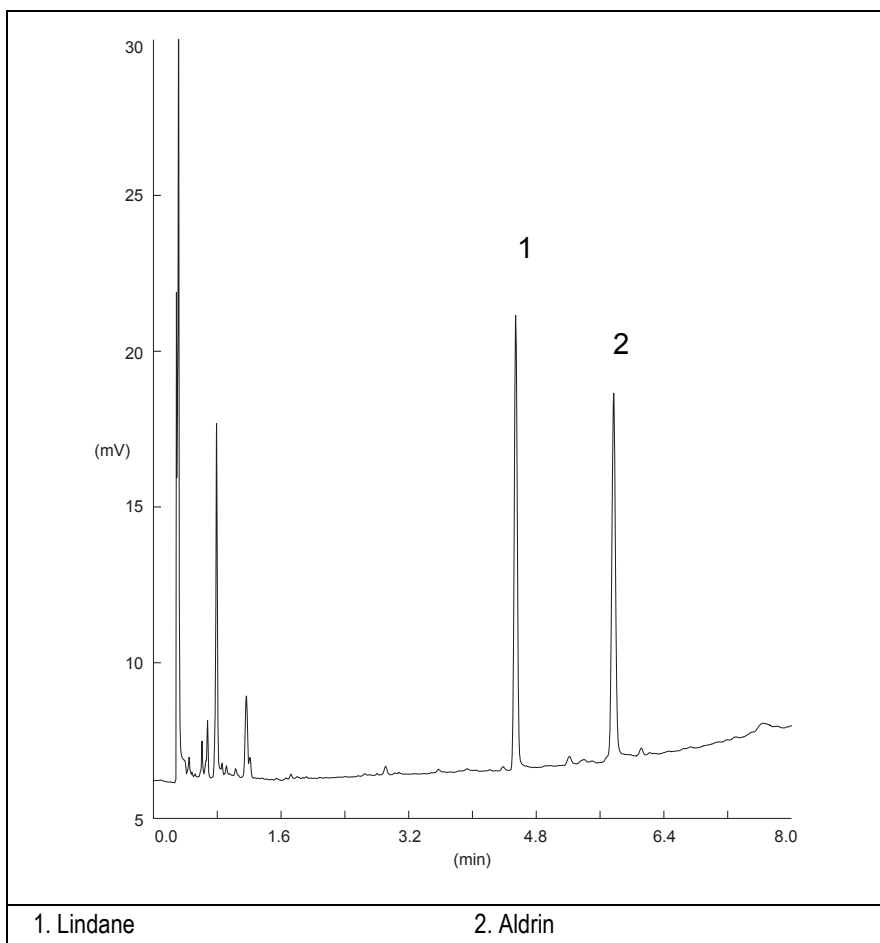


Figure 8-1. ECD-On-Column Injection

15. Establish the integration parameters and the peak table identifying the test mix components.
16. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Lindane component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Lindane.
- Repeat the procedure to calculate the signal-to-noise ratio also for Aldrin.
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Aldrin.

**NOTE**

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows:

Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:


- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.


**CAUTION**


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


17. The following criteria indicate successful completion of ECD-OCI checkout.
18. If these criteria are not met, repeat the test.

Table 8-3. ECD-OCI Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 40	< 400
	Wander (μV)	< 80	< 800
	Drift ($\mu\text{V/h}$)	< 200	< 2 000
	Analytical Results		
	Lindane Signal-to-Noise Ratio		> 3 000
	Aldrin Signal-to-Noise Ratio		> 3 000
			

CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 40
	Wander (μV)	< 80
	Drift ($\mu\text{V/h}$)	< 200
	Analytical Results	
	Lindane Signal-to-Noise Ratio	> 3 000
	Aldrin Signal-to-Noise Ratio	> 3 000
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 400
Wander (μV)	< 800
Drift ($\mu\text{V/h}$)	< 2 000
Analytical Results	
Lindane Signal-to-Noise Ratio	> 3 000
Aldrin Signal-to-Noise Ratio	> 3 000
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 4 000
Wander (Counts)	< 8 000
Drift (Counts/h)	< 20 000
Analytical Results	
Lindane Signal-to-Noise Ratio	> 3 000
Aldrin Signal-to-Noise Ratio	> 3 000
	

Analytical Acceptance Comments	
1	Using Chrom-Card, set the signal-to-noise report parameters as described in the current procedure.
2	Using ChromQuest, Atlas, Xcalibur or a Computing integrator (e.g. ChromJet), calculate the S/N ratio as <i>Peak Height (counts)/noise (counts)</i> .

Checkout Using ECD with PKD Injector

Chapter at a Glance...

SOP Number: P0310/07/E - 01 September 2009	120
Scope	120
Parts Referenced	120
Analytical Conditions Required for Packed Injector.....	122
Recommended Initial Operations	123

Operating Procedures

ECD-PKD Checkout	124
------------------------	-----

SOP Number: P0310/07/E - 01 September 2009

Scope

Use the following procedure to verify proper ECD operation with Packed Injector.

Parts Referenced

Table 9-1. ECD-PKD Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	2 mm ID tapered	453 220 50
Liner Seal	Graphite seal for glass liner	290 334 05
Pre-column	Fused Silica Capillary Column 0.5 mt long 0.53 mm ID	260 603 75
Retaining Nut	M4 capillary column retaining nut	350 324 23
Press-fit connections	For columns 0.53/0.32 mm ID	350 438 16
Injection side adapter	For Wide bore column	347 003 03
Retaining Nut	For Injection side adapter	350 024 04
Detector side adapter	For Wide bore column	347 103 04
Ferrule	6 mm ID double brass ferrule	290 341 37
Nut	1/4" G-6 mm ID nut	350 201 18
Graphite Ferrule	Graphite ferrule for 0.53 mm ID Column Graphite ferrule for 0.32 mm ID Column	290 134 86 290 134 87
Septum	Standard septum for Packed Injector	313 032 26
Syringe	10 µl size; 50 mm needle length	365 005 25

Table 9-1. ECD-PKD Parts Referenced (Continued)

Part	Description	Part Number	
Test Mixture	Two components in Iso-octane	338 190 11	
	<i>Component</i> <i>Concentration</i>		
	Lindane 0.030 µg/ml		
	Aldrin 0.030 µg/ml		
Gases	Chromatographic-grade purity		
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator		

Analytical Conditions Required for Packed Injector

Table 9-2. ECD-PKD Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Make-up Gas: Nitrogen = 30 ml/min
Oven Program	IsoTemperature = 140 °C Initial Time = 10 minute
Injector	Operating Mode = Packed Temperature = 180 °C
Detector	Base Temperature = 250 °C ECD temperature = 300 °C Reference Current = 1 nA Pulse Amplitude = 50 V Pulse Width = 1µs
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Insert the glass liner.
Remove the column and insert the 2 mm ID tapered glass liner, as required for the checkout, from the bottom of the injector with the appropriate liner seal. Fix the liner by using the appropriate adapter for capillary column. Refer to the TRACE GC Ultra Maintenance and Troubleshooting Manual.
2. Mount the adapter for capillary column on the detector base body.
3. Replace the septum
A new septum should be installed properly in your injector.
4. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
5. Install the precolumn.
Connect the precolumn to the injector.
6. Install the test column
Connect the test column to the precolumn by using the press fit connections provided.
7. Perform the manual leak check following the Operating Procedure “*Performing a Leak Check*” in Chapter 14 of the TRACE GC Ultra Operating Manual.
8. Connect the other end of the test column to the detector base body.
9. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

ECD-PKD Checkout

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	PKD
Left carrier or Right carrier	He (helium)
Left detector or Right detector	ECD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select con pres. Scrool to Pressure and set the pressure value to have the required carrier gas flow rate Col. flow.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	140	140
Initial Time		10.0
Ramp 1		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate PKD Injector Control Table. Set the required temperature setpoint Temp. Verify to operate in **Packed** mode. If not, scroll to Mode, press **MODE/TYPE** to access the selection menu, then select Packed.

LEFT INLET (PKD) ¹		
Temp	180	180
Pressure	30.0	30.0
Mode:	Packed	

1. These settings could also be for a right inlet.

- Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate ECD Detector Control Table. Set the required temperature Base Temp and the Mkup gas required setpoints.

LEFT DETECTOR (ECD) ¹		
Base temp	250	250
ECD Temp	300	300
Ref current nA		1.0
Freq kHz		(2.20)
Pulse amp V		50
Pulse width μ s		1.0
Mkup (N2)	30	30<

1. These settings could also be for a right detector.

- Set the Reference Current to 1.0 nA.
- Set the Pulse Amplitude to 50 V.
- Scroll to Pulse Width and press **ENTER** to open the menu selection. Select the pulse width to 1 μ s then press **ENTER**.
- Observe the ECD frequency value at the display. A base frequency value between 1 kHz and 3 kHz should be displayed. If the ECD frequency value is less than 1kHz, reduce gradually the pulse amplitude till the base frequency value is about 1 kHz (the pulse amplitude value should not be less than 20V). Then, if necessary modify the reference current in order to have a base frequency of about 1-1,5 kHz. Let the detector signal to stabilize.
- Activate your Data System and set the parameters required for the checkout.
- Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the ECD Detector Signal Control Table. Scroll to Auto zero? and turn it **YES**.

LEFT SIGNAL (ECD) ¹	
Output	(1000)
Offset	100<
Auto zero?	Y/N
Baseline comp?	Off

1. These settings could also be for a right signal.

11. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 9-3 according to the data handling in use.

12. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
13. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
14. When the GC is ready, inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 9.1*.

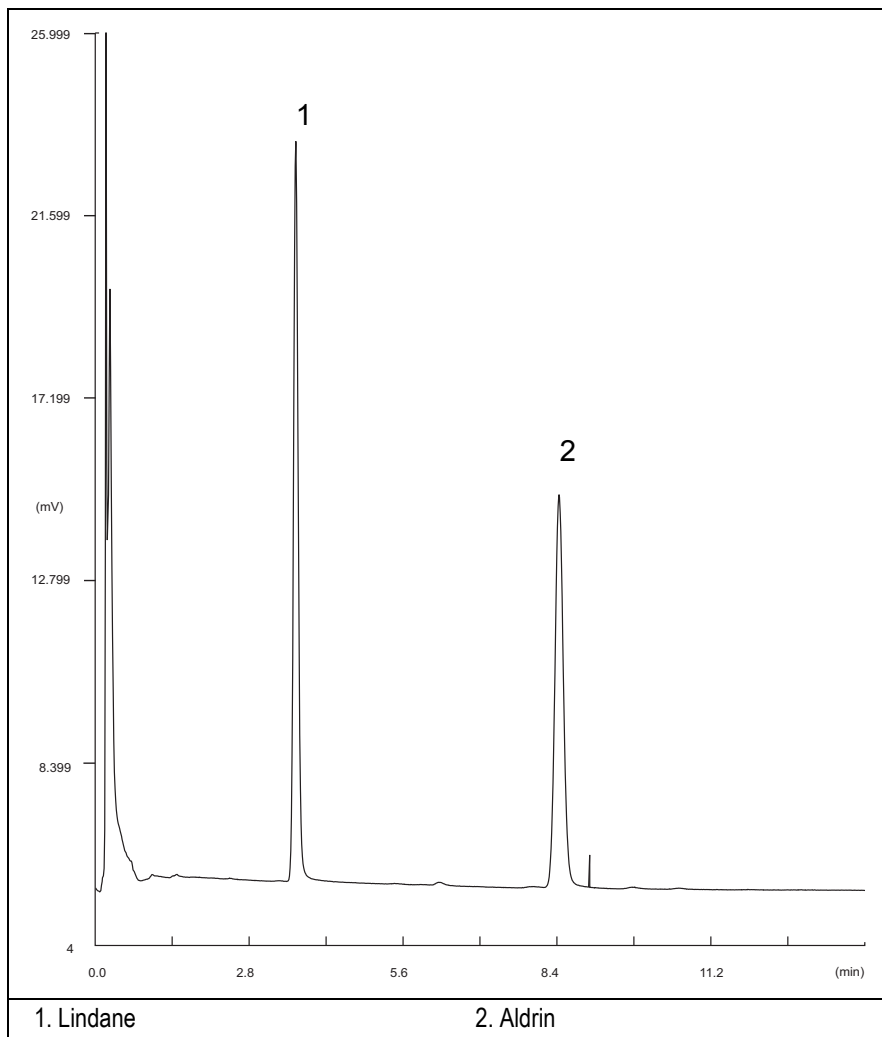


Figure 9-1. ECD-PKD Injection

15. Establish the integration parameters and the peak table identifying the test mix components.
16. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Lindane component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Lindane.
- Repeat the procedure to calculate the signal-to-noise ratio also for Aldrin.
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Aldrin.



NOTE

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows:

Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:

- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.




CAUTION


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


17. The following criteria indicate successful completion of ECD-PKD checkout.


18. If these criteria are not met, repeat the test.

Table 9-3. ECD-PKD Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 40	< 400
	Wander (μV)	< 80	< 800
	Drift ($\mu\text{V}/\text{h}$)	< 200	< 2000
	Analytical Results		
	Lindane Signal-to-Noise Ratio		> 2 000
	Aldrin Signal-to-Noise Ratio		> 1 000
			

CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 40
	Wander (μV)	< 80
	Drift ($\mu\text{V}/\text{h}$)	< 200
	Analytical Results	
	Lindane Signal-to-Noise Ratio	> 2 000
	Aldrin Signal-to-Noise Ratio	> 1 000
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 400
Wander (μV)	< 800
Drift ($\mu\text{V/h}$)	< 2 000
Analytical Results	
Lindane Signal-to-Noise Ratio	> 2 000
Aldrin Signal-to-Noise Ratio	> 1 000
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 4 000
Wander (Counts)	< 8 000
Drift (Counts/h)	< 20 000
Analytical Results	
Lindane Signal-to-Noise Ratio	> 2 000
Aldrin Signal-to-Noise Ratio	> 1 000
	

Analytical Acceptance Comments	
1	Using Chrom-Card, set the signal-to-noise report parameters as described in the current procedure.
2	Using ChromQuest, Atlas, Xcalibur or a Computing integrator (e.g. ChromJet), calculate the S/N ratio as <i>Peak Height (counts)/noise (counts)</i> .

Checkout Using ECD with PPKD Injector

Chapter at a Glance...

SOP Number: P0311/07/E - 01 September 2009	132
Scope	132
Parts Referenced.....	132
Analytical Conditions Required for Purged Packed Injector	134
Recommended Initial Operations	135

Operating Procedures

ECD-PPKD Checkout	136
-------------------------	-----

SOP Number: P0311/07/E - 01 September 2009

Scope

Use the following procedure to verify proper ECD operation with Purged Packed Injector.

Parts Referenced

Table 10-1. ECD-PPKD Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	2 mm ID tapered	453 220 50
Liner Seal	Graphite seal for glass liner	290 334 05
Pre-column	Fused Silica Capillary Column 0.5 mt long 0.53 mm ID	260 603 75
Retaining Nut	M4 capillary column retaining nut	350 324 23
Press-fit connections	For columns 0.53/0.32 mm ID	350 438 16
Injection side adapter	For Wide bore column	347 003 03
Retaining Nut	For Injection side adapter	350 024 04
Detector side adapter	For Wide bore column	347 103 04
Ferrule	6 mm ID double brass ferrule	290 341 37
Nut	1/4" G-6 mm ID nut	350 201 18
Graphite Ferrule	Graphite ferrule for 0.53 mm ID Column Graphite ferrule for 0.32 mm ID Column	290 134 86 290 134 87
Septum	Standard septum for Purged Packed Injector	313 032 26
Syringe	10 µl size; 50 mm needle length	365 005 25

Table 10-1. ECD-PPKD Parts Referenced (Continued)

Part	Description	Part Number	
Test Mixture	Two components in Iso-octane	338 190 11	
	<i>Component</i> <i>Concentration</i>		
	Lindane 0.030 µg/ml		
Aldrin 0.030 µg/ml			
Gases	Chromatographic-grade purity		
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator		

Analytical Conditions Required for Purged Packed Injector

Table 10-2. ECD-PPKD Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Make-up Gas: Nitrogen = 30 ml/min
Oven Program	Initial Temperature = 70 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 220 °C Final Time = 1 minute
Injector	Operating Mode = Wide bore Temperature = 200 °C
Detector	Base Temperature = 250 °C ECD temperature = 300 °C Reference Current = 1 nA Pulse Amplitude = 50 V Pulse Width = 1µs
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Insert the glass liner.
Remove the column and insert the 2 mm ID tapered glass liner, as required for the checkout, from the bottom of the injector with the appropriate liner seal. Fix the liner by using the appropriate adapter for capillary column. Refer to the TRACE GC Maintenance and Troubleshooting Manual.
2. Mount the adapter for capillary column on the detector base body.
3. Replace the septum
A new septum should be installed properly in your injector.
4. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
5. Install the precolumn.
Connect the precolumn to the injector.
6. Install the test column
Connect the test column to the precolumn by using the press fit connections provided.
7. Perform the manual leak check following the Operating Procedure “*Performing a Leak Check*” in Chapter 14 of the TRACE GC Ultra Operating Manual.
8. Connect the other end of the test column to the detector base body.
9. Perform Column Evaluation.
10. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

ECD-PPKD Checkout

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	PPKD
Left carrier or Right carrier	He (helium)
Left detector or Right detector	ECD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select con pres. Scrool to Pressure and set the pressure value to have the required carrier gas flow rate Col. flow.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		220
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate PPKD Injector Control Table. Set the required temperature setpoint Temp. Verify to operate

in **Wide bore** mode. If not, scroll to **Mode**, press **MODE/TYPE** to access the selection menu, then select **Wide bore**.

LEFT INLET (PPKD) ¹		
Temp	200	200
Pressure	30.0	30.0
Mode:	Wide bore<	
Constant sept purge?	Y	<

1. These settings could also be for a right inlet.

- Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate ECD Detector Control Table. Set the required temperature **Base Temp** and the **Mkup** gas required setpoints.

LEFT DETECTOR (ECD) ¹		
Base temp	250	250
ECD Temp	300	300
Ref current nA	1.0	
Freq kHz	(2.20)	
Pulse amp V	50	
Pulse width μ s	1.0	
Mkup (N2)	30	30<

1. These settings could also be for a right detector.

- Set the **Reference Current** to 1.0 nA.
- Set the **Pulse Amplitude** to 50 V.
- Scroll to **Pulse Width** and press **ENTER** to open the menu selection. Select the pulse width to 1 μ s then press **ENTER**.
- Observe the ECD frequency value at the display. A base frequency value between 1 kHz and 3 kHz should be displayed.
If the ECD frequency value is less than 1kHz, reduce gradually the pulse amplitude till the base frequency value is about 1 kHz (the pulse amplitude value should not be less than 20V). Then, if necessary modify the reference current in order to have a base frequency of about 1-1,5 kHz. Let the detector signal to stabilize.

9. Activate your Data System and set the parameters required for the checkout.
10. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the ECD Detector Signal Control Table. Scroll to `Auto zero?` and turn it **YES**.

LEFT SIGNAL (ECD) ¹	
Output	(1000)
Offset	100<
Auto zero?	Y/N
Baseline comp	Off

1. These settings could also be for a right signal.

11. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 10-3 according to the data handling in use.

12. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
13. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
14. When the GC is ready, inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 10.1*.

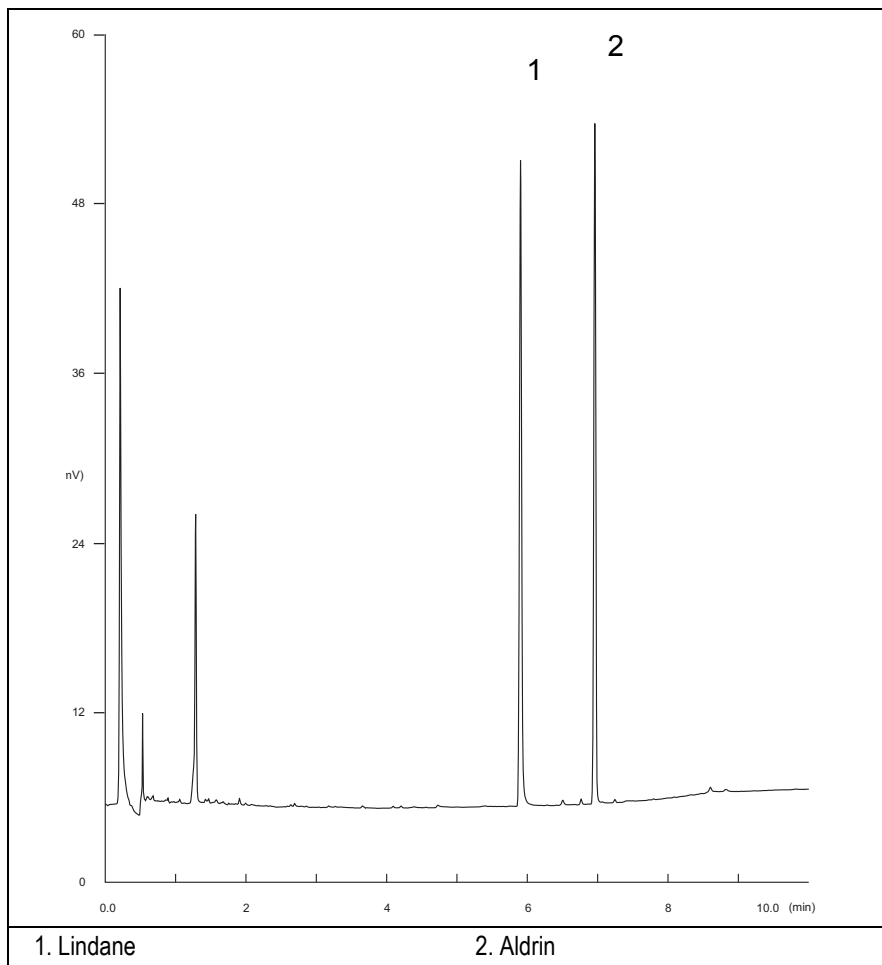


Figure 10-1. ECD-PPKD Injection

15. Establish the integration parameters and the peak table identifying the test mix components.
16. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Lindane component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Lindane.
- Repeat the procedure to calculate the signal-to-noise ratio also for Aldrin.
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Aldrin.



NOTE

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows:

Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:

- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.




CAUTION


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


17. The following criteria indicate successful completion of ECD-PPKD checkout.


18. If these criteria are not met, repeat the test.

Table 10-3. ECD-PPKD Acceptance Criteria

Acceptance Values	CHROM-CARD		
	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 40	< 400
	Wander (μV)	< 80	< 800
	Drift ($\mu\text{V}/\text{h}$)	< 200	< 2 000
	Analytical Results		
	Lindane Signal-to-Noise Ratio		> 3 000
	Aldrin Signal-to-Noise Ratio		> 3 000
			

Acceptance Values	CHROMQUEST	
	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 40
	Wander (μV)	< 80
	Drift ($\mu\text{V}/\text{h}$)	< 200
	Analytical Results	
	Lindane Signal-to-Noise Ratio	> 3 000
	Aldrin Signal-to-Noise Ratio	> 3 000
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 400
Wander (μV)	< 800
Drift ($\mu\text{V/h}$)	< 2 000
Analytical Results	
Lindane Signal-to-Noise Ratio	> 3 000
Aldrin Signal-to-Noise Ratio	> 3 000
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 4 000
Wander (Counts)	< 8 000
Drift (Counts/h)	< 20 000
Analytical Results	
Lindane Signal-to-Noise Ratio	> 3 000
Aldrin Signal-to-Noise Ratio	> 3 000
	

Analytical Acceptance Comments	
1	Using Chrom-Card, set the signal-to-noise report parameters as described in the current procedure.
2	Using ChromQuest, Atlas, Xcalibur or a Computing integrator (e.g. ChromJet), calculate the S/N ratio as <i>Peak Height (counts)/noise (counts)</i> .

Checkout Using ECD with PTV Injector

Chapter at a Glance...

SOP Number: P0312/07/E - 01 September 2009	144
Scope	144
Parts Referenced	144
Analytical Conditions Required for PTV Splitless Injection	145
Recommended Initial Operations	146

Operating Procedures

ECD-PTV Checkout in PTV Splitless Mode	147
--	-----

SOP Number: P0312/07/E - 01 September 2009

Scope

Use the following procedure to verify proper ECD operation with the Programmable Temperature Vaporizing Injector.

Parts Referenced

Table 11-1. ECD-PTV Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary ColumnTR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Liner	Silcosteel 2 mm ID (set of 2)	453 220 44
Liner Seal	Graphite seal for liner	290 034 17
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
Septum	Standard septum for PTV injector (set of 10)	313 132 25
Syringe	10 µl size; 50 mm needle length	365 005 25
Test Mixture	Two components in Iso-octane <i>Component</i> <i>Concentration</i> Lindane 0.030 µg/ml Aldrin 0.030 µg/ml	338 190 11
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for PTV Splitless Injection

Table 11-2 ECD-PTV Analytical Conditions

Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Make-up Gas: Nitrogen = 30 ml/min
Oven Program	Initial Temperature = 70 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 220 °C Final Time = 1 minute
Injector	Operating Mode = PTV splitless Splitless Time = 0.8 minutes Split Flow = 50 ml/min Constant Septum Purge = Yes Inject Temp = 50 °C Inject Time = 0.1 minute Transfer ramp = 10 °C/sec Transfer Temperature = 260 °C Transfer time = 1 minute
Detector	Base Temperature = 250 °C ECD temperature = 300 °C Reference Current = 1 nA Pulse Amplitude = 50 V Pulse Width = 1µs
Injected Volume	1 µl of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Replace the liner.
The liner currently installed in your injector should be carefully removed and replaced with the 2 mm ID Silcosteel liner, as required for the checkout, with the appropriate liner seal.
2. Replace the septum
A new septum should be installed properly in your injector.
3. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
4. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
5. Perform Column Evaluation and Leak Test.
6. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

ECD-PTV Checkout in PTV Splitless Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Right inlet	PTV
Right carrier	He (helium)
Right detector	ECD

1. Use **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to `Flow` mode, press **MODE/TYPE** to access the selection menu, then select `con pres`. Scroll to `Pressure` and set the pressure value to have the required carrier gas flow rate `Col.flow`.

RIGHT CARRIER ¹		
Pressure	30.0	30.0
Col.flow	3.00	
Lin. veloc.		(60.9) <

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		220
Final time 1		1.00 <
Ramp 2		Off

3. Use **RIGHT INLET** to display the appropriate Programmable Temperature Vaporizing Injector Control Table. Set the required temperature setpoint `Temp`. Verify to operate in **PTV splitless** mode. If not, scroll to `Mode`, press

MODE/TYPE to access the selection menu, then select PTV splitless. Scroll to Splitless time to set the required setpoint.

RIGHT INLET (PTV)		
Temp	70	70
Pressure	30.0	30.0
Mode:	PTV Splitless	
Total flow	(53.0)	
Split Flow	50.0	50.0
Splitless time	0.80	0.80
Constant sept purge?	Y	
Inject phase menu:	Y<	

4. Scroll to Inject phase menu. Press **MODE/TYPE** to enter the **PTV Phase Menu**.

PTV PHASE MENU	
Ramped pressure?	N
Inject temp	50
Inject time	0.1
Transfer ramp	10
Transfer temp	260
Transfer time	1.00<

5. Select Ramped pressure? **NO**.
Set the required Inject temp and Inject time setpoints as required. Then, set the Transfer ramp, the Transfer temp and the Transfer time required setpoints.
6. Use **RIGHT DETECTOR** to display the appropriate ECD Detector Control Table. Set the required temperature Base Temp and the Mkup gas required setpoints.

RIGHT DETECTOR (ECD)		
Base temp	250	250
ECD Temp	300	300
Ref current nA		1.0
Freq kHz		(2.20)
Pulse amp V		50
Pulse width μ s		1.0
Mkup (N2)	30	30<

7. Set the Reference Current to 1.0 nA.
8. Set the Pulse Amplitude to 50 V.
9. Scroll to Pulse Width and press **ENTER** to open the menu selection. Select the pulse width to 1 μ s then press **ENTER**.
10. Observe the ECD frequency value at the display. A base frequency value between 1 kHz and 3 kHz should be displayed.
If the ECD frequency value is less than 1kHz, reduce gradually the pulse amplitude till the base frequency value is about 1 kHz (the pulse amplitude value should not be less than 20V). Then, if necessary modify the reference current in order to have a base frequency of about 1-1,5 kHz. Let the detector signal to stabilize.
11. Activate your Data System and set the parameters required for the checkout.
12. Use **RIGHT SIGNAL** to display the ECD Detector Signal Control Table. Scroll to Auto zero? and turn it **YES**.

RIGHT SIGNAL (ECD)	
Output	(1000)
Offset	100<
Auto zero?	Y/N
Baseline com'	Off

13. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



Refer to the Acceptance Values reported in the Table 11-3 according to the data handling in use.

14. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
15. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
16. When the GC is ready, inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 11.1*.

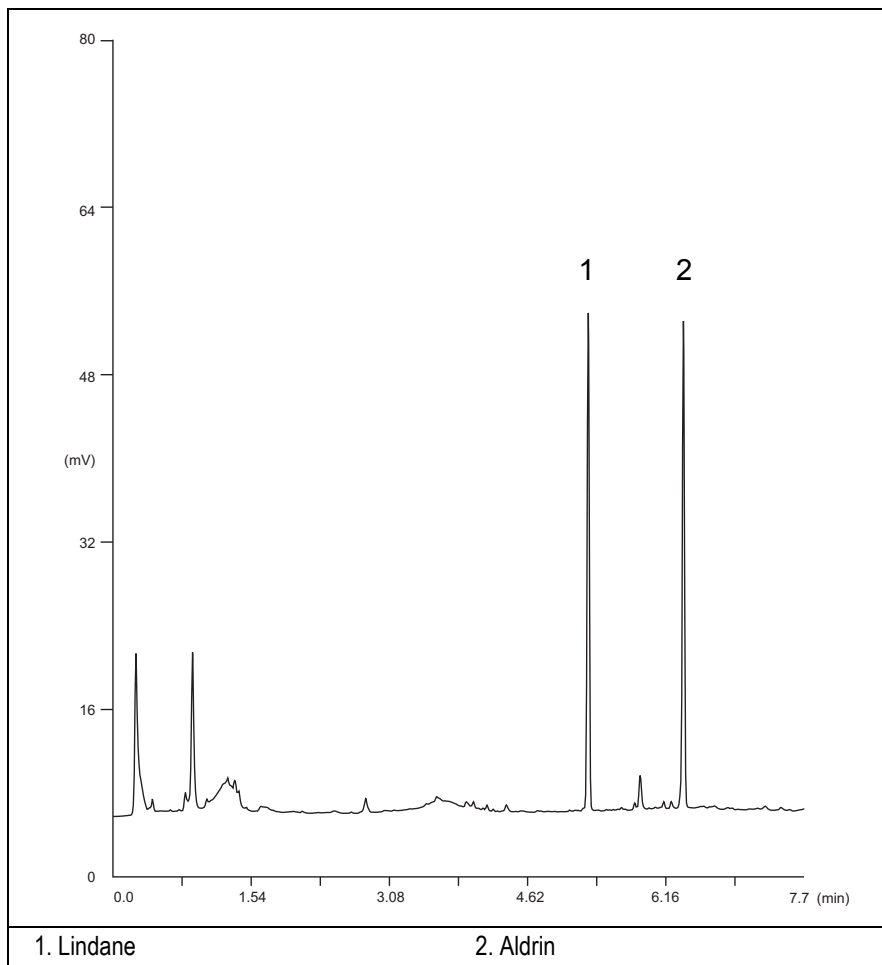


Figure 11-1. ECD-PTV Injection

17. Establish the integration parameters and the peak table identifying the test mix components.
18. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Lindane component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Lindane.
- Repeat the procedure to calculate the signal-to-noise ratio also for Aldrin.
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Aldrin.



NOTE

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows:

Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:

- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.




CAUTION


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


19. The following criteria indicate successful completion of ECD-PPKD checkout.


20. If these criteria are not met, repeat the test.

Table 11-3. ECD-PTV Acceptance Criteria

Acceptance Values	CHROM-CARD		
	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 40	< 400
	Wander (μV)	< 80	< 800
	Drift ($\mu\text{V}/\text{h}$)	< 200	< 2000
	Analytical Results		
	Lindane Signal-to-Noise Ratio		> 3 000
	Aldrin Signal-to-Noise Ratio		> 3 000
			

Acceptance Values	CHROMQUEST	
	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 40
	Wander (μV)	< 80
	Drift ($\mu\text{V}/\text{h}$)	< 200
	Analytical Results	
	Lindane Signal-to-Noise Ratio	> 3 000
	Aldrin Signal-to-Noise Ratio	> 3 000
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 400
Wander (μV)	< 800
Drift ($\mu\text{V/h}$)	< 2 000
Analytical Results	
Lindane Signal-to-Noise Ratio	> 3 000
Aldrin Signal-to-Noise Ratio	> 3 000
	

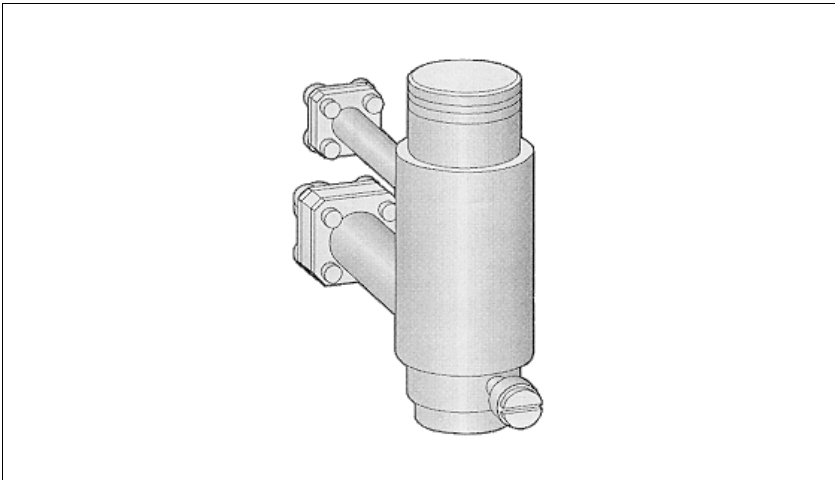
XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 4 000
Wander (Counts)	< 8 000
Drift (Counts/h)	< 20 000
Analytical Results	
Lindane Signal-to-Noise Ratio	> 3 000
Aldrin Signal-to-Noise Ratio	> 3 000
	

Analytical Acceptance Comments	
1	Using Chrom-Card, set the signal-to-noise report parameters as described in the current procedure.
2	Using ChromQuest, Atlas, Xcalibur or a Computing integrator (e.g. ChromJet), calculate the S/N ratio as <i>Peak Height (counts)/noise (counts)</i> .

SECTION

IV

SOPs Using NPD



The *SOPs Using NPD* section, contains the procedures to test the TRACE GC with the Nitrogen Phosphorus Detector (NPD) using different injectors.

Chapter 12, *Checkout Using NPD with S/SL Injector*.

Chapter 13, *Checkout Using NPD with OC Injector*.

Chapter 14, *Checkout Using NPD with PKD Injector*.

Chapter 15, *Checkout Using NPD with PPKD Injector*.

Chapter 16, *Checkout Using NPD with PTV Injector*.

Checkout Using NPD with S/SL Injector

Chapter at a Glance...

SOP Number: P0296/08/E - 01 September 2009	158
Scope	158
Parts Referenced	158
Analytical Conditions Required for Splitless Injection	159
Recommended Initial Operations	160

Operating Procedures

NPD-S/SL Checkout in Splitless Mode	161
---	-----

SOP Number: P0296/08/E - 01 September 2009

Scope

Use the following procedure to verify proper NPD operation with the Split/Splitless Injector.

Parts Referenced

Table 12-1. NPD-S/SL Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long; 0.32 mm ID; 0.25 µm film. thickness.	260 800 01
Glass Liner	3 mm ID for splitless injection	453 200 32
Liner Seal	Graphite seal for glass liner	290 334 06
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
Septum	Standard septum for S/SL injector	313 032 11
Syringe	10 µl size; 70 mm needle length	365 001 03
Test Mixture	Three components in Iso-Octane: <i>Component</i> <i>Concentration</i> Azobenzene 1 µg/ml Octadecane 1000 µg/ml Parathion methyl 1 µg/ml	338 190 06
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for Splitless Injection

Table 12-2. NPD-S/SL Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 2.3 ml/min Air = 60 ml/min Make-up: Nitrogen = 15 ml/min
Oven Program	Initial Temperature = 70 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 230 °C Final Time = 1 minute
Injector	Operating Mode = Splitless Temperature = 230 °C Splitless Time = 0.8 minutes Split Flow = 60 ml/min Constant Septum Purge = Yes
Detector	Base Temperature = 300 °C Source Current = Refer to <i>Source Ignition</i> Polarizer voltage = 3.5 V Detector Signal Range = 10 ⁰
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Replace the glass liner.
The glass liner currently installed in your injector should be carefully removed and replaced with the 3 mm ID glass liner for splitless application, as required for the checkout, with the appropriate graphite seal.
2. Replace the septum
A new septum should be installed properly in your injector.
3. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
4. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
5. Perform Column Evaluation and Leak Test
6. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

NPD-S/SL Checkout in Splitless Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	S/SL
Left carrier or Right carrier	He (helium)
Left detector or Right detector	NPD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select con pres. Scrool to Pressure and set the pressure value to have the required carrier gas flow rate Col.flow.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col.flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		230
Final time 1		1.00 <
Ramp 2		Off

- Use **LEFT INLET** or **RIGHT INLET** to display the appropriate Split/Splitless Injector Control Table and set the required temperature setpoint *Temp*. Verify to operate in **Splitless** mode. If not, scroll to Mode, press **MODE/TYPE** to access the selection menu, then select Splitless. Scroll to Splitless time to set the required setpoint.

LEFT INLET (S/SL) ¹		
Temp	230	230
Pressure	30.0	30.0
Mode:	Splitless	
Total flow	(63.0)	
Split Flow	60.0	60.0
Splitless time	0.80	0.80
Constant sept purge?	Y<	

1. These settings could also be for a right inlet.

- Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate NPD Detector Control Table. .

LEFT DETECTOR (NPD) ¹		
Source cur,A	OFF	
Base temp	300	300
Signal pA	(10.4)	
Target curr. pA	(X.XX)	
Auto adjust	No	
Polarizer V	3.5	
H2 delay time	Off	
H2	2.3	2.3
Air	60	60
Mkup N2	15	15

1. These settings could also be for a right detector.

- Scroll to Polarizer V and set 3.5 V.

Source Ignition

6. Switch on the source operating as follows:
 - a. Open the detector gases H₂, Air and M_{kup} and set the gas flow rates as follows:
 - H₂ = 2.3 ml/min
 - Air = 60 ml/min
 - M_{kup} N₂ = 15 ml/min)
 - b. Increase the Base Temp to 300°C and wait that the NPD cell reaches the correct set temperature.
 - c. Be sure that the backoff signal is between 0 and 0.5 pA.
 - d. Switch on the source with an initial current of 2.50 A. The backoff signal can slightly increase, but should remain within 0 and 1.5 pA.
 - e. Monitor the signal through the keypad or through the data system, increase the current value by steps of 0.002 A, until an immediate and strong increase of the signal is observed.
 - f. Wait five minutes to let the source stabilizes.
 - g. Check that source is correctly switched on decreasing hydrogen flow to 0.5 ml/min until signal decreases down to zero, then increase again to original value.
 - If the signal remains around zero, it means that the source is not switched on and it is necessary to increase further the current, accordingly to the procedure just described.
 - If the signal rises back to original value, it means that source is correctly switched on
 - h. Increase the current value of 2% of the actual ignition current. Let the signal stabilizes until its level drops below 20 pA.

7. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate NPD Detector Signal Control Table. Scroll to `Range` and set the electrometer amplifier input range required.

LEFT SIGNAL (NPD) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline comp	Off

1. These settings could also be for a right signal.

8. Activate your Data System and set the parameters required for the checkout.
9. In the NPD Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
10. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 12-3 according to the data handling in use.

11. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
12. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
13. Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 12.1*.



CAUTION

Verify the peak shape. If any peak distortion is visible, change the analytical test column.

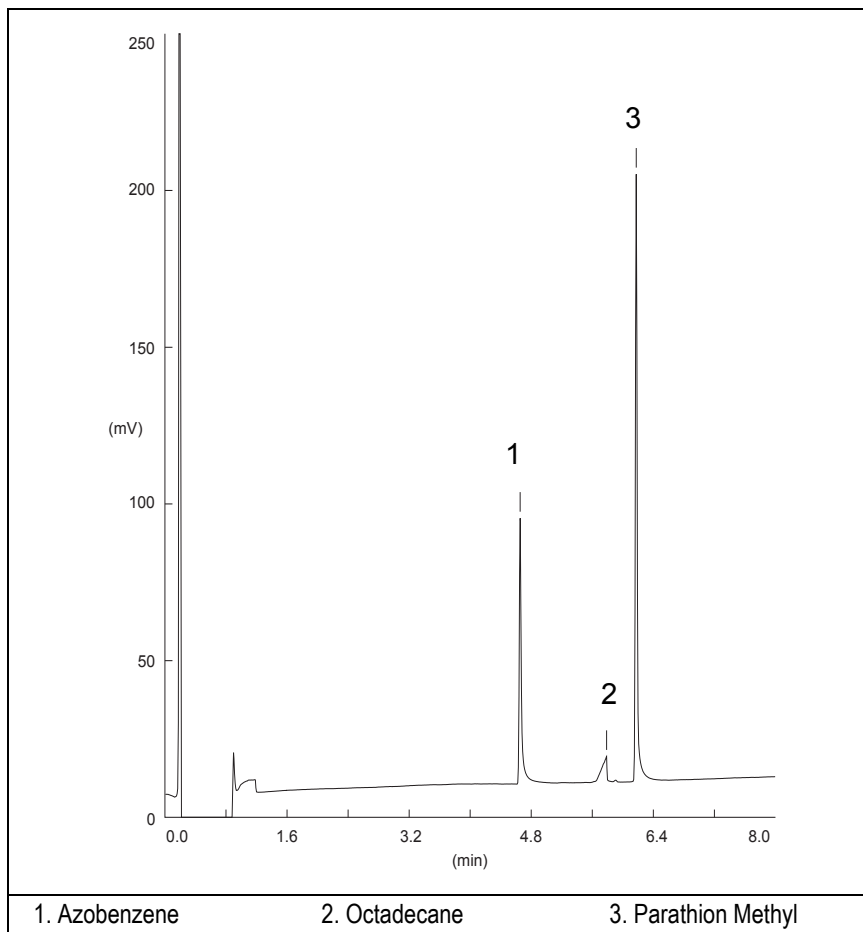


Figure 12-1. NPD-Splitless Injection

14. Establish the integration parameters and the table identifying the test mix components.
15. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Azobenzene component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Azobenzene.
- Repeat the procedure to calculate the signal-to-noise ratio also for Parathion Methyl.
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Parathion Methyl.



NOTE

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows:

Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:

- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.





CAUTION


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


16. The following criteria indicate successful completion of NPD-S/SL checkout.
17. If these criteria are not met, repeat the test.


Table 12-3. NPD-S/SL Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 40	< 400
	Wander (μV)	< 100	< 1 000
	Drift ($\mu\text{V/h}$)	< 300	<3 000
	Analytical Results		
	Azobenzene Signal-to-Noise Ratio	> 800	
	Parathion Methyl Signal-to-Noise Ratio	> 2 500	
	Octadecane Signal-to-Noise Ratio	Negligible	
			

Computing-integrator (e.g. ChromJet)		
		

CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 40
	Wander (μV)	< 100
	Drift ($\mu\text{V/h}$)	< 300
	Analytical Results	
	Azobenzene Signal-to-Noise Ratio	> 800
	Parathion Methyl Signal-to-Noise Ratio	> 2 500
	Octadecano Signal-to-Noise Ratio	Negligible
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 400
Wander (μV)	< 1 000
Drift ($\mu\text{V/h}$)	< 3 000
Analytical Results	
Azobenzene Signal-to-Noise Ratio	> 800
Parathion Methyl Signal-to-Noise Ratio	> 2 500
Octadecano Signal-to-Noise Ratio	Negligible
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 4 000
Wander Counts)	< 10 000
Drift (Counts/h)	< 30 000
Analytical Results	
Azobenzene Signal-to-Noise Ratio	> 800
Parathion Methyl Signal-to-Noise Ratio	> 2 500
Octadecano Signal-to-Noise Ratio	Negligible
	

Checkout Using NPD with OC Injector

Chapter at a Glance...

SOP Number: P0297/08/E - 01 September 2009	170
Scope	170
Parts Referenced	170
Analytical Conditions Required for On-Column Injection	171
Recommended Initial Operations	172

Operating Procedures

NPD-OCI Checkout in On-Column Mode	173
--	-----

SOP Number: P0297/08/E - 01 September 2009

Scope

Use the following procedure to verify proper NPD operation with the On-Column Injector.

Parts Referenced

Table 13-1. NPD-OCI Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long; 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Graphite Ferrule	Graphite Ferrule for 0.32 mm ID Column	290 134 87
Vespel Ferrule	Vespel Ferrule for 0.32 mm ID Column	290 134 60
Retaining Nut	M4 capillary column retaining nut	350 324 23
Syringe	10 µl size; 75 mm needle length	365 020 07
Test Mixture	Three components in Iso-Octane: <i>Component</i> <i>Concentration</i> Azobenzene 1 µg/ml Octadecane 1000 µg/ml Parathion methyl 1 µg/ml	338 190 06
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	
Syringe	10 µl size; 80 mm needle length	365 020 19
Pre-column	2 m long; 0.53 mm ID	260 603 75
Press-fit set	Set of Press-fit connectors for TRACE OC	350 038 45

Analytical Conditions Required for On-Column Injection

Table 13-2. NPD-OCI Analytical Conditions

Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 2.3 ml/min Air = 60 ml/min Make-up: Nitrogen = 15 ml/min
Oven Program	Initial Temperature = 85 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 230 °C Final Time = 1 minute
Injector	Secondary cooling = 0.2 minutes
Detector	Base Temperature = 300 °C Source Current = Refer to <i>Source Ignition</i> Polarizer voltage = 3.5 V Detector Signal Range = 10 ⁰
Injected Volume	1 µl of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
2. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
In case of automatic On-column for TriPlus sampler, install the pre-column and connect it to the test column by press-fit connector.
3. Install and connect the TriPlus sampler and its components.
4. Perform Column Evaluation and Leak Test.
5. Connect your data handling.
Verify that your data handling is properly connected to your GC system.
6. Verify the opening/closing of the OC injector actuator by using the proper commands.
7. Verify the alignment of the syringe on the OC injector.

OPERATING PROCEDURE

NPD-OCI Checkout in On-Column Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	OCI
Left carrier or Right carrier	He (helium)
Left detector or Right detector	NPD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select con pres. Scrool to Pressure and set the pressure value to have the required carrier gas flow rate Col. flow.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	85.0	85.0
Initial Time		1.00
Ramp 1		20.0
Final temp		230
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate On-Column Injector Control Table. Scroll to *Sec. cool time* and set the required secondary cooling time.

LEFT INLET (OCI) ¹		
Pressure	30.0	30.0
Sec. Cool Time		0.2<

1. These settings could also be for a right inlet.

4. Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate NPD Detector Control Table. Set the required temperature *Base Temp* and the detector gases *H2, Air* and *Mkup* required setpoints.

LEFT DETECTOR (NPD) ¹		
Source cur, A		Off
Base temp	300	300
Signal pA		(10.4)
Target curr. pA		(X.XX)
Auto adjust		No
Polarizer V		3.5
H2 delay time		Off
H2	2.3	2.3
Air	60	60
Mkup N2	15	15

1. These settings could also be for a right detector.

5. Scroll to *Polarizer V* and set 3.5 V.

Source Ignition

6. Switch on the source operating as follows:
 - a. Open the detector gases *H2, Air* and *Mkup* and set the gas flow rates as follows:

- $H_2 = 2.3$ ml/min
 - $Air = 60$ ml/min
 - $Mkup N_2 = 15$ ml/min
- b. Increase the `Base Temp` to 300°C and wait that the NPD cell reaches the correct set temperature.
 - c. Be sure that the backoff signal is between 0 and 0.5 pA.
 - d. Switch on the source with an initial current of 2.50 A. The backoff signal can slightly increase, but should remain within 0 and 1.5 pA.
 - e. Monitor the signal through the keypad or through the data system, increase the current value by steps of 0.002 A, until an immediate and strong increase of the signal is observed.
 - f. Wait five minutes to let the source stabilizes.
 - g. Check that source is correctly switched on decreasing hydrogen flow to 0.5 ml/min until signal decreases down to zero, then increase again to original value.
 - If the signal remains around zero, it means that the source is not switched on and it is necessary to increase further the current, accordingly to the procedure just described.
 - If the signal rises back to original value, it means that source is correctly switched on
 - h. Increase the current value of 2% of the actual ignition current. Let the signal stabilizes until its level drops below 20 pA.
7. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate NPD Detector Signal Control Table. Scroll to `Range` and set the electrometer amplifier input range required.

LEFT SIGNAL (NPD) ¹	
Output	(1000)
Offset	100
Autozero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline comp	Off

1. These settings could also be for a right signal.

8. Activate your Data System and set the parameters required for the checkout.
9. In the NPD Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
10. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 13-3 according to the data handling in use.

11. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
12. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
13. Perform the analysis.

Manual injection

- Inject the test mixture and press **START** on the GC to begin the checkout run.

Automatic injection with TriPlus sampler

- Fill a vial with the standard mix and place that vial in the sample tray.

- Load the method for OC and perform the sampling.

The resulting chromatogram should look like the one shown in *Figure 7.1*.



CAUTION Verify the peak shape. If any peak distortion is visible, change the analytical test column.

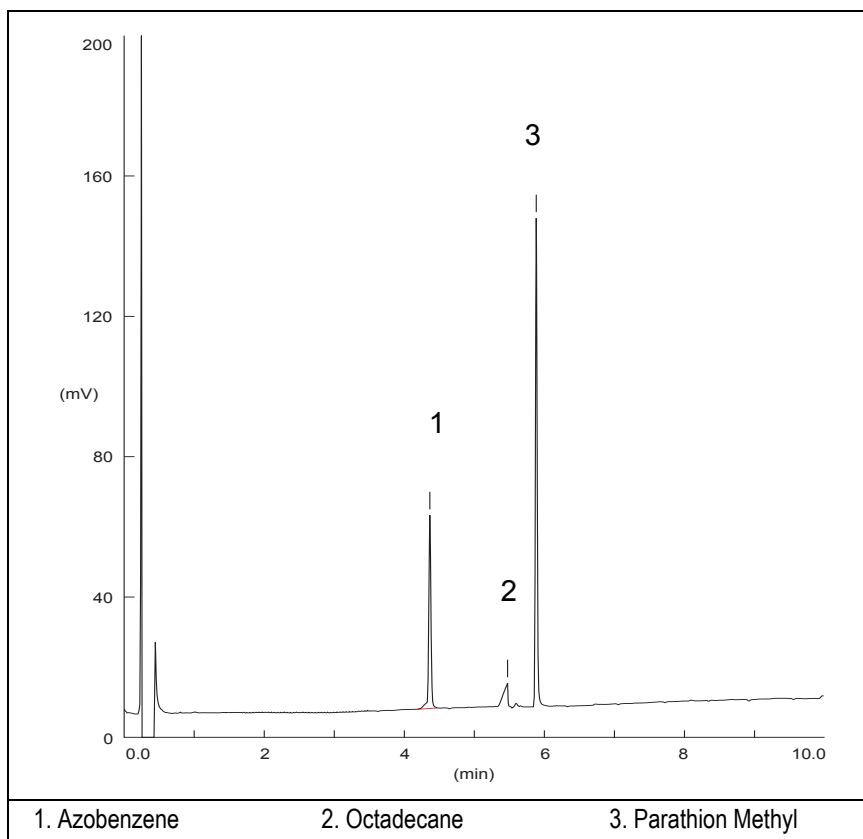


Figure 13-1. NPD-On-Column Injection

14. Establish the integration parameters and the peak table identifying the test mix components.
15. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Azobenzene component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Azobenzene.
- Repeat the procedure to calculate the signal-to-noise ratio also for Parathion Methyl.
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Parathion Methyl.



NOTE

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows:

Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:

- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.





CAUTION


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


16. The following criteria indicate successful completion of NPD-OCI checkout. If these criteria are not met, repeat the test.


Table 13-3. NPD-OCI Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 40	< 400
	Wander (μV)	< 100	< 1 000
	Drift ($\mu\text{V}/\text{h}$)	< 300	<3 000
	Analytical Results		
	Azobenzene Signal-to-noise ratio	> 550	
	Parathion Methyl Signal-to-noise ratio	> 1 500	
	Octadecane Signal-to-noise ratio	Negligible	
			

Computing-integrator (e.g. ChromJet)	
	

CHROMQUEST			
Acceptance Values	Baseline Parameters (1V Full Scale)		
	Noise (μV)	< 40	
	Wander (μV)	< 100	
	Drift ($\mu\text{V}/\text{h}$)	< 300	
	Analytical Results		
	Azobenzene Signal-to-noise ratio	> 550	
	Parathion Methyl Signal-to-noise ratio	> 1 500	
	Octadecane Signal-to-noise ratio	Negligible	
			

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 400
Wander (μV)	< 1 000
Drift ($\mu\text{V/h}$)	< 3 000
Analytical Results	
Azobenzene Signal-to-noise ratio	> 550
Parathion Methyl Signal-to-noise ratio	> 1 500
Octadecane Signal-to-noise ratio	Negligible
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 4 000
Wander Counts)	< 10 000
Drift (Counts/h)	< 30 000
Analytical Results	
Azobenzene	> 550
Parathion Methyl	> 1 500
Octadecane	Negligible
	

Checkout Using NPD with PKD Injector

Chapter at a Glance...

SOP Number: P0313/08/E - 01 September 2009	182
Scope	182
Parts Referenced	182
Analytical Conditions Required for Packed Injector.....	184
Recommended Initial Operations	185

Operating Procedures

NPD-PKD Checkout	186
------------------------	-----

SOP Number: P0313/08/E - 01 September 2009

Scope

Use the following procedure to verify proper NPD operation with Packed Injector.

Parts Referenced

Table 14-1. NPD-PKD Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	2 mm ID tapered	453 220 50
Liner Seal	Graphite seal for glass liner	290 334 05
Pre-column	Fused Silica Capillary Column 0.5 mt long 0.53 mm ID	260 603 75
Retaining Nut	M4 capillary column retaining nut	350 324 23
Press-fit connections	For columns 0.53/0.32 mm ID	350 438 16
Injection side adapter	For Wide bore column	347 003 03
Retaining Nut	For Injection side adapter	350 024 04
Detector side adapter	For Wide bore column	347 103 04
Ferrule	6 mm ID double brass ferrule	290 341 37
Nut	1/4" G-6 mm ID nut	350 201 18
Graphite Ferrule	Graphite ferrule for 0.53 mm ID Column Graphite ferrule for 0.32 mm ID Column	290 134 86 290 134 87
Septum	Standard septum for Packed Injector	313 032 26
Syringe	10 µl size; 50 mm needle length	365 005 25

Table 14-1. NPD-PKD Parts Referenced (Continued)

Part	Description	Part Number
Test Mixture	Three components in Iso-Octane: <i>Component</i> <i>Concentration</i> Azobenzene 1 µg/ml Octadecane 1000 µg/ml Parathion methyl 1 µg/ml	338 190 06
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for Packed Injector

Table 14-2. NPD-PKD Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 2.3 ml/min Air = 60 ml/min Make-up Gas: Nitrogen = 15 ml/min
Oven Program	IsoTemperature = 50 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 200 °C Final Time = 1 minute
Injector	Operating Mode = Packed Temperature = 200 °C
Detector	Base Temperature = 300 °C Source Current = Refer to <i>Source Ignition</i> Polarizer voltage = 3.5 V Detector Signal Range = 10 ⁰
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Insert the glass liner.
Remove the column and insert the 2 mm ID tapered glass liner, as required for the checkout, from the bottom of the injector with the appropriate liner seal. Fix the liner by using the appropriate adapter for capillary column. Refer to the TRACE GC Ultra Maintenance and Troubleshooting Manual.
2. Mount the adapter for capillary column on the detector base body.
3. Replace the septum
A new septum should be installed properly in your injector.
4. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
5. Install the precolumn.
Connect the precolumn to the injector.
6. Install the test column
Connect the test column to the precolumn by using the press fit connections provided.
7. Perform the manual leak check following the Operating Procedure “*Performing a Leak Check*” in Chapter 14 of the TRACE GC Ultra Operating Manual.
8. Connect the other end of the test column to the detector base body.
9. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

NPD-PKD Checkout

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	PKD
Left carrier or Right carrier	He (helium)
Left detector or Right detector	NPD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select con pres. Scrool to Pressure and set the pressure value to have the required carrier gas flow rate Col.flow.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col.flow	3.00	
Lin. veloc.		(60.9)<

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	50.0	50.0
Initial Time		1.00
Ramp 1		20.0
Final temp		200
Final time 1		1.00<
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate PKD Injector Control Table. Set the required temperature setpoint Temp. Verify to operate

in **Packed** mode. If not, scroll to Mode, press **MODE/TYPE** to access the selection menu, then select Packed.

LEFT INLET (PKD) ¹		
Temp	200	200
Pressure	30.0	30.0
Mode:		Packed

1. These settings could also be for a right inlet.

4. Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate NPD Detector Control Table. Set the required temperature Base Temp and the detector gases H₂, Air and Mkup required setpoints.

LEFT DETECTOR (NPD) ¹		
Source cur, A		Off
Base temp	300	300
Signal pA		(10.4)
Target curr. pA		(X.XX)
Auto adjust		No
Polarizer V		3.5
H ₂ delay time		Off
H ₂	2.3	2.3
Air	60	60
Mkup N ₂	15	15

1. These settings could also be for a right detector.

5. Scroll to Polarizer V and set 3.5 V.

Source Ignition

6. Switch on the source operating as follows:
 - a. Open the detector gases H₂, Air and Mkup and set the gas flow rates as follows:

- H₂ = 2.3 ml/min
 - Air = 60 ml/min
 - Makeup N₂ = 15 ml/min)
- b. Increase the Base Temp to 300°C and wait that the NPD cell reaches the correct set temperature.
 - c. Be sure that the backoff signal is between 0 and 0.5 pA.
 - d. Switch on the source with an initial current of 2.50 A. The backoff signal can slightly increase, but should remain within 0 and 1.5 pA.
 - e. Monitor the signal through the keypad or through the data system, increase the current value by steps of 0.002 A, until an immediate and strong increase of the signal is observed.
 - f. Wait five minutes to let the source stabilizes.
 - g. Check that source is correctly switched on decreasing hydrogen flow to 0.5 ml/min until signal decreases down to zero, then increase again to original value.
 - If the signal remains around zero, it means that the source is not switched on and it is necessary to increase further the current, accordingly to the procedure just described.
 - If the signal rises back to original value, it means that source is correctly switched on
 - h. Increase the current value of 2% of the actual ignition current. Let the signal stabilizes until its level drops below 20 pA.
7. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate NPD Detector Signal Control Table. Scroll to Range and set the electrometer amplifier input range required.

LEFT SIGNAL (NPD) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline comp	Off

1. These settings could also be for a right signal.

8. Activate your Data System and set the parameters required for the checkout.
9. In the NPD Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
10. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 14-3 according to the data handling in use.

11. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
12. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
13. Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 14.1*.



CAUTION

Verify the peak shape. If any peak distortion is visible, change the analytical test column.

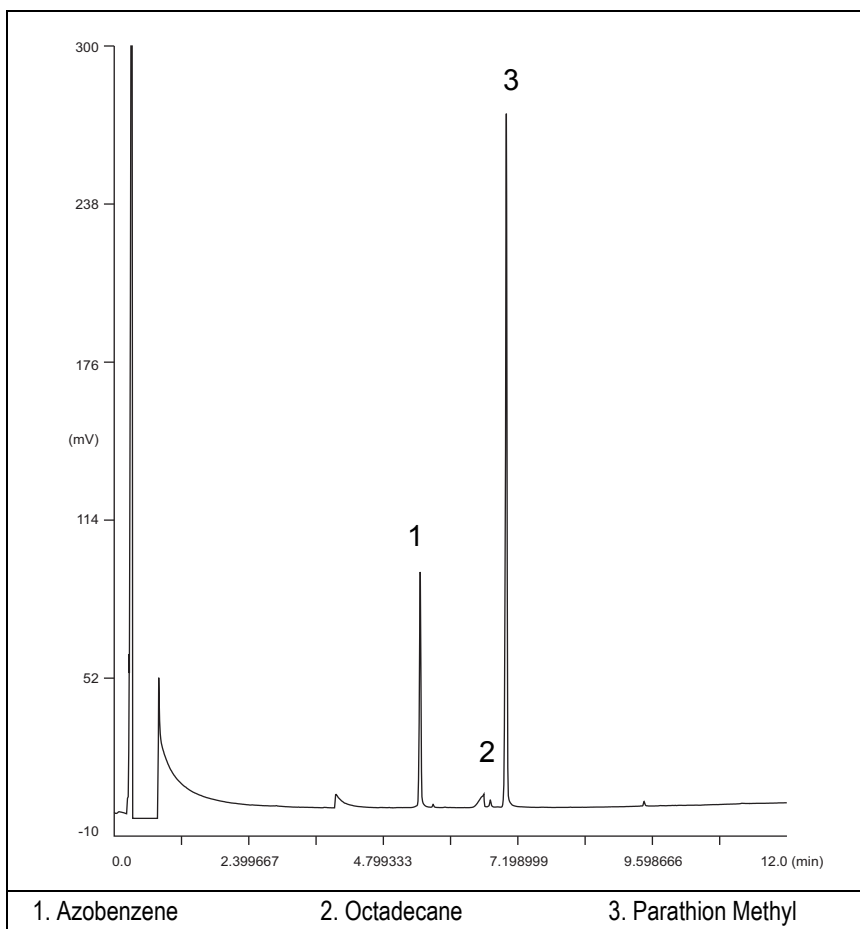


Figure 14-1. NPD-Packed Injection

14. Establish the integration parameters and the peak table identifying the test mix components.
15. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Azobenzene component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Azobenzene.
- Repeat the procedure to calculate the signal-to-noise ratio also for Parathion Methyl.
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Parathion Methyl.

**NOTE**

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows:

Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:


- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.


**CAUTION**


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


16. The following criteria indicate successful completion of NPD-PKD checkout.
17. If these criteria are not met, repeat the test.


Table 14-3. NPD-PKD Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 40	< 400
	Wander (μV)	< 100	< 1 000
	Drift ($\mu\text{V/h}$)	< 300	<3 000
	Analytical Results		
	Azobenzene Signal-to-noise ratio		600
	Parathion Methyl Signal-to-noise ratio		1 700
	Octadecano Signal-to-noise ratio		Negligible
			

Computing-integrator (e.g. ChromJet)	
	

CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 40
	Wander (μV)	< 100
	Drift ($\mu\text{V/h}$)	< 300
	Analytical Results	
	Azobenzene Signal-to-noise ratio	> 600
	Parathion Methyl Signal-to-noise ratio	> 1700
	Octadecane Signal-to-noise ratio	Negligible
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 400
Wander (μV)	< 1 000
Drift ($\mu\text{V/h}$)	< 3 000
Analytical Results	
Azobenzene Signal-to-noise ratio	> 600
Parathion Methyl Signal-to-noise ratio	> 1 700
Octadecane Signal-to-noise ratio	Negligible
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 4 000
Wander Counts)	< 10 000
Drift (Counts/h)	< 30 000
Analytical Results	
Azobenzene Signal-to-noise ratio	> 600
Parathion Methyl Signal-to-noise ratio	> 1 700
Octadecane Signal-to-noise ratio	Negligible
	

Checkout Using NPD with PPKD Injector

SOP Number: P0314/08/E - 01 September	196
Scope	196
Parts Referenced	196
Recommended Initial Operations	199
<i>Operating Procedures</i>	
NPD-PPKD Checkout	200

SOP Number: P0314/08/E - 01 September

Scope

Use the following procedure to verify proper NPD operation with Purged Packed Injector.

Parts Referenced

Table 15-1. NPD-PPKD Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	2 mm ID tapered	453 220 50
Liner Seal	Graphite seal for glass liner	290 334 05
Pre-column	Fused Silica Capillary Column 0.5 mt long 0.53 mm ID	260 603 75
Retaining Nut	M4 capillary column retaining nut	350 324 23
Press-fit connections	For columns 0.53/0.32 mm ID	350 438 16
Injection side adapter	For Wide bore column	347 003 03
Retaining Nut	For Injection side adapter	350 024 04
Detector side adapter	For Wide bore column	347 103 04
Ferrule	6 mm ID double brass ferrule	290 341 37
Nut	1/4" G-6 mm ID nut	350 201 18
Graphite Ferrule	Graphite ferrule for 0.53 mm ID Column Graphite ferrule for 0.32 mm ID Column	290 134 86 290 134 87
Septum	Standard septum for Purged Packed Injector	313 032 26
Syringe	10 µl size; 50 mm needle length	365 005 25

Table 15-1. NPD-PPKD Parts Referenced (Continued)

Part	Description	Part Number
Test Mixture	Three components in Iso-Octane: <i>Component</i> <i>Concentration</i> Azobenzene 1 µg/ml Octadecane 1000 µg/ml Parathion methyl 1 µg/ml	338 190 06
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for Purged Packed Injector

Table 15-2. NPD-PPKD Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 2.3 ml/min Air = 60 ml/min Make-up Gas: Nitrogen = 15 ml/min
Oven Program	Initial Temperature = 50 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 200 °C Final Time = 1 minute
Injector	Operating Mode = Wide bore Temperature = 200 °C
Detector	Base Temperature = 300 °C Source Current = Refer to <i>Source Ignition</i> Polarizer voltage = 3.5 V Detector Signal Range = 10 ⁰
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Insert the glass liner.
Remove the column and insert the 2 mm ID tapered glass liner, as required for the checkout, from the bottom of the injector with the appropriate liner seal. Fix the liner by using the appropriate adapter for capillary column. Refer to the TRACE GC Ultra Maintenance and Troubleshooting Manual.
2. Mount the adapter for capillary column on the detector base body.
3. Replace the septum
A new septum should be installed properly in your injector.
4. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
5. Install the precolumn.
Connect the precolumn to the injector.
6. Install the test column
Connect the test column to the precolumn by using the press fit connections provided.
7. Perform the manual leak check following the Operating Procedure “*Performing a Leak Check*” in Chapter 14 of the TRACE GC Ultra Operating Manual.
8. Connect the other end of the test column to the detector base body.
9. Perform Column Evaluation.
10. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

NPD-PPKD Checkout

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	PPKD
Left carrier or Right carrier	He (helium)
Left detector or Right detector	NPD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to Pressure and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	50.0	50.0
Initial Time		1.00
Ramp 1		20.0
Final temp		200
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate PPKD Injector Control Table. Set the required temperature setpoint **Temp**. Verify to operate in

Wide bore mode. If not, scroll to *Mode*, press **MODE/TYPE** to access the selection menu, then select *Widebore*.

LEFT INLET (PPKD) ¹		
Temp	200	200
Pressure	30.0	30.0
Mode:	Wide bore<	
Constant sept purge?		Y<

1. These settings could also be for a right inlet.

4. Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate NPD Detector Control Table. Set the required temperature *Base Temp* and the detector gases *H2*, *Air* and *Mkup* required setpoints.

LEFT DETECTOR (NPD) ¹		
Source cur, A		Off
Base temp	300	300
Signal pA		(10.4)
Target curr. pA		(X.XX)
Auto adjust		No
Polarizer V		3.5
H2 delay time		Off
H2	2.3	2.3
Air	60	60
Mkup N2	15	15

1. These settings could also be for a right detector.

5. Scroll to *Polarizer V* and set 3.5 V.

Source Ignition

6. Switch on the source operating as follows:
 - a. Open the detector gases *H2*, *Air* and *Mkup* and set the gas flow rates as follows:

- H₂ = 2.3 ml/min
 - Air = 60 ml/min
 - Makeup N₂ = 15 ml/min)
- b. Increase the Base Temp to 300°C and wait that the NPD cell reaches the correct set temperature.
 - c. Be sure that the backoff signal is between 0 and 0.5 pA.
 - d. Switch on the source with an initial current of 2.50 A. The backoff signal can slightly increase, but should remain within 0 and 1.5 pA.
 - e. Monitor the signal through the keypad or through the data system, increase the current value by steps of 0.002 A, until an immediate and strong increase of the signal is observed.
 - f. Wait five minutes to let the source stabilizes.
 - g. Check that source is correctly switched on decreasing hydrogen flow to 0.5 ml/min until signal decreases down to zero, then increase again to original value.
 - If the signal remains around zero, it means that the source is not switched on and it is necessary to increase further the current, accordingly to the procedure just described.
 - If the signal rises back to original value, it means that source is correctly switched on
 - h. Increase the current value of 2% of the actual ignition current. Let the signal stabilizes until its level drops below 20 pA.
7. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate NPD Detector Signal Control Table. Scroll to Range and set the electrometer amplifier input range required.

LEFT SIGNAL (NPD) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline comp	Off

1. These settings could also be for a right signal.

8. Activate your Data System and set the parameters required for the checkout.
9. In the NPD Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
10. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 15-3 according to the data handling in use.

11. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
12. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
13. Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 15.1*.



CAUTION

Verify the peak shape. If any peak distortion is visible, change the analytical test column.

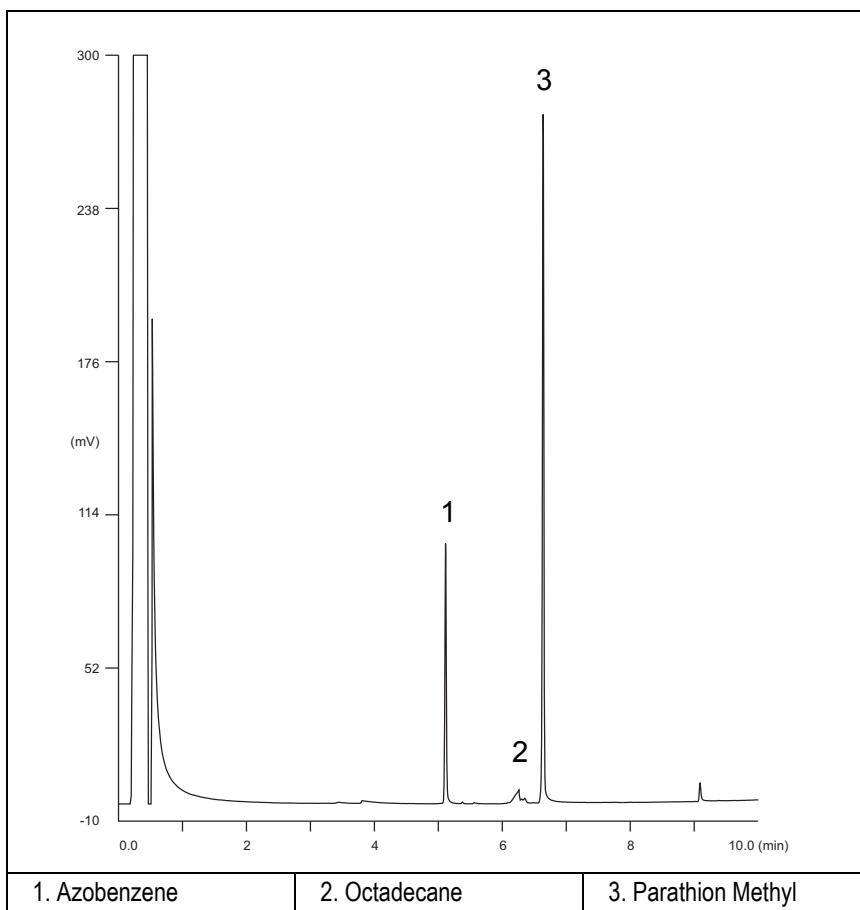


Figure 15-1. NPD-PPKD Injection

14. Establish the integration parameters and the peak table identifying the test mix components.
15. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Azobenzene component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Azobenzene.
- Repeat the procedure to calculate the signal-to-noise ratio also for Parathion Methyl.
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Parathion Methyl.

**NOTE**

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows: Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:


- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.


**CAUTION**


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


16. The following criteria indicate successful completion of NPD-PPKD checkout.
17. If these criteria are not met, repeat the test.


Table 15-3. NPD-PPKD Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 40	< 400
	Wander (μV)	< 100	< 1 000
	Drift ($\mu\text{V}/\text{h}$)	< 300	<3 000
	Analytical Results		
	Azobenzene Signal-to-noise ratio		600
	Parathion Methyl Signal-to-noise ratio		1 700
	Octadecane Signal-to-noise ratio		Negligible
			

Computing-integrator (e.g. ChromJet)	
	

CHROMQUEST			
Acceptance Values	Baseline Parameters (1V Full Scale)		
	Noise (μV)	< 40	
	Wander (μV)	< 100	
	Drift ($\mu\text{V}/\text{h}$)	< 300	
	Analytical Results		
	Azobenzene Signal-to-noise ratio		600
	Parathion Methyl Signal-to-noise ratio		1 700
	Octadecane Signal-to-noise ratio		Negligible
			

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 400
Wander (μV)	< 1 000
Drift ($\mu\text{V/h}$)	< 3 000
Analytical Results	
Azobenzene Signal-to-noise ratio	> 600
Parathion Methyl Signal-to-noise ratio	> 1 700
Octadecane Signal-to-noise ratio	Negligible
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 4 000
Wander Counts)	< 10 000
Drift (Counts/h)	< 30 000
Analytical Results	
Azobenzene Signal-to-noise ratio	> 600
Parathion Methyl Signal-to-noise ratio	> 1 700
Octadecane Signal-to-noise ratio	Negligible
	

Checkout Using NPD with PTV Injector

Chapter at a Glance...

SOP Number: P0315/08/E - 01 September 2009	210
Scope	210
Parts Referenced	210
Analytical Conditions Required for PTV Splitless Injection	211
Recommended Initial Operations	212

Operating Procedures

NPD-PTV Checkout in PTV Splitless Mode	213
--	-----

SOP Number: P0315/08/E - 01 September 2009

Scope

Use the following procedure to verify proper NPD operation with the Programmable Temperature Vaporizing Injector.

Parts Referenced

Table 16-1. NPD-PTV Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Liner	Silcosteel 2 mm ID (set of 2)	453 220 44
Liner Seal	Graphite seal for liner	290 034 17
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
Septum	Standard septum for PTV injector (set of 10)	313 132 25
Syringe	10 µl size; 50 mm needle length	365 005 25
Test Mixture	Three components in Iso-Octane: <i>Component</i> <i>Concentration</i> Azobenzene 1 µg/ml Octadecane 1000 µg/ml Parathion methyl 1 µg/ml	338 190 06
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for PTV Splitless Injection

Table 16-2. NPD-PTV Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 2.3 ml/min Air = 60 ml/min Make-up Gas: Nitrogen = 30 ml/min
Oven Program	Initial Temperature = 70 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 230 °C Final Time = 1 minute
Injector	Operating Mode = PTV Splitless Splitless Time = 0.8 minutes Split Flow = 50 ml/min Constant Septum Purge = Yes Inject Temp = 50 °C Inject Time = 0.1 minute Transfer ramp = 10 °C/sec Transfer Temperature = 260 °C Transfer time = 1 minutes
Detector	Base Temperature = 300 °C Source Current = Refer to <i>Source Ignition</i> Polarizer voltage = 3.5 V Detector Signal Range = 10 ⁰
Injected Volume	1 µl of Test Mixture

Table 16-2. NPD-PTV Analytical Conditions (Continued)

Parameters Setting	
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Replace the liner.
The liner currently installed in your injector should be carefully removed and replaced with the 2 mm ID Silcosteel glass liner, as required for the checkout, with the appropriate liner seal.
2. Replace the septum
A new septum should be installed properly in your injector.
3. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
4. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
5. Perform Column Evaluation and Leak Test.
6. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

NPD-PTV Checkout in PTV Splitless Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Right inlet	PTV
Right carrier	He (helium)
Right detector	NPD

1. Use **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to `Flow` mode, press **MODE/TYPE** to access the selection menu, then select `con pres`. Scroll to `Pressure` and set the pressure value to have the required carrier gas flow rate `Col.flow`.

RIGHT CARRIER ¹		
Pressure	30.0	30.0
Col.flow	3.00	
Lin. veloc.		(60.9)<

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		230
Final time 1		1.00<
Ramp 2		Off

3. Use **RIGHT INLET** to display the appropriate Programmable Temperature Vaporizing Injector Control Table. Set the required temperature setpoint

Temp. Verify to operate in **PTV splitless** mode. If not, scroll to Mode, press **MODE/TYPE** to access the selection menu, then select PTV splitless. Scroll to Splitless time to set the required setpoint.

RIGHT INLET (PTV)		
Temp	70	70
Pressure	30.0	30.0
Mode:	PTV Splitless	
Total flow	(53.0)	
Split Flow	50.0	50.0
Splitless time	0.80	0.80
Constant sept purge?		Y
Inject phase menu:		Y<

4. Scroll to Inject phase menu. Press **MODE/TYPE** to enter the **PTV Phase Menu**.

PTV PHASE MENU	
Ramped pressure?	N
Inject temp	50
Inject time	0.1
Transfer ramp	10
Transfer temp	260
Transfer time	1.00<

5. Select Ramped pressure? **NO**. Set the required Inject temp and *Inject time* setpoints as required. Then, set the Transfer ramp, the Transfer temp and the Transfer time required setpoints.
6. Use **RIGHT DETECTOR** to display the appropriate NPD Detector Control Table. Set the required temperature Base Temp and the detector gases H2, Air and Mkup required setpoints.

RIGHT DETECTOR (NPD)		
Source cur, A		Off
Base temp	300	300
Signal pA		(10.4)
Target curr. pA		(X.XX)
Auto adjust		No
Polarizer V		3.5
H2 delay time		Off
H2	2.3	2.3
Air	60	60
Mkup N2	15	15

7. Scroll to Polarizer V and set 3.5 V.

Source Ignition

8. Switch on the source operating as follows:
 - a. Open the detector gases H₂, Air and Mkup and set the gas flow rates as follows:
 - H₂ = 2.3 ml/min
 - Air = 60 ml/min
 - Mkup N₂ = 15 ml/min)
 - b. Increase the Base Temp to 300°C and wait that the NPD cell reaches the correct set temperature.
 - c. Be sure that the backoff signal is between 0 and 0.5 pA.
 - d. Switch on the source with an initial current of 2.50 A. The backoff signal can slightly increase, but should remain within 0 and 1.5 pA.

- e. Monitor the signal through the keypad or through the data system, increase the current value by steps of 0.002 A, until an immediate and strong increase of the signal is observed.
 - f. Wait five minutes to let the source stabilize.
 - g. Check that source is correctly switched on decreasing hydrogen flow to 0.5 ml/min until signal decreases down to zero, then increase again to original value.
 - If the signal remains around zero, it means that the source is not switched on and it is necessary to increase further the current, accordingly to the procedure just described.
 - If the signal rises back to original value, it means that source is correctly switched on
 - h. Increase the current value of 2% of the actual ignition current. Let the signal stabilize until its level drops below 20 pA.
9. Use **RIGHT SIGNAL** to display the appropriate NPD Detector Signal Control Table. Scroll to `Range` and set the electrometer amplifier input range required.

RIGHT SIGNAL (NPD)	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range $10^{(0..3)}$	0<
Analog filter	Off
Baseline comp	Off

10. Activate your Data System and set the parameters required for the checkout.
11. In the NPD Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
12. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 16-3 according to the data handling in use.

13. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
14. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
15. Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 16.1*.



CAUTION

Verify the peak shape. If any peak distortion is visible, change the analytical test column.

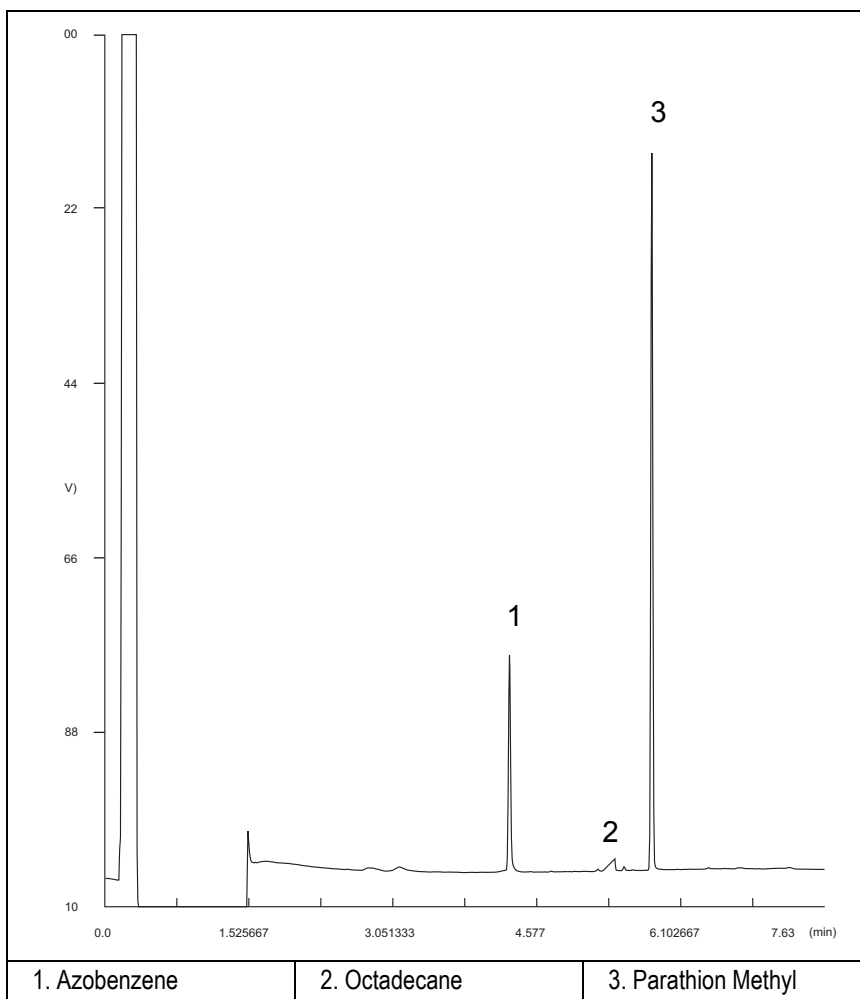


Figure 16-1. NPD-PTV Injection

16. Establish the integration parameters and the peak table identifying the test mix components.
17. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Azobenzene component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Azobenzene.
- Repeat the procedure to calculate the signal-to-noise ratio also for Parathion Methyl.
- Generate a report showing the chromatogram, peak area and signal-to-noise information for Parathion Methyl.



NOTE

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows: Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:

- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.




CAUTION


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


18. The following criteria indicate successful completion of NPD-PTV checkout.


19. If these criteria are not met, repeat the test.


Table 16-3. NPD-PTV Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 40	< 400
	Wander (μV)	< 100	< 1 000
	Drift ($\mu\text{V/h}$)	< 300	<3 000
	Analytical Results		
	Azobenzene Signal-to-noise ratio	> 550	
	Parathion Methyl Signal-to-noise ratio	> 1500	
	Octadecane Signal-to-noise ratio	Negligible	
			

Computing-integrator (e.g. ChromJet)	
	

CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 40
	Wander (μV)	< 100
	Drift ($\mu\text{V/h}$)	< 300
	Analytical Results	
	Azobenzene Signal-to-noise ratio	> 550
	Parathion Methyl Signal-to-noise ratio	> 1500
	Octadecane Signal-to-noise ratio	Negligible
		

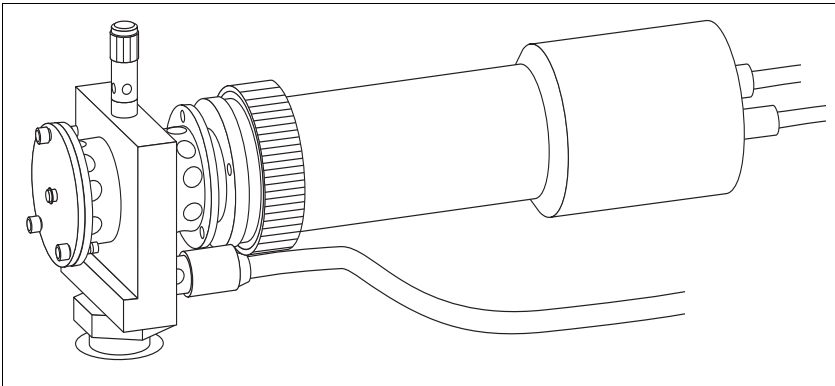
ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 400
Wander (μV)	< 1 000
Drift ($\mu\text{V}/\text{h}$)	< 3 000
Analytical Results	
Azobenzene Signal-to-noise ratio	> 550
Parathion Methyl Signal-to-noise ratio	> 1500
Octadecane Signal-to-noise ratio	Negligible
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 4 000
Wander Counts)	< 10 000
Drift (Counts/h)	< 30 000
Analytical Results	
Azobenzene Signal-to-noise ratio	> 550
Parathion Methyl Signal-to-noise ratio	> 1 500
Octadecane Signal-to-noise ratio	Negligible
	

SECTION

V

SOPs Using FPD



The *SOPs Using Fast FPD* section, contains the procedures to test the TRACE GC Ultra with the fast Flame Photometric Detector (FPD) using different injectors.

Chapter 17, *Checkout Using FPD with S/SL Injector*.

Chapter 18, *Checkout Using FPD with OC Injector*.

Chapter 19, *Checkout Using FPD with PKD Injector*.

Chapter 20, *Checkout Using FPD with PPKD Injector*.

Chapter 21, *Checkout Using FPD with PTV Injector*.

Checkout Using FPD with S/SL Injector

Chapter at a Glance...

SOP Number: P0316/08/E - 01 September 2009	226
Scope	226
Parts Referenced	226
Analytical Conditions Required for Splitless Injection	227
Recommended Initial Operations	228

Operating Procedures

FPD-S/SL Checkout in Splitless Mode	229
---	-----

SOP Number: P0316/08/E - 01 September 2009

Scope

Use the following procedure to verify proper FPD operation with the Split/Splitless Injector. This SOP is applicable both for the control card labeled **FPD** and for the control card labeled **FPD/F**.

Parts Referenced

Table 17-1. FPD-S/SL Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long; 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	3 mm ID for splitless injection	453 200 32
Liner Seal	Graphite seal for glass liner	290 334 06
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
Septum	Standard septum for S/SL injector	313 032 11
Syringe	10 µl size; 70 mm needle length	365 001 03
Interferential Filter	526 nm for phosphorus 394 nm for sulphur	281 071 00 281 070 00
Test Mixture	Three components in Iso-Octane: <i>Component</i> <i>Concentration</i> Azobenzene 1 µg/ml Octadecane 1000 µg/ml Parathion methyl 1 µg/ml	338 190 06
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for Splitless Injection

Table 17-2. FPD-S/SL Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 90 ml/min Air = 115 ml/min
Oven Program	Initial Temperature = 70 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 230 °C Final Time = 1 minute
Injector	Operating Mode = Splitless Temperature = 230 °C Splitless Time = 0.8 minutes Split Flow = 60 ml/min Constant Septum Purge = Yes
Detector	Base Temperature = 300 °C FPD Temperature = 150 °C High voltage mode = No Detector Signal Range = 10 ⁰ (See Note)
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz



NOTE

In the case your GC is equipped with the previous non-fast FPD control card, labeled **FPD**, set Detector Signal Range to 10¹.

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Replace the glass liner.
The glass liner currently installed in your injector should be carefully removed and replaced with the 3 mm ID glass liner for splitless application, as required for the checkout, with the appropriate graphite seal.
2. Replace the septum
A new septum should be installed properly in your injector.
3. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
4. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
5. Perform Column Evaluation and Leak Test
6. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

FPD-S/SL Checkout in Splitless Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	S/SL
Left carrier or Right carrier	He (helium)
Left detector or Right detector	FPD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to Pressure and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		230
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate Split/Splitless Injector Control Table and set the required temperature setpoint **Temp**.

Verify to operate in **Splitless** mode. If not, scroll to Mode, press **MODE/TYPE** to access the selection menu, then select Splitless. Scroll to Splitless time to set the required setpoint.

LEFT INLET (S/SL) ¹		
Temp	230	230
Pressure	30.0	30.0
Mode:	Splitless	
Total flow	(63.0)	
Split Flow	60.0	60.0
Splitless time	0.80	0.80
Constant sept purge?	Y	<

1. These settings could also be for a right inlet.

4. Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate FPD Detector Control Table. Set the required temperatures Base Temp and FPD Temp. Then, set the detector gases H2 and Air required setpoints.

LEFT DETECTOR (FPD) ¹		
Flame	Off	
Base temp	300	300
FPD temp	150	150
Signal pA	(1.4)	
High voltage mode?	N	
H2	90	90
Air	115	115
Mkup N2	00	00

1. These settings could also be for a right detector.

5. Verify that High voltage mode is set to **NO**.
6. Scroll to Flame and press **ON**. This start the ignition sequence. When ignition is confirmed, the photomultiplier tube is energized.

The baseline level `Signal pA`, will suddenly increase meaning that the flame is lit inside the detector. After a few seconds, the baseline should stabilize to the standing current of the system.

- Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate FPD Detector Signal Control Table. Scroll to `Range` and set the electrometer amplifier input range required.

LEFT SIGNAL (FPD) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Baseline comp	Off

1. These settings could also be for a right signal.

- Activate your Data System and set the parameters required for the checkout.
- In the FPD Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
- Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 17-3 according to the data handling in use.

- With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
- After the baseline evaluation has been completed, set-up the data system to acquire a single run.
- When the GC is ready, inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 17.1* or *17.2*.

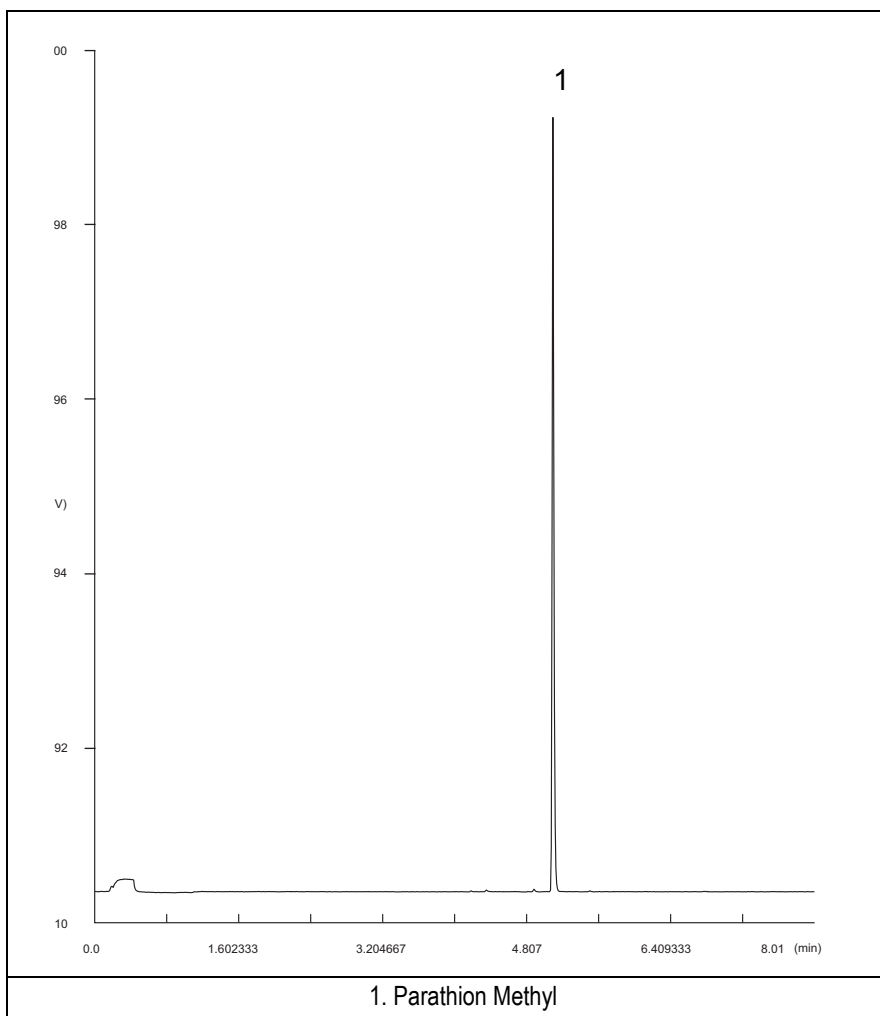


Figure 17-1. FPD-Splitless Injection with 526 nm filter

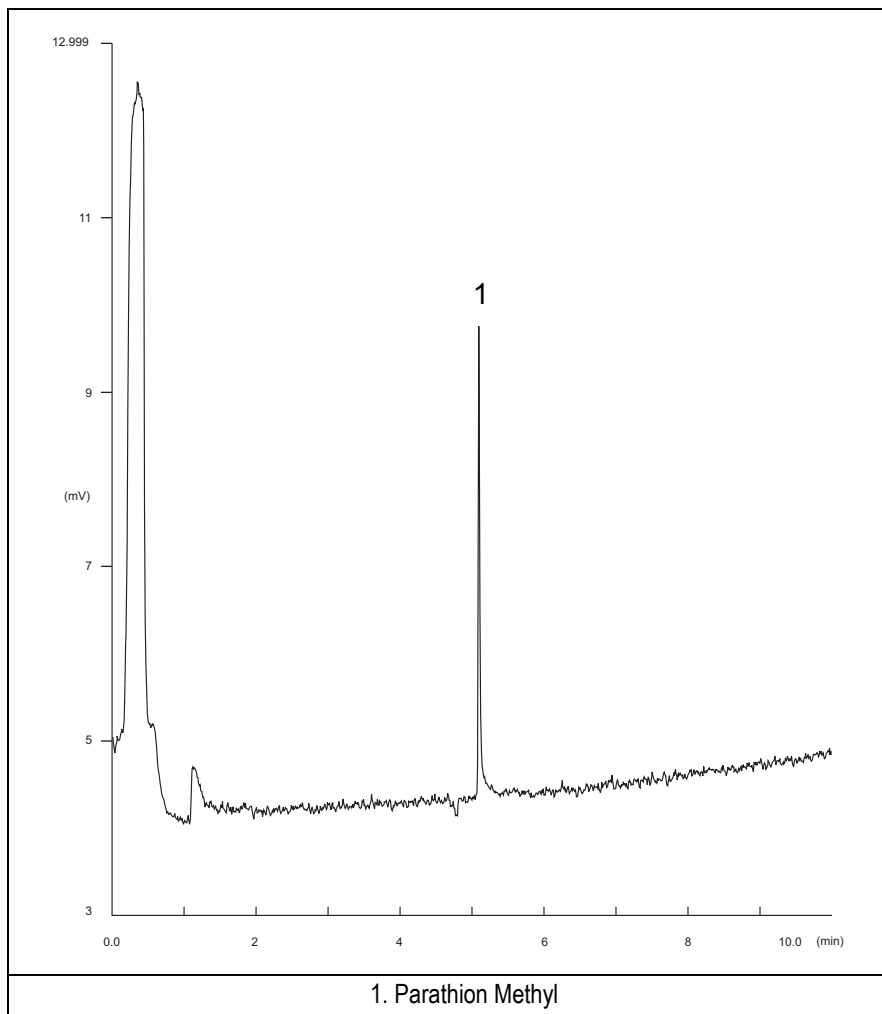


Figure 17-2. FPD-Splitless Injection with 394 nm

14. Establish the integration parameters and the peak table identifying the test mix components.
15. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Parathion Methyl component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for the component.



NOTE

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows: Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:

- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.





CAUTION


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


16. The following criteria indicate successful completion of FPD-S/SL checkout.
17. If these criteria are not met, repeat the test.

Table 17-3. FPD-S/SL Acceptance Criteria

Acceptance Values	CHROM-CARD				
	Baseline Parameters	Analog (1V Full Scale)		Digital (10V Full Scale)	
		394 nm (S) Filter	526 nm (P) Filter	394 nm (S) Filter	526 nm (P) Filter
	Noise (μV)	< 50	< 50	< 500	< 500
	Wander (μV)	< 100	< 100	< 1 000	< 1 000
Drift ($\mu\text{V/h}$)	< 100	< 100	< 1 000	< 1 000	
Analytical Results					
Parathion Methyl Signal to Noise Ratio			394 nm (S) Filter	526 nm (P) Filter	
			> 40	> 2 000	
					

Acceptance Values	CHROMQUEST		
	Baseline Parameters (1V Full Scale)		
			526 nm (P) Filter
	Noise (μV)	< 50	< 50
	Wander (μV)	< 100	< 100
Drift ($\mu\text{V/h}$)	< 100	< 100	
Analytical Results			
Parathion Methyl Signal to Noise Ratio		394 nm (S) Filter	526 nm (P) Filter
		> 40	> 2 000
			

ATLAS		
Baseline Parameters (10V Full Scale)		
		526 nm (P) Filter
Noise (μV)	< 500	< 500
Wander (μV)	< 1 000	< 1 000
Drift ($\mu\text{V}/\text{h}$)	< 1 000	< 1 000
Analytical Results		
Parathion Methyl Signal to Noise Ratio	394 nm (S) Filter	526 nm (P) Filter
	> 40	> 2 000
		

XCALIBUR			
Baseline Parameters (Acquisition Frequency = 10 Hz)			
Acceptance Values		526 nm (P) Filter	
	Noise (Counts)	< 5 000	< 5 000
	Wander (Counts)	< 10 000	< 10 000
	Drift (Counts/h)	< 10 000	< 10 000
Analytical Results			
Parathion Methyl Signal to Noise Ratio	394 nm (S) Filter	526 nm (P) Filter	
	> 40	> 2 000	
			

Analytical Acceptance Comments	
1	Using Chrom-Card, set the signal-to-noise report parameters as described in the current procedure.
2	Using ChromQuest, Atlas, Xcalibur or a Computing integrator (e.g. ChromJet), calculate the S/N ratio as <i>Peak Height (counts)/noise (counts)</i> .

Checkout Using FPD with OC Injector

Chapter at a Glance...

SOP Number: P0317/09/E - 01 September 2009	238
Scope	238
Parts Referenced	238
Analytical Conditions Required for On-Column Injection	239
Recommended Initial Operations	240

Operating Procedures

FPD-OCI Checkout in On-Column Mode	241
--	-----

SOP Number: P0317/09/E - 01 September 2009

Scope

Use the following procedure to verify proper FPD operation with the On-Column Injector. This SOP is applicable both for the control card labeled **FPD** and for the control card labeled **FPD/F**.

Parts Referenced

Table 18-1. FPD-OCI Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long; 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Graphite Ferrule	Graphite Ferrule for 0.32 mm ID Column	290 134 87
Vespel Ferrule	Vespel Ferrule for 0.32 mm ID Column	290 134 60
Retaining Nut	M4 capillary column retaining nut	350 324 23
Syringe	10 µl size; 75 mm needle length	365 020 07
Interferential Filter	526 nm for phosphorus 394 nm for sulphur	281 071 00 281 070 00
Test Mixture	Three components in Iso-Octane: <i>Component</i> <i>Concentration</i> Azobenzene 1 µg/ml Octadecane 1000 µg/ml Parathion methyl 1 µg/ml	338 190 06
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	
In case of automatic On-Column for TriPlus Sampler AS		
Syringe	10 µl size; 80 mm needle length	365 020 19

Table 18-1. FPD-OCI Parts Referenced (Continued)

Part	Description	Part Number
Pre-column	2 m long; 0.53 mm ID	260 603 75
Press-fit set	Set of Press-fit connectors for TRACE OC	350 038 45

Analytical Conditions Required for On-Column Injection

Table 18-2. FPD-OCI Analytical Conditions

Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 90 ml/min Air = 115 ml/min
Oven Program	Initial Temperature = 85 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 230 °C Final Time = 1 minute
Injector	Secondary cooling = 10 seconds
Detector	Base Temperature = 300 °C FPD Temperature = 150 °C High voltage mode = No Detector Signal Range = 10 ⁰ (see Note)
Injected Volume	1 µl of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

**NOTE**

In the case your GC is equipped with the previous non-fast FPD control card, labeled **FPD**, set Detector Signal Range to 10¹.

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
2. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
In case of automatic On-column for TriPlus sampler, install the pre-column and connect it to the test column by press-fit connector.
3. Install and connect the TriPlus sampler and its components.
4. Perform Column Evaluation and Leak Test.
5. Connect your data handling.
Verify that your data handling is properly connected to your GC system.
6. Verify the opening/closing of the OC injector actuator by using the proper commands.
7. Verify the alignment of the syringe on the OC injector.

OPERATING PROCEDURE

FPD-OCI Checkout in On-Column Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	OCI
Left carrier or Right carrier	He (helium)
Left detector or Right detector	FPD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to Pressure and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	85.0	85.0
Initial Time		1.00
Ramp 1		20.0
Final temp		230
Final time 1		1.00 <
Ramp 2		Off

- Use **LEFT INLET** or **RIGHT INLET** to display the appropriate On-Column Injector Control Table. Scroll to *Sec. cool time* and set the required secondary cooling time.

LEFT INLET (OCI) ¹		
Pressure	30.0	30.0
Sec. Cool Time		10.00<

1. These settings could also be for a right inlet.

- Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate FPD Detector Control Table. Set the required temperatures *Base Temp* and *FPD Temp*. Then, set the detector gases *H2* and *Air* required setpoints.

LEFT DETECTOR (FPD) ¹		
Flame		Off
Base temp	300	300
FPD temp	150	150
Signal pA		(1.4)
High voltage mode?		N
H2	90	90
Air	115	115
Mkup N2	00	00

1. These settings could also be for a right detector.

- Verify that *High voltage mode* is set to **NO**.
- Scroll to *Flame* and press **ON**. This start the ignition sequence. When ignition is confirmed, the photomultiplier tube is energized. The baseline level *Signal pA*, will suddenly increase meaning that the flame is lit inside the detector. After a few seconds, the baseline should stabilize to the standing current of the system.
- Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate FPD Detector Signal Control Table. Scroll to *Range* and set the electrometer amplifier input range required.

LEFT SIGNAL (FPD) ¹	
Output	(1000)
Offset	100
Autozero?	Y/N
Range 10^(0..3)	0<
Baseline comp	Off

1. These settings could also be for a right signal.

8. Activate your Data System and set the parameters required for the checkout.
9. In the FPD Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
10. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 18-3 according to the data handling in use.

11. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
12. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
13. Perform the analysis.

Manual injection

- Inject the test mixture and press **START** on the GC to begin the checkout run.

Automatic injection with TriPlus sampler

- Fill a vial with the standard mix and place that vial in the sample tray.
- Load the method for OC and perform the sampling..

The resulting chromatogram should look like the one shown in *Figure 18.1* or *18.2*.

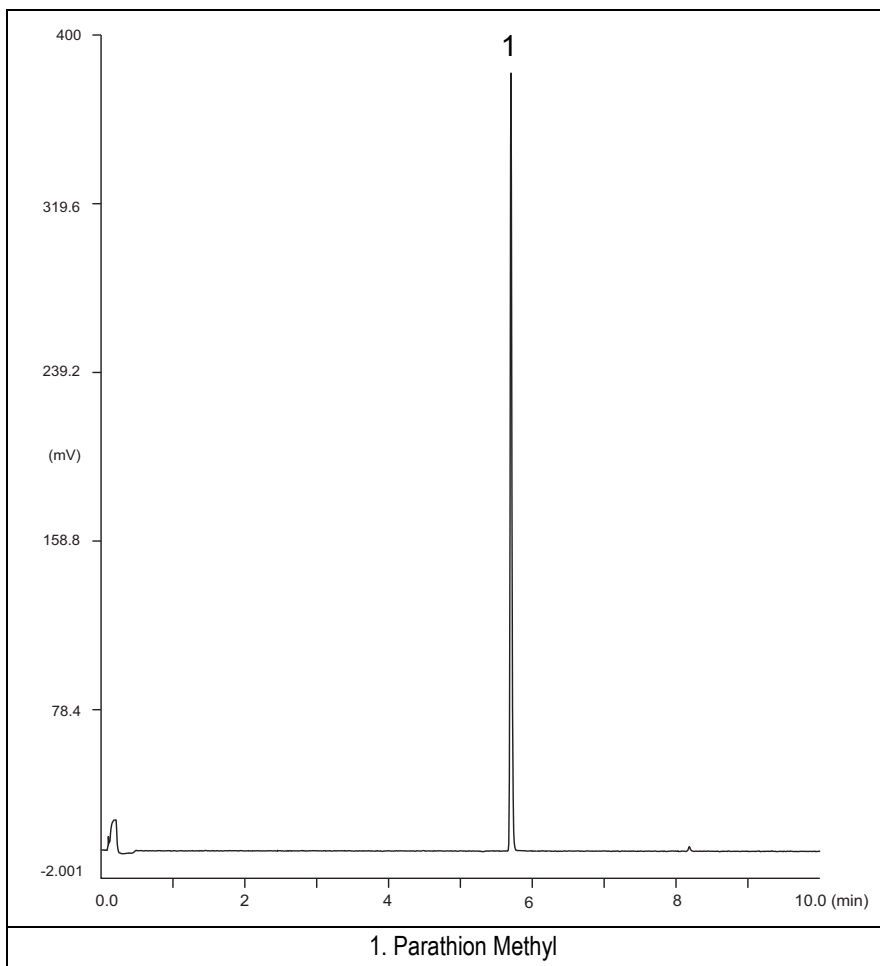


Figure 18-1. FPD-On-Column Injection with 526 nm filter

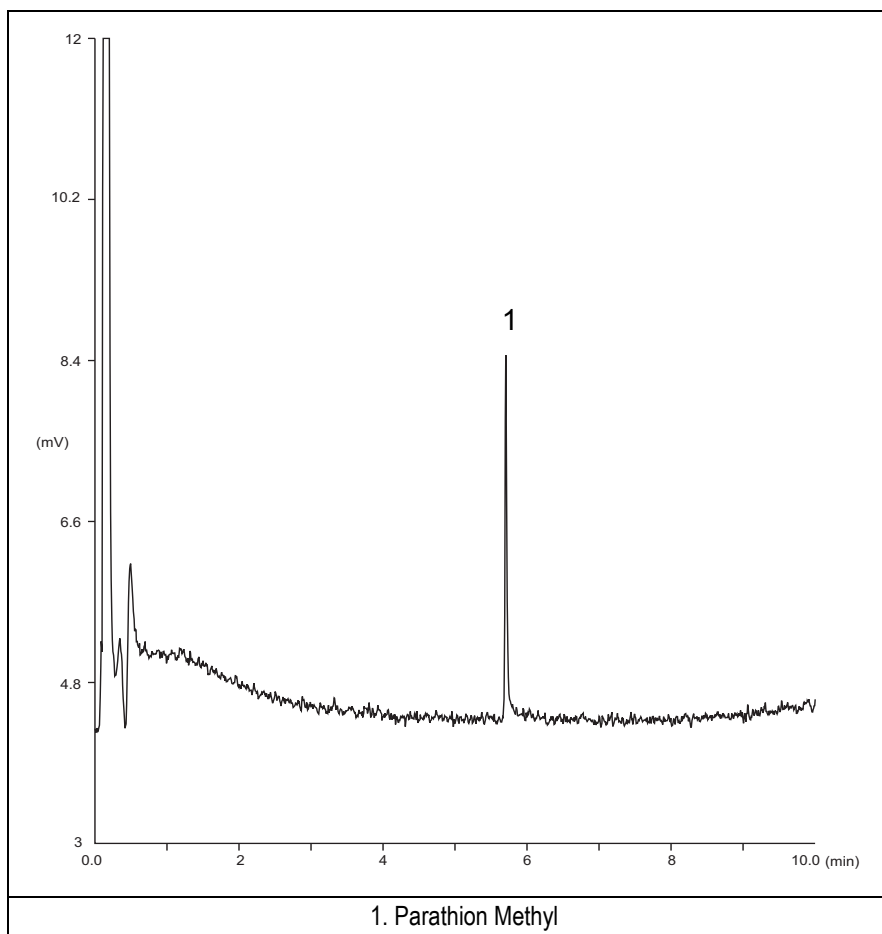


Figure 18-2. FPD-On-Column Injection with 394 nm

14. Establish the integration parameters and the peak table identifying the test mix components.
15. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Methylparathion component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for the component.



NOTE

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows: Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:

- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.





CAUTION


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


16. The following criteria indicate successful completion of FPD-OCI checkout.
17. If these criteria are not met, repeat the test.

Table 18-3. FPD-OCI Acceptance Criteria

Acceptance Values	CHROM-CARD				
	Baseline Parameters	Analog (1V Full Scale)		Digital (10V Full Scale)	
		394 nm (S) Filter	526 nm (P) Filter	394 nm (S) Filter	526 nm (P) Filter
	Noise (μV)	< 50	< 50	< 500	< 500
	Wander (μV)	< 100	< 100	< 1 000	< 1 000
Drift ($\mu\text{V/h}$)	< 100	< 100	< 1 000	< 1 000	
Analytical Results					
Parathion Methyl Signal to Noise Ratio			394 nm (S) Filter	526 nm (P) Filter	
			> 20	> 1 000	
					

Acceptance Values	CHROMQUEST		
	Baseline Parameters (1V Full Scale)		
			526 nm (P) Filter
	Noise (μV)	< 50	< 50
	Wander (μV)	< 100	< 100
Drift ($\mu\text{V/h}$)	< 100	< 100	
Analytical Results			
Parathion Methyl Signal to Noise Ratio		394 nm (S) Filter	526 nm (P) Filter
		> 20	> 1 000
			

ATLAS		
Baseline Parameters (10V Full Scale)		
		526 nm (P) Filter
Noise (μV)	< 500	< 500
Wander (μV)	< 1 000	< 1 000
Drift ($\mu\text{V}/\text{h}$)	< 1 000	< 1 000
Analytical Results		
Parathion Methyl Signal to Noise Ratio	394 nm (S) Filter	526 nm (P) Filter
	> 20	> 1 000
		

XCALIBUR			
Baseline Parameters (Acquisition Frequency = 10 Hz)			
Acceptance Values		526 nm (P) Filter	
	Noise (Counts)	< 5 000	< 5 000
	Wander (Counts)	< 10 000	< 10 000
	Drift (Counts/h)	< 10 000	< 10 000
Analytical Results			
Parathion Methyl Signal to Noise Ratio	394 nm (S) Filter	526 nm (P) Filter	
	> 20	> 1 000	
			

Analytical Acceptance Comments	
1	Using Chrom-Card, set the signal-to-noise report parameters as described in the current procedure.
2	Using ChromQuest, Atlas, Xcalibur or a Computing integrator (e.g. ChromJet), calculate the S/N ratio as <i>Peak Height (counts)/noise (counts)</i> .

Checkout Using FPD with PKD Injector

Chapter at a Glance...

SOP Number: P0318/08/E - 01 September 2009	250
Scope	250
Parts Referenced	250
Analytical Conditions Required for Packed Injector.....	252
Recommended Initial Operations	253

Operating Procedures

FPD-PKD Checkout	254
------------------------	-----

SOP Number: P0318/08/E - 01 September 2009

Scope

Use the following procedure to verify proper FPD operation with Packed Injector. This SOP is applicable both for the control card labeled **FPD** and for the control card labeled **FPD/F**.

Parts Referenced

Table 19-1. FPD-PKD Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	2 mm ID tapered	453 220 50
Liner Seal	Graphite seal for glass liner	290 334 05
Pre-column	Fused Silica Capillary Column 0.5 mt long 0.53 mm ID	260 603 75
Retaining Nut	M4 capillary column retaining nut	350 324 23
Press-fit connections	For columns 0.53/0.32 mm ID	350 438 16
Injection side adapter	For Wide bore column	347 003 03
Retaining Nut	For Injection side adapter	350 024 04
Detector side adapter	For Wide bore column	347 103 04
Ferrule	6 mm ID double brass ferrule	290 341 37
Nut	1/4" G-6 mm ID nut	350 201 18
Graphite Ferrule	Graphite ferrule for 0.53 mm ID Column Graphite ferrule for 0.32 mm ID Column	290 134 86 290 134 87
Septum	Standard septum for Packed Injector	313 032 26
Syringe	10 µl size; 50 mm needle length	365 005 25

Table 19-1. FPD-PKD Parts Referenced (Continued)

Part	Description	Part Number
Interferential Filter	526 nm for phosphorus 394 nm for sulphur	281 071 00 281 070 00
Test Mixture	Three components in Iso-Octane: <i>Component</i> <i>Concentration</i> Azobenzene 1 µg/ml Octadecane 1000 µg/ml Parathion methyl 1 µg/ml	338 190 06
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for Packed Injector

Table 19-2. FPD-PKD Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 90 ml/min Air = 115 ml/min
Oven Program	IsoTemperature = 70 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 230 °C Final Time = 1 minute
Injector	Operating Mode = Packed Temperature = 200 °C
Detector	Base Temperature = 300 °C FPD Temperature = 150 °C High voltage mode = No Detector Signal Range = 10 ⁰ (see Note)
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz



NOTE

In the case your GC is equipped with the previous non-fast FPD control card, labeled **FPD**, set Detector Signal Range to 10¹.

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Insert the glass liner.
Remove the column and insert the 2 mm ID tapered glass liner, as required for the checkout, from the bottom of the injector with the appropriate liner seal. Fix the liner by using the appropriate adapter for capillary column. Refer to the TRACE GC Ultra Maintenance and Troubleshooting Manual.
2. Mount the adapter for capillary column on the detector base body.
3. Replace the septum
A new septum should be installed properly in your injector.
4. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
5. Install the precolumn.
Connect the precolumn to the injector.
6. Install the test column
Connect the test column to the precolumn by using the press fit connections provided.
7. Perform the manual leak check following the Operating Procedure “*Performing a Leak Check*” in Chapter 14 of the TRACE GC Ultra Operating Manual.
8. Connect the other end of the test column to the detector base body.
9. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

FPD-PKD Checkout

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	PKD
Left carrier or Right carrier	He (helium)
Left detector or Right detector	FPD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select con pres. Scrool to Pressure and set the pressure value to have the required carrier gas flow rate Col. flow.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		230
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate PKD Injector Control Table. Set the required temperature setpoint Temp. Verify to operate in

Packed mode. If not, scroll to Mode, press **MODE/TYPE** to access the selection menu, then select Packed.

LEFT INLET (PKD) ¹		
Temp	200	200
Pressure	30.0	30.0
Mode:	Packed	

1. These settings could also be for a right inlet.

- Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate FPD Detector Control Table. Set the required temperatures Base Temp and FPD Temp. Then, set the detector gases H₂ and Air required setpoints.

LEFT DETECTOR (FPD) ¹		
Flame		Off
Base temp	300	300
FPD temp	150	150
Signal pA		(1.4)
High voltage mode?		N
H ₂	90	90
Air	115	115
Mkup N ₂	00	00

1. These settings could also be for a right detector.

- Verify that High voltage mode is set to **NO**.
- Scroll to Flame and press **ON**. This start the ignition sequence. When ignition is confirmed, the photomultiplier tube is energized. The baseline level Signal pA, will suddenly increase meaning that the flame is lit inside the detector. After a few seconds, the baseline should stabilize to the standing current of the system.

7. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate FPD Detector Signal Control Table. Scroll to `Range` and set the electrometer amplifier input range required.

LEFT SIGNAL (FPD) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Baseline comp	Off

1. These settings could also be for a right signal.

8. Activate your Data System and set the parameters required for the checkout.
9. In the FPD Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
10. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 19-3 according to the data handling in use.

11. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
12. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
13. When the GC is ready, inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 19.1* or *19.2*.

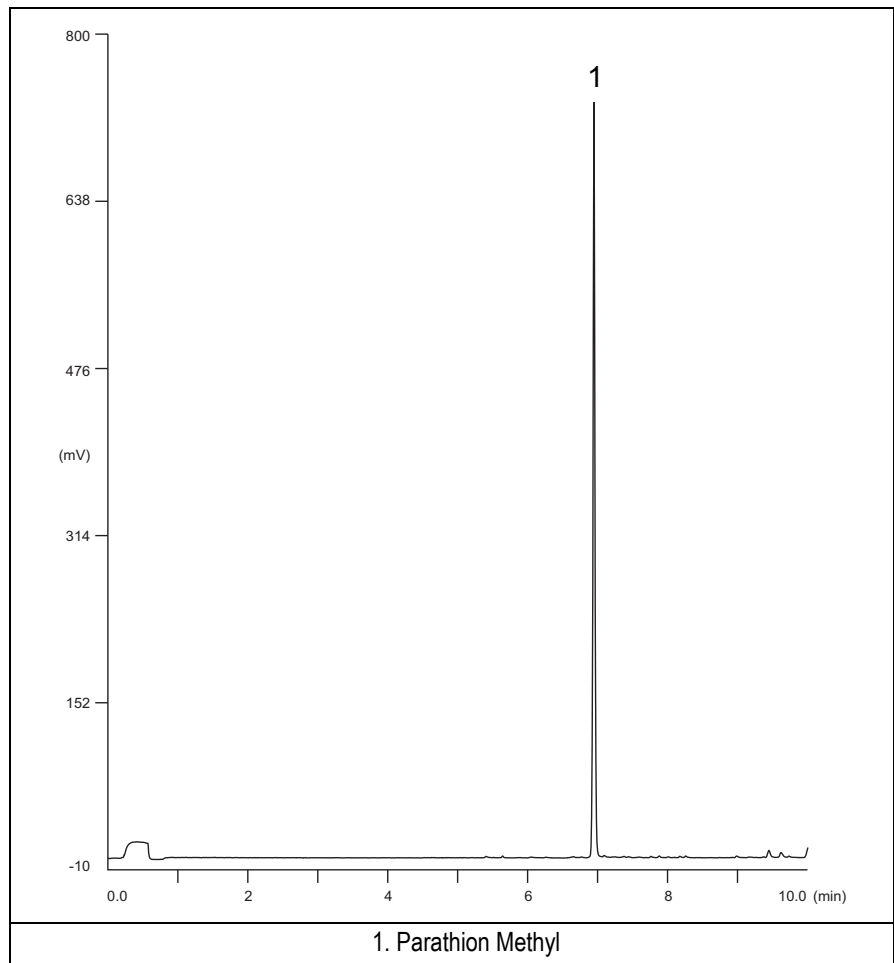


Figure 19-1. FPD-PKD Injection with 526 nm filter

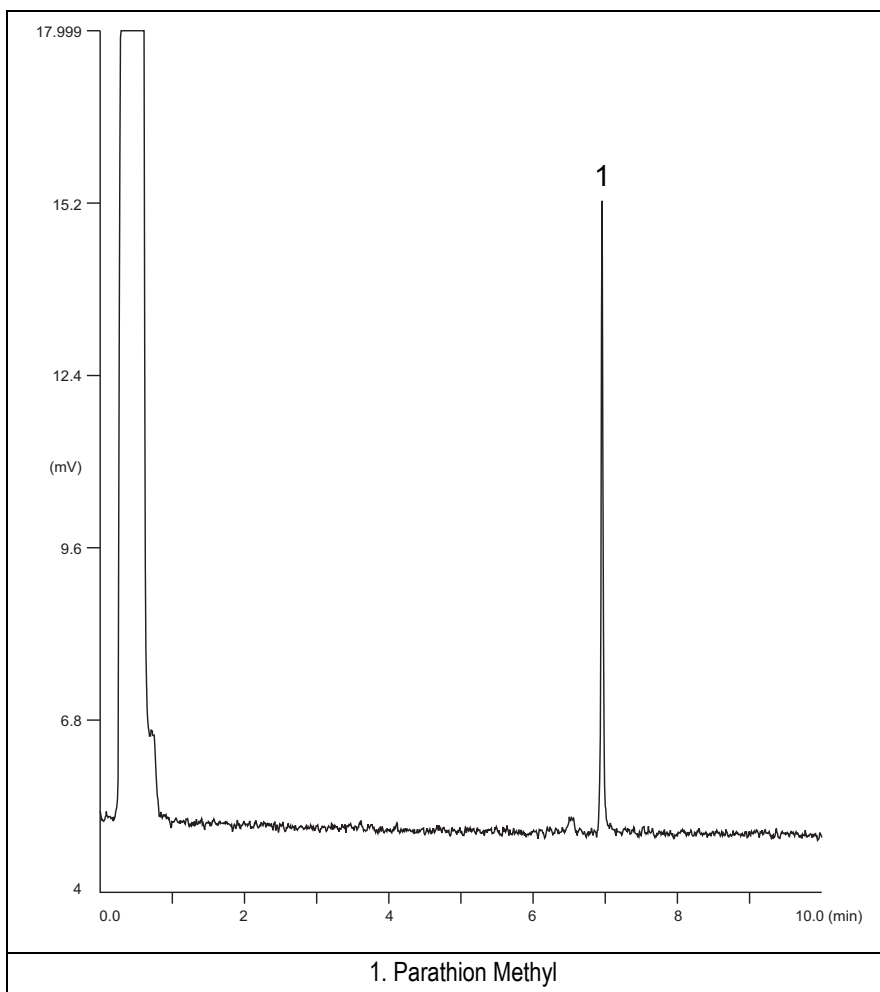


Figure 19-2. FPD-PKD Injection with 394 nm

14. Establish the integration parameters and the peak table identifying the test mix components.
15. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Methylparathion component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for the component.



NOTE

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows: Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:

- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.





CAUTION


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


16. The following criteria indicate successful completion of FPD-PKD checkout.
17. If these criteria are not met, repeat the test.

Table 19-3. FPD-PKD Acceptance Criteria

CHROM-CARD					
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)		Digital (10V Full Scale)	
		394 nm (S) Filter	526 nm (P) Filter	394 nm (S) Filter	526 nm (P) Filter
	Noise (μV)	< 50	< 50	< 500	< 500
	Wander (μV)	< 100	< 100	< 1 000	< 1 000
	Drift ($\mu\text{V/h}$)	< 100	< 100	< 1 000	< 1 000
Analytical Results					
Parathion Methyl Signal to Noise Ratio			394 nm (S) Filter	526 nm (P) Filter	
			> 30	> 1 500	
					

CHROMQUEST			
Acceptance Values	Baseline Parameters (1V Full Scale)		
		526 nm (P) Filter	
	Noise (μV)	< 50	< 50
	Wander (μV)	< 100	< 100
	Drift ($\mu\text{V/h}$)	< 100	< 100
Analytical Results			
Parathion Methyl Signal to Noise Ratio		394 nm (S) Filter	
		> 30	
			

ATLAS		
Baseline Parameters (10V Full Scale)		
		526 nm (P) Filter
Noise (μV)	< 500	< 500
Wander (μV)	< 1 000	< 1 000
Drift ($\mu\text{V/h}$)	< 1 000	< 1 000
Analytical Results		
Parathion Methyl Signal to Noise Ratio	394 nm (S) Filter	526 nm (P) Filter
	> 30	> 1 500
		

XCALIBUR			
Baseline Parameters (Acquisition Frequency = 10 Hz)			
Acceptance Values		526 nm (P) Filter	
	Noise (Counts)	< 5 000	< 5 000
	Wander (Counts)	< 10 000	< 10 000
	Drift (Counts/h)	< 10 000	< 10 000
	Analytical Results		
Parathion Methyl Signal to Noise Ratio	394 nm (S) Filter	526 nm (P) Filter	
	> 30	> 1 500	
			

Analytical Acceptance Comments	
1	Using Chrom-Card, set the signal-to-noise report parameters as described in the current procedure.
2	Using ChromQuest, Atlas, Xcalibur or a Computing integrator (e.g. ChromJet), calculate the S/N ratio as <i>Peak Height (counts)/noise (counts)</i> .

Checkout Using FPD with PPKD Injector

Chapter at a Glance...

SOP Number: P0319/08/E - 01 September 2009	264
Scope	264
Parts Referenced.....	264
Analytical Conditions Required for Purged Packed Injector	266
Recommended Initial Operations	267

Operating Procedures

FPD-PPKD Checkout.....	268
------------------------	-----

SOP Number: P0319/08/E - 01 September 2009

Scope

Use the following procedure to verify proper FPD operation with Purged Packed Injector. This SOP is applicable both for the control card labeled **FPD** and for the control card labeled **FPD/F**.

Parts Referenced

Table 20-1. FPD-PPKD Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	2 mm ID tapered	453 220 50
Liner Seal	Graphite seal for glass liner	290 334 05
Pre-column	Fused Silica Capillary Column 0.5 mt long 0.53 mm ID	260 603 75
Retaining Nut	M4 capillary column retaining nut	350 324 23
Press-fit connections	For columns 0.53/0.32 mm ID	350 438 16
Injection side adapter	For Wide bore column	347 003 03
Retaining Nut	For Injection side adapter	350 024 04
Detector side adapter	For Wide bore column	347 103 04
Ferrule	6 mm ID double brass ferrule	290 341 37
Nut	1/4" G-6 mm ID nut	350 201 18
Graphite Ferrule	Graphite ferrule for 0.53 mm ID Column Graphite ferrule for 0.32 mm ID Column	290 134 86 290 134 87
Septum	Standard septum for Purged Packed Injector	313 032 26
Syringe	10 µl size; 50 mm needle length	365 005 25

Table 20-1. FPD-PPKD Parts Referenced (Continued)

Part	Description	Part Number
Interferential Filter	526 nm for phosphorus	281 071 00
	394 nm for sulphur	281 070 00
Test Mixture	Three components in Iso-Octane: <i>Component</i> <i>Concentration</i> Azobenzene 1 µg/ml Octadecane 1000 µg/ml Parathion methyl 1 µg/ml	338 190 06
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for Purged Packed Injector

Table 20-2. FPD-PPKD Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 90 ml/min Air = 115 ml/min
Oven Program	Initial Temperature = 50 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 200 °C Final Time = 1 minute
Injector	Operating Mode = Wide bore Temperature = 200 °C
Detector	Base Temperature = 300 °C FPD Temperature = 150 °C High voltage mode = No Detector Signal Range = 10 ⁰ (see Note)
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz



NOTE

In the case your GC is equipped with the previous non-fast FPD control card, labeled **FPD**, set Detector Signal Range to 10¹.

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Insert the glass liner.
Remove the column and insert the 2 mm ID tapered glass liner, as required for the checkout, from the bottom of the injector with the appropriate liner seal. Fix the liner by using the appropriate adapter for capillary column. Refer to the TRACE GC Ultra Maintenance and Troubleshooting Manual.
2. Mount the adapter for capillary column on the detector base body.
3. Replace the septum
A new septum should be installed properly in your injector.
4. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
5. Install the precolumn.
Connect the precolumn to the injector.
6. Install the test column
Connect the test column to the precolumn by using the press fit connections provided.
7. Perform the manual leak check following the Operating Procedure “*Performing a Leak Check*” in Chapter 14 of the TRACE GC Ultra Operating Manual.
8. Connect the other end of the test column to the detector base body.
9. Perform Column Evaluation.
10. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

FPD-PPKD Checkout

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	PPKD
Left carrier or Right carrier	He (helium)
Left detector or Right detector	FPD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select con pres. Scrool to Pressure and set the pressure value to have the required carrier gas flow rate Col.flow.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col.flow	3.00	
Lin. veloc.		(60.9) <

These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		250
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate PPKD Injector Control Table. Set the required temperature setpoint Temp. Verify to operate in

Widebore mode. If not, scroll to **Mode**, press **MODE/TYPE** to access the selection menu, then select **Widebore**.

LEFT INLET (PPKD) ¹		
Temp	200	200
Pressure	30.0	30.0
Mode:	Wide bore<	
Constant sept purge?		Y<

These settings could also be for a right inlet.

- Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate FPD Detector Control Table. Set the required temperatures **Base Temp** and **FPD Temp**. Then, set the detector gases **H2** and **Air** required setpoints.

LEFT DETECTOR (FPD) ¹		
Flame		Off
Base temp	300	300
FPD temp	150	150
Signal pA		(1.4)
High voltage mode?		N
H2	90	90
Air	115	115
Mkup N2	00	00

These settings could also be for a right detector.

- Verify that **High voltage mode** is set to **NO**.
- Scroll to **Flame** and press **ON**. This start the ignition sequence. When ignition is confirmed, the photomultiplier tube is energized. The baseline level **Signal pA**, will suddenly increase meaning that the flame is lit inside the detector. After a few seconds, the baseline should stabilize to the standing current of the system.

7. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate FPD Detector Signal Control Table. Scroll to `Range` and set the electrometer amplifier input range required.

LEFT SIGNAL (FPD) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Baseline comp	Off

These settings could also be for a right signal.

8. Activate your Data System and set the parameters required for the checkout.
9. In the FPD Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
10. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 20-3 according to the data handling in use.

11. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
12. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
13. When the GC is ready, inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 20.1* or *20.2*.

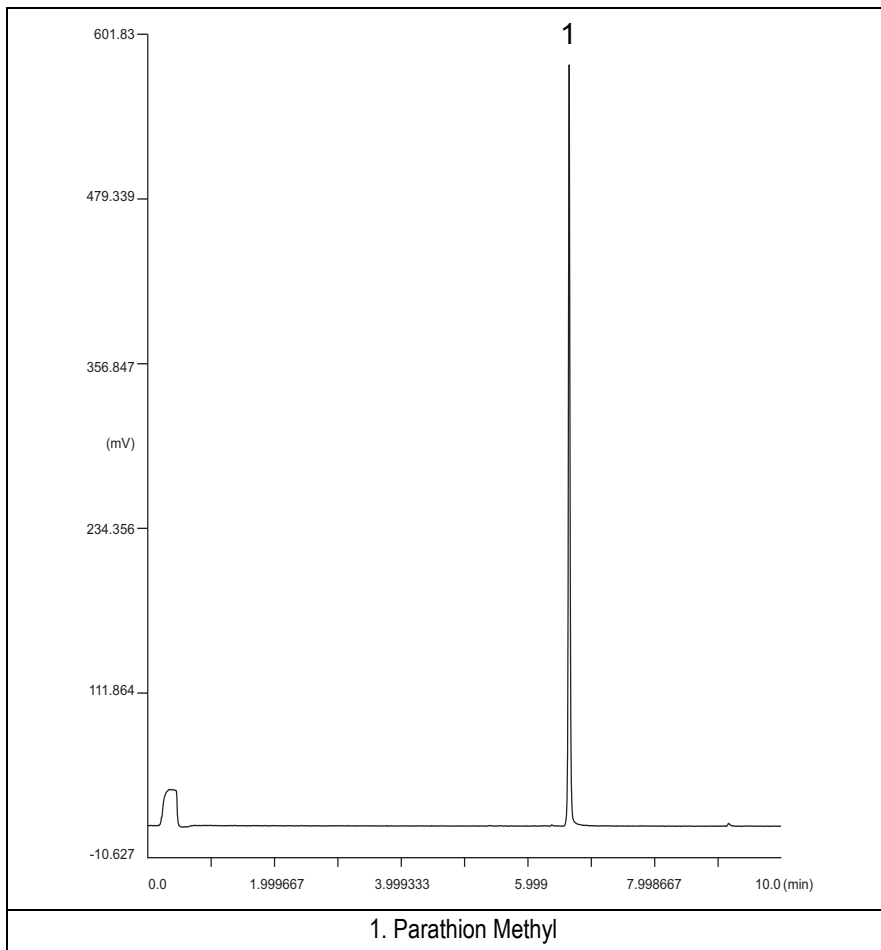


Figure 20-1. FPD-PPKD Injection with 526 nm filter

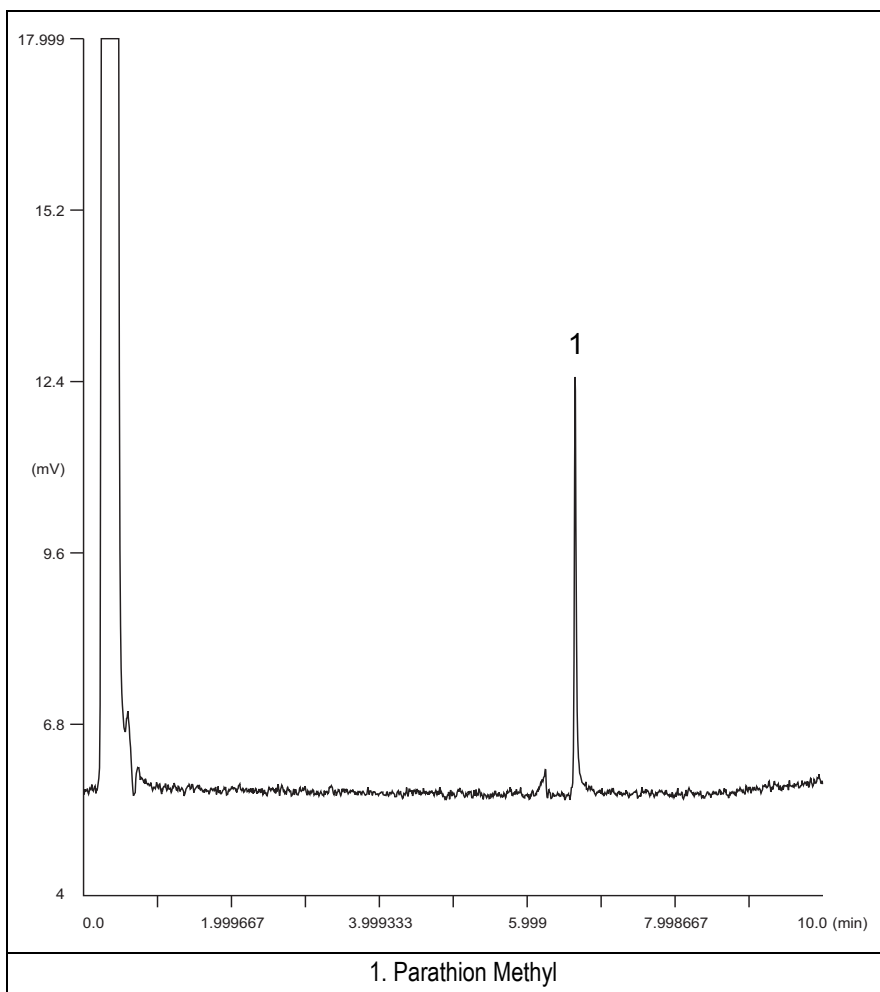


Figure 20-2. FPD-PPKD Injection with 394 nm

14. Establish the integration parameters and the peak table identifying the test mix components.
15. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Methylparathion component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for the component.



NOTE

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows: Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:

- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.





CAUTION


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


16. The following criteria indicate successful completion of FPD-PPKD checkout.
17. If these criteria are not met, repeat the test.

Table 20-3. FPD-PPKD Acceptance Criteria

CHROM-CARD					
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)		Digital (10V Full Scale)	
		394 nm (S) Filter	526 nm (P) Filter	394 nm (S) Filter	526 nm (P) Filter
	Noise (μV)	< 50	< 50	< 500	< 500
	Wander (μV)	< 100	< 100	< 1 000	< 1 000
	Drift ($\mu\text{V}/\text{h}$)	< 100	< 100	< 1 000	< 1 000
Analytical Results					
Parathion Methyl Signal to Noise Ratio			394 nm (S) Filter	526 nm (P) Filter	
			> 30	> 1 500	
					

CHROMQUEST			
Acceptance Values	Baseline Parameters (1V Full Scale)		
		526 nm (P) Filter	
	Noise (μV)	< 50	< 50
	Wander (μV)	< 100	< 100
	Drift ($\mu\text{V}/\text{h}$)	< 100	< 100
Analytical Results			
Parathion Methyl Signal to Noise Ratio		394 nm (S) Filter	
		> 30	
			

ATLAS		
Baseline Parameters (10V Full Scale)		
		526 nm (P) Filter
Noise (μV)	< 500	< 500
Wander (μV)	< 1 000	< 1 000
Drift ($\mu\text{V}/\text{h}$)	< 1 000	< 1 000
Analytical Results		
Parathion Methyl Signal to Noise Ratio	394 nm (S) Filter	526 nm (P) Filter
	> 30	> 1 500
		

XCALIBUR			
Baseline Parameters (Acquisition Frequency = 10 Hz)			
Acceptance Values		526 nm (P) Filter	
	Noise (Counts)	< 5 000	< 5 000
	Wander (Counts)	< 10 000	< 10 000
	Drift (Counts/h)	< 10 000	< 10 000
	Analytical Results		
Parathion Methyl Signal to Noise Ratio	394 nm (S) Filter	526 nm (P) Filter	
	> 30	> 1 500	
			

Analytical Acceptance Comments	
1	Using Chrom-Card, set the signal-to-noise report parameters as described in the current procedure.
2	Using ChromQuest, Atlas, Xcalibur or a Computing integrator (e.g. ChromJet), calculate the S/N ratio as <i>Peak Height (counts)/noise (counts)</i> .

Checkout Using FPD with PTV Injector

Chapter at a Glance...

SOP Number: P0320/08/E - 01 September 2009	278
Scope	278
Parts Referenced	278
Analytical Conditions Required for PTV Splitless Injection	279
Recommended Initial Operations	280

Operating Procedures

FPD-PTV Checkout in PTV Splitless Mode	281
--	-----

SOP Number: P0320/08/E - 01 September 2009

Scope

Use the following procedure to verify proper FPD operation with the Programmable Temperature Vaporizing Injector. This SOP is applicable both for the control card labeled **FPD** and for the control card labeled **FPD/F**.

Parts Referenced

Table 21-1. FPD-PTV Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Liner	Silcosteel 2 mm ID (set of 2)	453 220 44
Liner Seal	Graphite seal for liner	290 034 17
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
Septum	Standard septum for PTV injector (set of 10)	313 132 25
Syringe	10 µl size; 50 mm needle length	365 005 25
Interferential Filter	526 nm for phosphorus 394 nm for sulphur	281 071 00 281 070 00
Test Mixture	Three components in Iso-Octane: <i>Component</i> <i>Concentration</i> Azobenzene 1 µg/ml Octadecane 1000 µg/ml Parathion methyl 1 µg/ml	338 190 06
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for PTV Splitless Injection

Table 21-2. FPD-PTV Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 90 ml/min Air = 115 ml/min
Oven Program	Initial Temperature = 70 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 230 °C Final Time = 1 minute
Injector	Operating Mode = PTV splitless Splitless Time = 0.8 minutes Split Flow = 50 ml/min Constant Septum Purge = Yes Inject Temp = 50 °C Inject Time = 0.1 minute Transfer ramp = 10 °C/sec Transfer Temperature = 260 °C Transfer time = 1 minute
Detector	Base Temperature = 300 °C FPD Temperature = 150 °C High voltage mode = No Detector Signal Range = 10 ⁰ (see Note)
Injected Volume	1 µl of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz



NOTE

In the case your GC is equipped with the previous non-fast FPD control card, labeled **FPD**, set Detector Signal Range to 10^1 .

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Replace the liner.
The liner currently installed in your injector should be carefully removed and replaced with the 2 mm ID Silcolsteel liner, as required for the checkout, with the appropriate liner seal.
2. Replace the septum
A new septum should be installed properly in your injector.
3. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
4. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
5. Perform Column Evaluation and Leak Test.
6. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

FPD-PTV Checkout in PTV Splitless Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Right inlet	PTV
Right carrier	He (helium)
Right detector	FPD

1. Use **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to `Flow` mode, press **MODE/TYPE** to access the selection menu, then select `con pres`. Scroll to `Pressure` and set the pressure value to have the required carrier gas flow rate `Col.flow`.

RIGHT CARRIER ¹		
Pressure	30.0	30.0
Col.flow	3.00	
Lin. veloc.		(60.9)<

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		230
Final time 1		1.00<
Ramp 2		Off

3. Use **RIGHT INLET** to display the appropriate Programmable Temperature Vaporizing Injector Control Table. Set the required temperature setpoint

Temp. Verify to operate in **PTV splitless** mode. If not, scroll to Mode, press **MODE/TYPE** to access the selection menu, then select PTV splitless. Scroll to Splitless time to set the required setpoint.

RIGHT INLET (PTV)		
Temp	70	70
Pressure	30.0	30.0
Mode:	PTV Splitless	
Total flow	(53.0)	
Split Flow	50.0	50.0
Splitless time	0.80	0.80
Constant sept purge?		Y
Inject phase menu:		Y<

4. Scroll to Inject phase menu. Press **MODE/TYPE** to enter the **PTV Phase Menu**.

PTV PHASE MENU	
Ramped pressure?	N
Inject temp	50
Inject time	0.1
Transfer ramp	10
Transfer temp	260
Transfer time	1.00<

5. Select Ramped pressure? **NO**. Set the required Inject temp and *Inject time* setpoints as required. Then, set the Transfer ramp, the Transfer temp and the Transfer time required setpoints.
6. Use **RIGHT DETECTOR** to display the appropriate FPD Detector Control Table. Set the required temperatures Base Temp and FPD Temp. Then, set the detector gases H2 and Air required setpoints.

RIGHT DETECTOR (FPD)		
Flame		Off
Base temp	300	300
FPD temp	150	150
Signal pA		(1.4)
High voltage mode?		N
H2	90	90
Air	115	115
Mkup N2	00	00

7. Verify that High voltage mode is set to **NO**.
8. Scroll to Flame and press **ON**. This start the ignition sequence. When ignition is confirmed, the photomultiplier tube is energized. The baseline level Signal pA, will suddenly increase meaning that the flame is lit inside the detector. After a few seconds, the baseline should stabilize to the standing current of the system.
9. Use **RIGHT SIGNAL** to display the appropriate FPD Detector Signal Control Table. Scroll to Range and set the electrometer amplifier input range required.

RIGHT SIGNAL (FPD)	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Baseline comp	Off

10. Activate your Data System and set the parameters required for the checkout.
11. In the FPD Detector Signal Control Table scroll to Auto zero? and turn it **YES**.

12. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



Refer to the Acceptance Values reported in the Table 21-3 according to the data handling in use.

13. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
14. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
15. When the GC is ready, inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 21.1* or *21.2*.

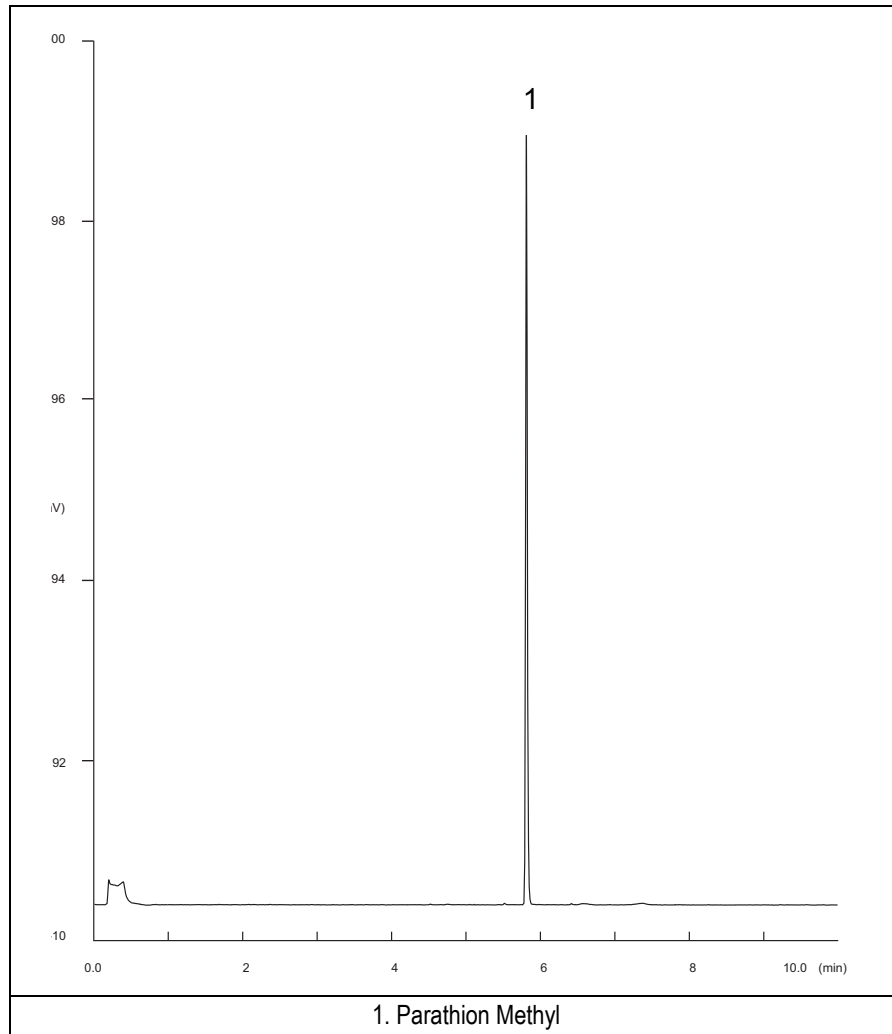


Figure 21-1. FPD-PTV Injection with 526 nm filter

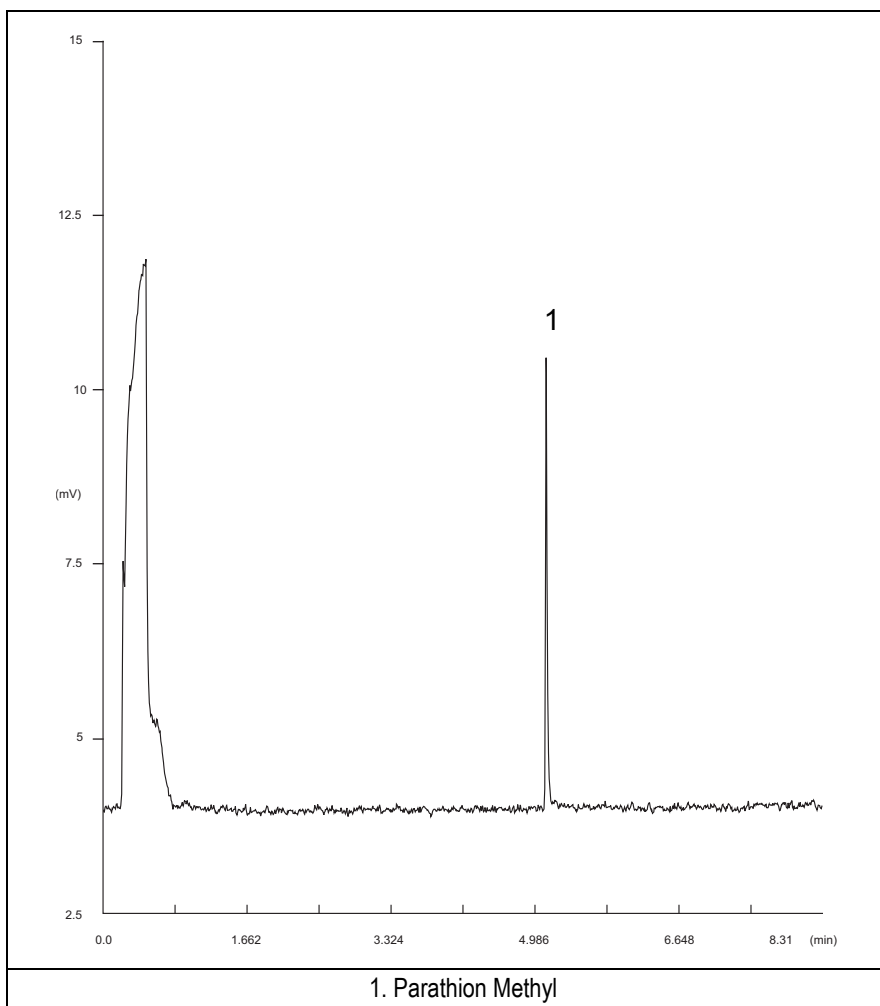


Figure 21-2. FPD-PTV Injection with 394 nm filter

16. Establish the integration parameters and the peak table identifying the test mix components.
17. Set up the data system to calculate the signal-to-noise ratio.

Using Chrom-Card Data System

Operate as follows:

- Open the Method Editor and include the signal-to-noise Report into the Report Parameters Page.
- By clicking on the side icon, open the signal-to-noise calculation and set Methylparathion component as signal peak ID.
- Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).
- Generate a report showing the chromatogram, peak area and signal-to-noise information for the component.



NOTE

If it is not possible to find a suitable part of baseline, to have a complete Chrom-Card final report proceed as follows: Start the acquisition on the data system. Let the system to acquire for about 1 minute and then inject the test mixture. The noise can be calculated during the first minute of acquisition. Since the retention time will shift by a delayed time, the Component Table has to be updated.

Using non-Chrom-Card Data System

Operate as follows:

- Set the parameters to calculate the signal-to-noise ratio according to the instruction reported in the relevant data system manual.




CAUTION


Choose a part of baseline without peaks or interference signals and calculate the noise for 0.1 min. (Verify that the noise value is comparable with the one of the previous *Baseline Acquisition and Analysis*).


18. The following criteria indicate successful completion of FPD-PTV checkout.


19. If these criteria are not met, repeat the test.

Table 21-3. FPD-PTV Acceptance Criteria

CHROM-CARD					
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)		Digital (10V Full Scale)	
		394 nm (S) Filter	526 nm (P) Filter	394 nm (S) Filter	526 nm (P) Filter
	Noise (μV)	< 50	< 50	< 500	< 500
	Wander (μV)	< 100	< 100	< 1 000	< 1 000
	Drift ($\mu\text{V/h}$)	< 100	< 100	< 1 000	< 1 000
Analytical Results					
Parathion Methyl Signal to Noise Ratio			394 nm (S) Filter	526 nm (P) Filter	
			> 20	> 1 000	
					

CHROMQUEST			
Acceptance Values	Baseline Parameters (1V Full Scale)		
		526 nm (P) Filter	
	Noise (μV)	< 50	< 50
	Wander (μV)	< 100	< 100
	Drift ($\mu\text{V/h}$)	< 100	< 100
Analytical Results			
Parathion Methyl Signal to Noise Ratio		394 nm (S) Filter	
		> 20	
			

ATLAS		
Baseline Parameters (10V Full Scale)		
		526 nm (P) Filter
Noise (μV)	< 500	< 500
Wander (μV)	< 1 000	< 1 000
Drift ($\mu\text{V/h}$)	< 1 000	< 1 000
Analytical Results		
Parathion Methyl Signal to Noise Ratio	394 nm (S) Filter	526 nm (P) Filter
	> 20	> 1 000
		

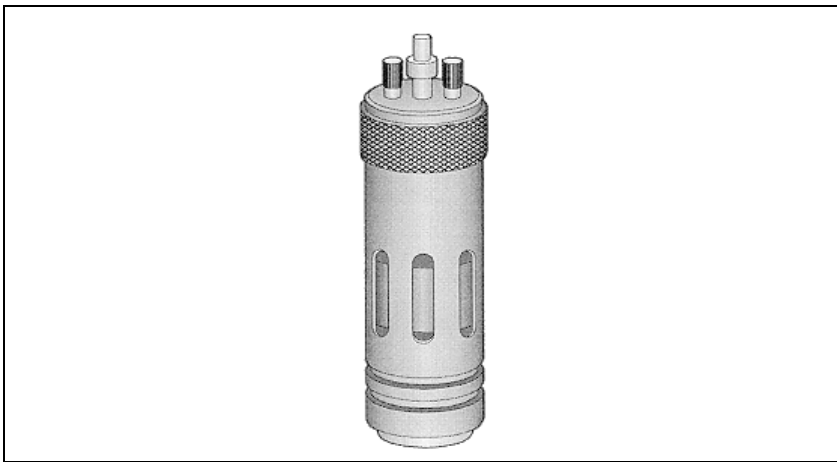
XCALIBUR			
Baseline Parameters (Acquisition Frequency = 10 Hz)			
Acceptance Values		526 nm (P) Filter	
	Noise (Counts)	< 5 000	< 5 000
	Wander (Counts)	< 10 000	< 10 000
	Drift (Counts/h)	< 10 000	< 10 000
	Analytical Results		
Parathion Methyl Signal to Noise Ratio	394 nm (S) Filter	526 nm (P) Filter	
	> 20	> 1 000	
			

Analytical Acceptance Comments	
1	Using Chrom-Card, set the signal-to-noise report parameters as described in the current procedure.
2	Using ChromQuest, Atlas, Xcalibur or a Computing integrator (e.g. ChromJet), calculate the S/N ratio as <i>Peak Height (counts)/noise (counts)</i> .

SECTION

VI

SOPs Using PID



The *SOPs Using PID* section, contains the procedures to test the TRACE GC Ultra with the Photoionization Detector (PID) using different injectors.

Chapter 22, *Checkout Using PID with S/SL Injector*.

Chapter 23, *Checkout Using PID with OC Injector*.

Chapter 24, *Checkout Using PID with PTV Injector*.

Checkout Using PID with S/SL Injector

Chapter at a Glance...

SOP Number: P0321/09/E - 01 September 2009	294
Scope	294
Parts Referenced	294
Analytical Conditions Required for Splitless Injection	296
Recommended Initial Operations	297

Operating Procedures

PID-S/SL Checkout in Splitless Mode	298
---	-----

SOP Number: P0321/09/E - 01 September 2009

Scope

Use the following procedure to verify proper PID operation with the Split/Splitless Injector.

Parts Referenced

Table 22-1. PID-S/SL Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long; 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	3 mm ID for splitless injection	453 200 32
Liner Seal	Graphite seal for glass liner	290 334 06
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
Septum	Standard septum for S/SL injector	313 032 11
Syringe	10 µl size; 70 mm needle length	365 001 03
Test Mixture	Three components in Iso-Octane: <i>Component</i> <i>Concentration</i> Azobenzene 1 µg/ml Octadecane 1000 µg/ml Parathion Methyl 1 µg/ml	338 190 06
Detector UV Lamp	8.4 eV 9.6 eV 10.6 eV 11.8 eV	305 030 13 305 030 14 305 030 15 305 030 16
Gases	Chromatographic-grade purity	

Table 22-1. PID-S/SL Parts Referenced (Continued)

Part	Description	Part Number
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for Splitless Injection

Table 22-2. PID-S/SL Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Make-up: Nitrogen = 7 ml/min Sheath Gas: Nitrogen = 40 ml/min
Oven Program	Initial Temperature = 70 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 230 °C Final Time = 1 minute
Injector	Operating Mode = Splitless Temperature = 230 °C Splitless Time = 0.8 minutes Split Flow = 60 ml/min Constant Septum Purge = Yes
Detector	Base Temperature = 230 °C High Current = No Detector Signal Range = 10 ⁰
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Replace the glass liner.
The glass liner currently installed in your injector should be carefully removed and replaced with the 3 mm ID glass liner for splitless application, as required for the checkout, with the appropriate graphite seal.
2. Replace the septum
A new septum should be installed properly in your injector.
3. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
4. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
5. Perform Column Evaluation and Leak Test
6. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

PID-S/SL Checkout in Splitless Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	S/SL
Left carrier or Right carrier	He (helium)
Left detector or Right detector	PID

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to Pressure and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		230
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate Split/Splitless Injector Control Table and set the required temperature setpoint *Temp*. Verify

to operate in **Splitless** mode. If not, scroll to *Mode*, press **MODE/TYPE** to access the selection menu, then select *Splitless*. Scroll to *Splitless time* to set the required setpoint.

LEFT INLET (S/SL) ¹		
Temp	230	230
Pressure	30.0	30.0
Mode:	Splitless	
Total flow	(63.0)	
Split Flow	60.0	60.0
Splitless time	0.80	0.80
Constant sept purge?	Y<	

1. These settings could also be for a right inlet.

- Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate PID Detector Control Table. Set the required temperature *Base Temp* and the detector gases *Mkup* and *Sheath Gas* required setpoints.

LEFT DETECTOR (PID) ¹		
Lamp	On	
Base temp	230	230
High current	N	
Signal pA	(15.4)	
Mkup N2	7.0	7.0
Sheath Gas	40	40

1. These settings could also be for a right detector.

- Verify that *High current* is set to **NO**.
- Scroll to *Lamp* and press **ON**. This start the ignition sequence. The baseline level *Signal pA*, will suddenly increase meaning that the lamp is lit inside the detector. A short period of conditioning is required in order to obtain a stable baseline.

7. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate PID Detector Signal Control Table. Scroll to `Range` and set the electrometer amplifier input range required.

LEFT SIGNAL (PID) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline comp	Off

1. These settings could also be for a right signal.

8. Activate your Data System and set the parameters required for the checkout.
9. In the PID Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
10. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 22-3 according to the data handling in use.

11. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
12. After the baseline the baseline evaluation has been completed, set-up the data system to acquire a single run.
13. Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatograms should look like the one shown in *Figures 22.1, 22.2, 22.3 or 22.4*.

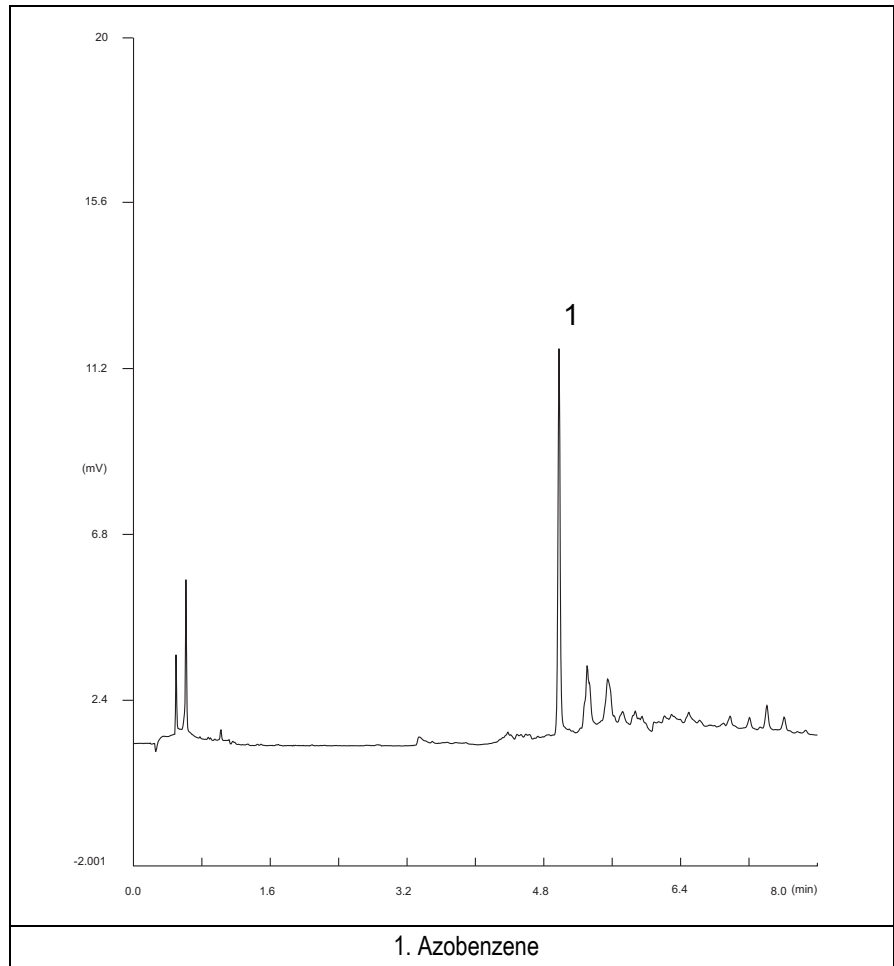


Figure 22-1. PID-Splitless Injection with 8.4 eV UV Lamp

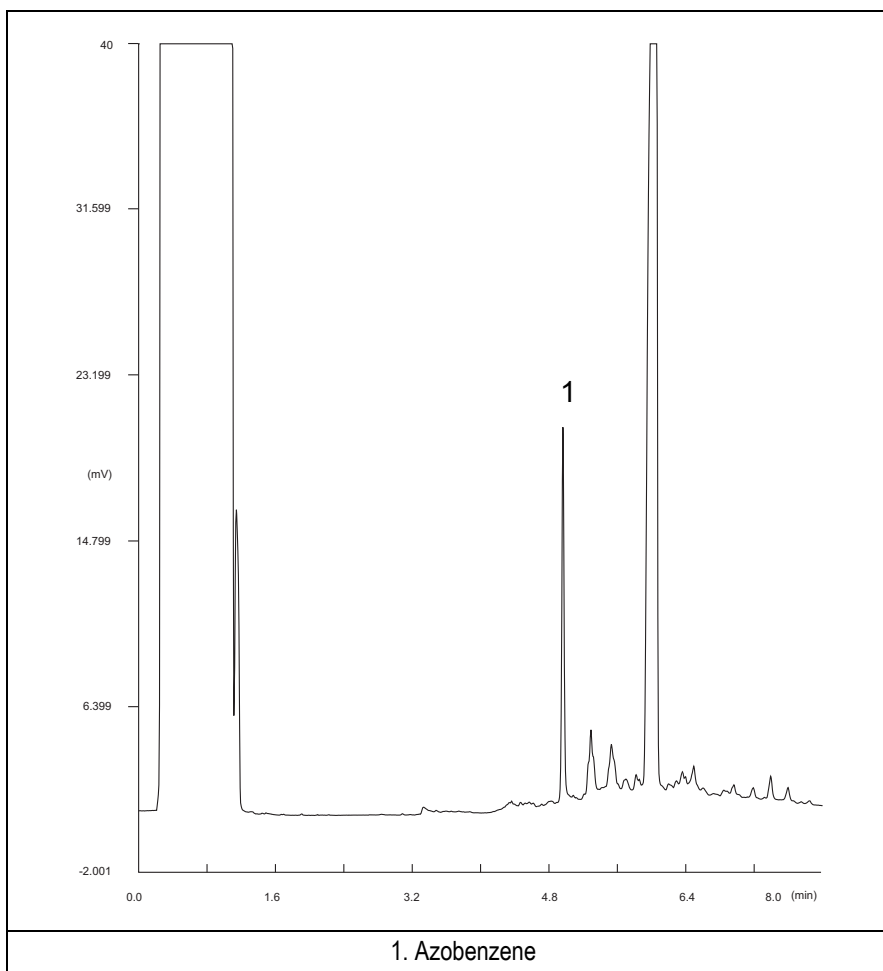


Figure 22-2. PID-Splitless Injection with 9.6 eV UV Lamp

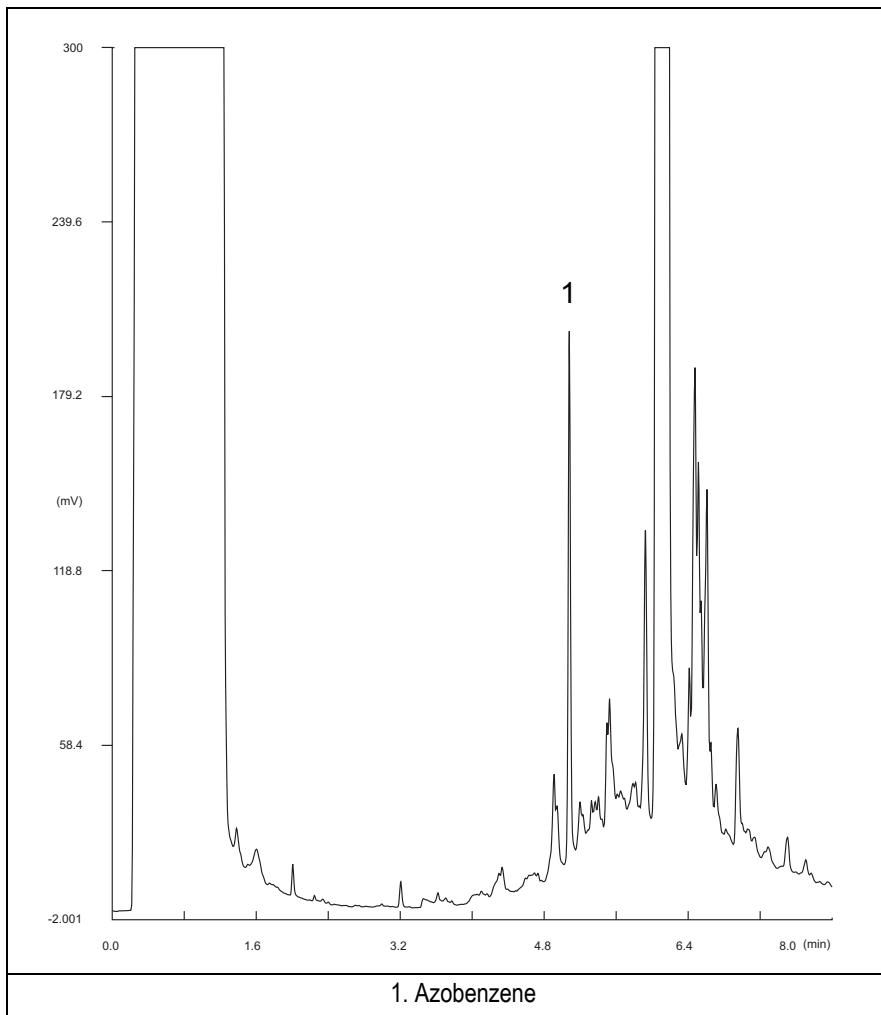


Figure 22-3. PID-Splitless Injection with 10.6 eV UV Lamp

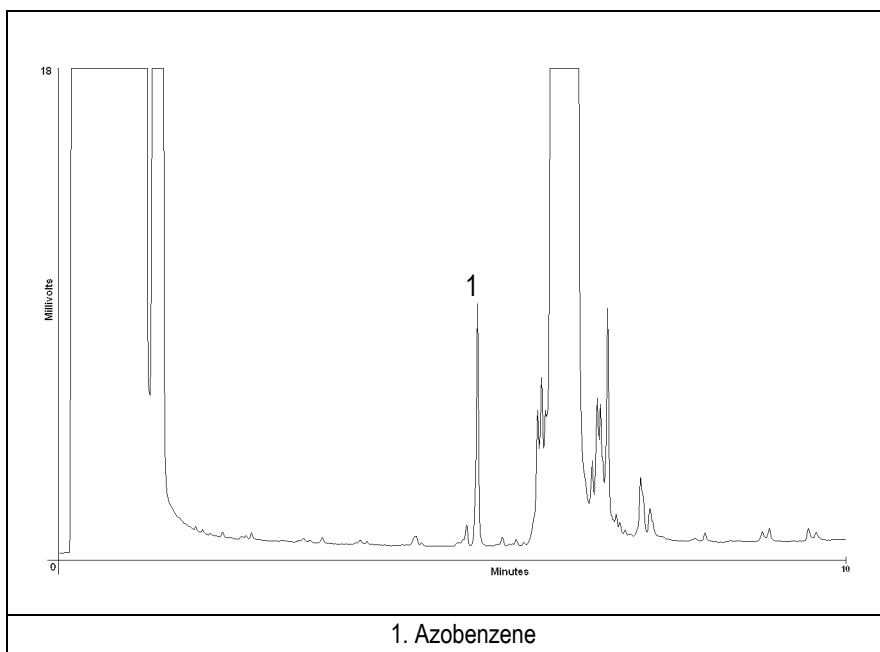





Figure 22-4. PID-Splitless Injection with 11.8 eV UV Lamp


14. The following criteria indicate successful completion of PID-S/SL checkout.
15. If these criteria are not met, repeat the test.


Table 22-3. PID-S/SL Acceptance Criteria

CHROM-CARD										
Acceptance Values	Baseline Parameters		Analog (1V Full Scale)				Digital (10V Full Scale)			
	Lamp (eV)	8.4	9.6	10.6	11.8	8.4	9.6	10.6	11.8	
	Noise (µV)	< 100	< 50	< 150	< 150	< 1 000	< 500	< 1 500	< 1 500	
	Wander (µV)	< 150	< 100	< 100	< 100	< 1 500	< 1 000	< 1 000	< 1 000	
	Drift (µV/h)	< 100	< 100	< 100	< 100	< 1 000	< 1 000	< 1 000	< 1 000	
	Analytical Results		Analog (1V Full Scale) Area Counts (0.1 µVs)				Digital (10V Full Scale) Area Counts (0.1 µVs)			
	Lamp (eV)	8.4	9.6	10.6	11.8	8.4	9.6	10.6	11.8	
	Azobenzene	> 100*	> 160*	> 2 000*	> 40*	> 1 000*	> 1 600*	> 20 000*	> 400*	
										

Computing-integrator (e.g. ChromJet)									
									

CHROMQUEST									
Acceptance Values	Baseline Parameters (1V Full Scale)								
	Lamp (eV)	8.4	9.6	10.6	11.8				
	Noise (µV)	< 100	< 50	< 150	< 150				
	Wander (µV)	< 150	< 100	< 100	< 100				
	Drift (µV/h)	< 100	< 100	< 100	< 100				
	Analytical Results (1V Full Scale) - Area Counts (0.01 µVs)								
	Lamp (eV)	8.4	9.6	10.6	11.8				
	Azobenzene	> 1 000 000	> 1 600 000	> 20 000 000	> 400 000				
									

ATLAS				
Baseline Parameters (10V Full Scale)				
Lamp (eV)	8.4	9.6	10.6	11.8
Noise (μV)	< 1 000	< 500	< 1 500	< 1 500
Wander (μV)	< 1 500	< 1 000	< 1 000	< 1 000
Drift ($\mu\text{V}/\text{h}$)	< 1 000	< 1 000	< 1 000	< 1 000
Analytical Results (10V Full Scale) - Area Counts (μVs)				
Lamp (eV)	8.4	9.6	10.6	11.8
Azobenzene	> 100 000	> 160 000	> 2 000 000	> 40 000
				

XCALIBUR				
Baseline Parameters (Acquisition Frequency = 10 Hz)				
Lamp (eV)	8.4	9.6	10.6	11.8
Noise (Counts)	< 10 000	< 5 000	< 15 000	< 15 000
Wander Counts)	< 15 000	< 10 000	< 10 000	< 10 000
Drift (Counts/h)	< 10 000	< 10 000	< 10 000	< 10 000
Analytical Results Area Counts (Cts*s)				
Lamp (eV)	8.4	9.6	10.6	11.8
Azobenzene	> 1 000 000	> 1 600 000	> 20 000 000	> 400 000
				

Checkout Using PID with OC Injector

Chapter at a Glance...

SOP Number: P0322/09/E - 01 September 2009	308
Scope	308
Parts Referenced	308
Analytical Conditions Required for On-Column Injection	309
Recommended Initial Operations	310

Operating Procedures

PID-OCI Checkout in On-Column Mode.....	311
---	-----

SOP Number: P0322/09/E - 01 September 2009

Scope

Use the following procedure to verify proper PID operation with the On-Column Injector.

Parts Referenced

Table 23-1. PID-OCI Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long; 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Graphite Ferrule	Graphite Ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
Vespel Ferrule	Vespel Ferrule for 0.32 mm ID Column	290 134 60
Syringe	10 µl size; 75 mm needle length	365 020 07
Test Mixture	Three components in Iso-Octane: <i>Component</i> <i>Concentration</i> Azobenzene 1 µg/ml Octadecane 1000 µg/ml Parathion methyl 1 µg/ml	338 190 06
Detector UV Lamp	8.4 eV 9.6 eV 10.6 eV 11.8 eV	305 030 13 305 030 14 305 030 15 305 030 16
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Table 23-1. PID-OCI Parts Referenced (Continued)

Part	Description	Part Number
In case of automatic On-Column for TriPlus Sampler AS		
Syringe	10 µl size; 80 mm needle length	365 020 19
Pre-column	2 m long; 0.53 mm ID	260 603 75
Press-fit set	Set of Press-fit connectors for TRACE OC	350 038 45

Analytical Conditions Required for On-Column Injection

Table 23-2. PID-OCI Analytical Conditions

Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Make-up: Nitrogen = 7 ml/min Sheath Gas: Nitrogen = 40 ml/min
Oven Program	Initial Temperature = 85 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 230 °C Final Time = 1 minute
Injector	Secondary cooling = 0.2 minutes
Detector	Base Temperature = 230 °C High Current = No Detector Signal Range = 10 ⁰
Injected Volume	1 µl of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
2. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
In case of automatic On-column for TriPlus sampler, install the pre-column and connect it to the test column by press-fit connector.
3. Install and connect the TriPlus sampler and its components.
4. Perform Column Evaluation and Leak Test.
5. Connect your data handling.
Verify that your data handling is properly connected to your GC system.
6. Verify the opening/closing of the OC injector actuator by using the proper commands.
7. Verify the alignment of the syringe on the OC injector.

OPERATING PROCEDURE

PID-OCI Checkout in On-Column Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	OCI
Left carrier or Right carrier	He (helium)
Left detector or Right detector	PID

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to Pressure and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	85.0	85.0
Initial Time		1.00
Ramp 1		20.0
Final temp		230
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate On-Column Injector Control Table. Scroll to `Sec. cool time` and set the required secondary cooling time.

LEFT INLET (OCI) ¹		
Pressure	30.0	30.0
Sec. Cool Time		10.00<

1. These settings could also be for a right inlet.

4. Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate PID Detector Control Table. Set the required temperatures `Base Temp` and the detector gases `Mkup` and `Sheath Gas` required setpoints.

LEFT DETECTOR (PID) ¹		
Lamp		On
Base temp	230	230
High current		N
Signal pA		(15.4)
Mkup N2	7.0	7.0
Sheath Gas	40	40

1. These settings could also be for a right detector.

5. Verify that `High current` is set to **NO**.
6. Scroll to `Lamp` and press **ON**. This start the ignition sequence. The baseline level `Signal pA`, will suddenly increase meaning that the lamp is lit inside the detector. A short period of conditioning is required in order to obtain a stable baseline.
7. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate PID Detector Signal Control Table. Scroll to `Range` and set the electrometer amplifier input range required.

LEFT SIGNAL (PID) ¹	
Output	(1000)
Offset	100
Autozero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline comp	Off

1. These settings could also be for a right signal.

8. Activate your Data System and set the parameters required for the checkout.
9. In the PID Detector Signal Control Table scroll to *Auto zero?* and turn it **YES**.
10. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 23-3 according to the data handling in use.

11. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
12. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
13. Perform the analysis.

Manual injection

- Inject the test mixture and press **START** on the GC to begin the checkout run.

Automatic injection with TriPlus sampler

- Fill a vial with the standard mix and place that vial in the sample tray.

- Load the method for OC and perform the sampling.

The resulting chromatograms should look like the one shown in *Figures 23.1, 23.2, 23.3 or 23.4.*

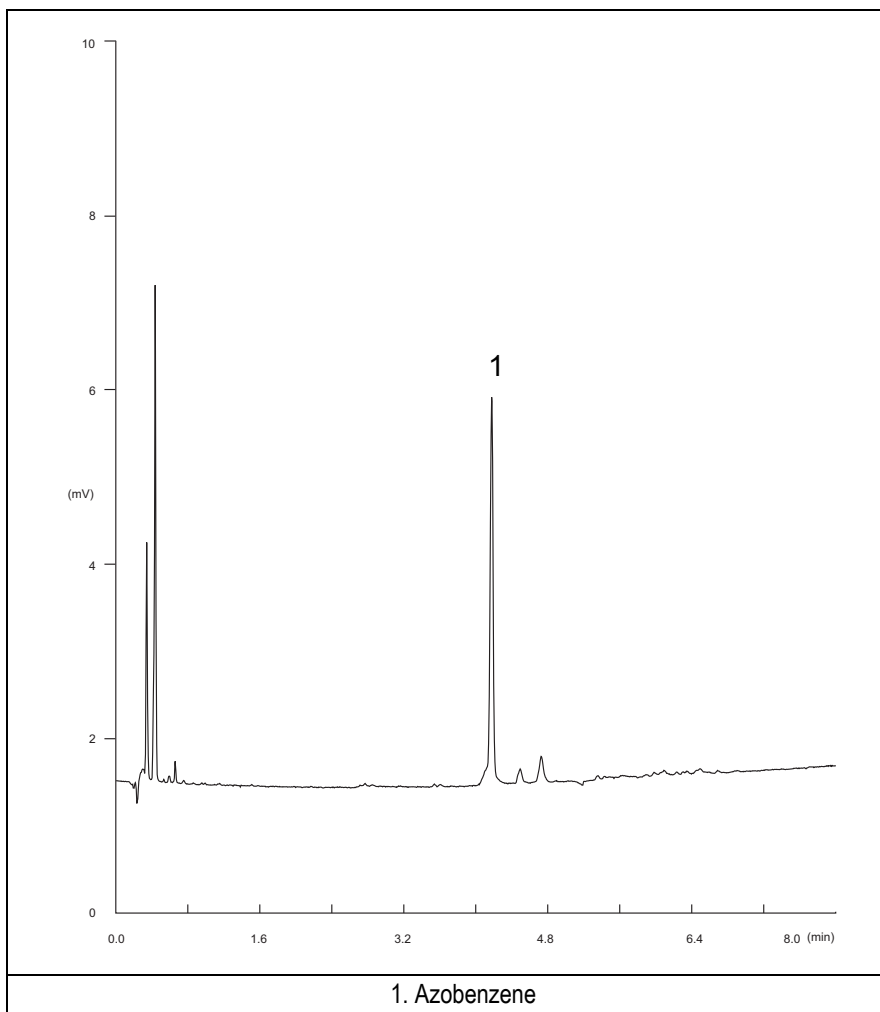


Figure 23-1. PID-On-Column Injection with 8.4 eV UV Lamp

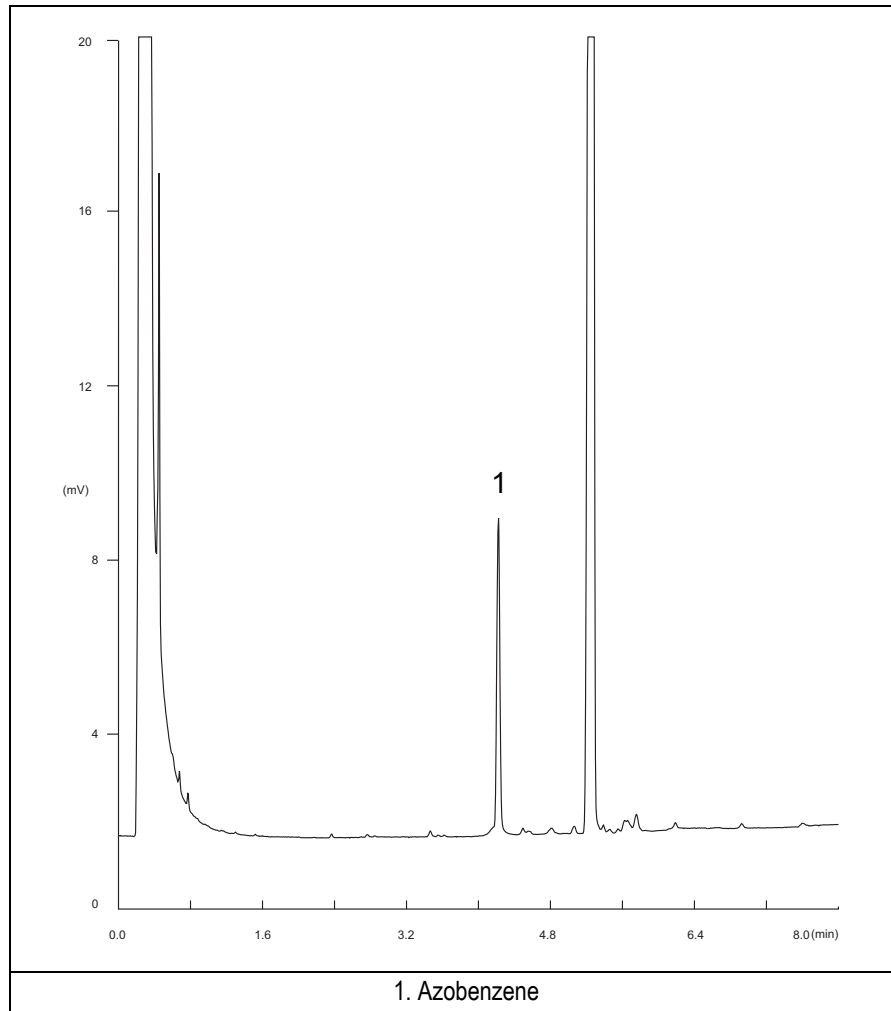


Figure 23-2. PID-On-Column Injection with 9.6 eV UV Lamp

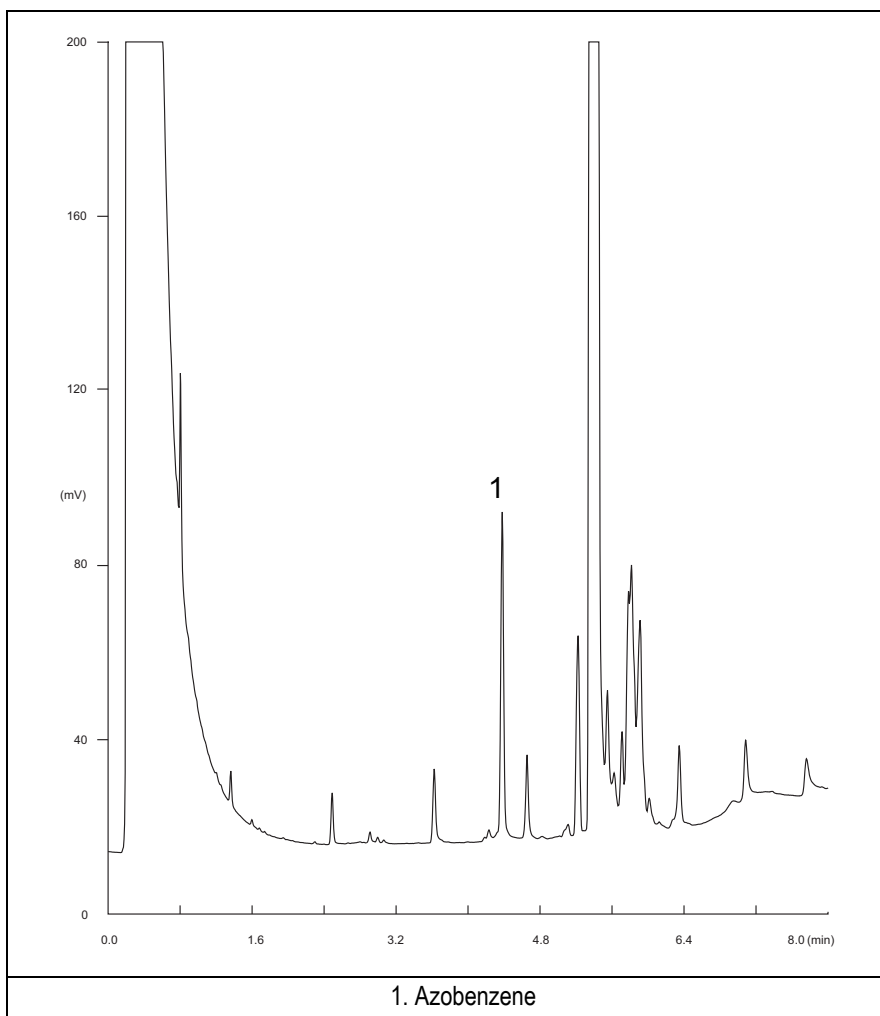


Figure 23-3. PID-On-Column Injection with 10.6 eV UV Lamp

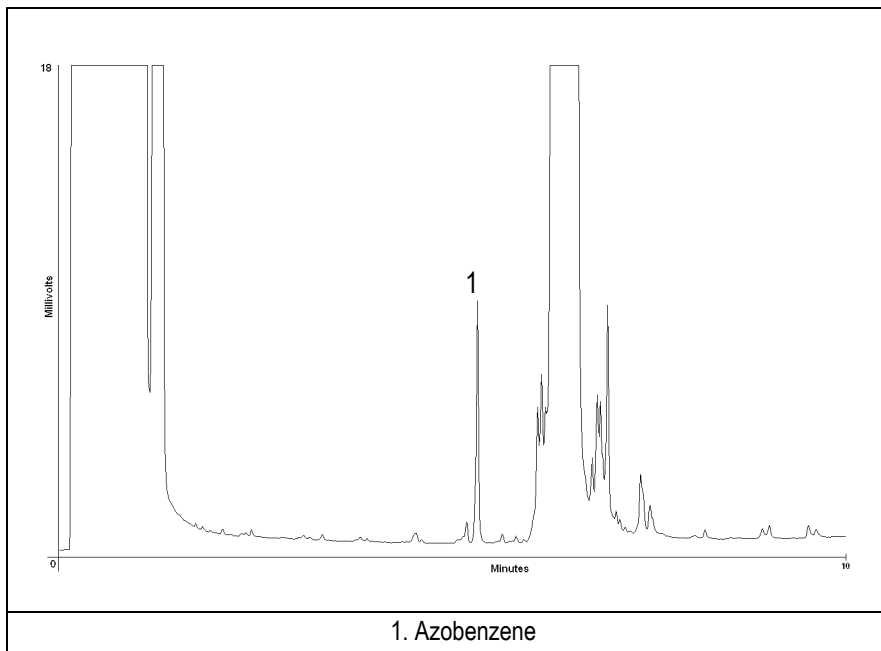


Figure 23-4. PID-On-Column Injection with 11.8 eV UV Lamp


- 14. The following criteria indicate successful completion of PID-OCI checkout.
- 15. If these criteria are not met, repeat the test.


Table 23-3. PID-OCI Acceptance Criteria

CHROM-CARD									
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)				Digital (10V Full Scale)			
	Lamp (eV)	8.4	9.6	10.6	11.8	8.4	9.6	10.6	11.8
	Noise (μV)	< 100	< 50	< 150	< 150	< 1 000	< 500	< 1 500	< 1 500
	Wander (μV)	< 150	< 100	< 100	< 100	< 1 500	< 1 000	< 1 000	< 1 000
	Drift (μV/h)	< 100	< 100	< 100	< 100	< 1 000	< 1 000	< 1 000	< 1 000
	Analytical Results	Analog (1V Full Scale) Area Counts (0.1 μVs)				Digital (10V Full Scale) Area Counts (0.1 μVs)			
	Lamp (eV)	8.4	9.6	10.6	11.8	8.4	9.6	10.6	11.8
	Azobenzene	> 60*	> 110*	> 1 300*	> 24*	> 600*	> 1 100*	> 13 000*	> 240*
	⚠								

Computing-integrator (e.g. ChromJet)	
⚠	

CHROMQUEST					
Acceptance Values	Baseline Parameters (1V Full Scale)				
	Lamp (eV)	8.4	9.6	10.6	11.8
	Noise (μV)	< 100	< 50	< 150	< 150
	Wander (μV)	< 150	< 100	< 100	< 100
	Drift (μV/h)	< 100	< 100	< 100	< 100
	Analytical Results (1V Full Scale) - Area Counts (0.01 μVs)				
	Lamp (eV)	8.4	9.6	10.6	11.8
	Azobenzene	> 600 000	> 1 100 000	> 13 000 000	> 240 000
	⚠				

ATLAS				
Baseline Parameters (10V Full Scale)				
Lamp (eV)	8.4	9.6	10.6	11.8
Noise (μV)	< 1 000	< 500	< 1 500	< 1 500
Wander (μV)	< 1 500	< 1 000	< 1 000	< 1 000
Drift ($\mu\text{V}/\text{h}$)	< 1 000	< 1 000	< 1 000	< 1 000
Analytical Results (10V Full Scale) - Area Counts (μVs)				
Lamp (eV)	8.4	9.6	10.6	11.8
Azobenzene	> 60 000	> 110 000	> 1 300 000	> 24 000
				

XCALIBUR				
Baseline Parameters (Acquisition Frequency = 10 Hz)				
Lamp (eV)	8.4	9.6	10.6	11.8
Noise (Counts)	< 10 000	< 5 000	< 15 000	< 15 000
Wander Counts)	< 15 000	< 10 000	< 10 000	< 10 000
Drift (Counts/h)	< 10 000	< 10 000	< 10 000	< 10 000
Analytical Results Area Counts (Cts*s)				
Lamp (eV)	8.4	9.6	10.6	11.8
Azobenzene	> 600 000	> 1 100 000	> 13 000 000	> 240 000
				

Checkout Using PID with PTV Injector

Chapter at a Glance...

SOP Number: P0325/08/E - 01 September 2009	322
Scope	322
Parts Referenced	322
Analytical Conditions Required for PTV Splitless Injection	324
Recommended Initial Operations	325

Operating Procedures

PID-PTV Checkout in PTV Splitless Mode	326
--	-----

SOP Number: P0325/08/E - 01 September 2009

Scope

Use the following procedure to verify proper PID operation with the Programmable Temperature Vaporizing Injector.

Parts Referenced

Table 24-1. PID-PTV Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Liner	Silcosteel 2 mm ID (set of 2)	453 220 44
Liner Seal	Graphite seal for liner	290 034 17
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
Septum	Standard septum for PTV injector (set of 10)	313 132 25
Syringe	10 µl size; 50 mm needle length	365 005 25
Test Mixture	Three components in Iso-Octane: <i>Component</i> <i>Concentration</i> Azobenzene 1 µg/ml Octadecane 1000 µg/ml Parathion methyl 1 µg/ml	338 190 06
Detector UV Lamp	8.4 eV 9.6 eV 10.6 eV 11.8 eV	305 030 13 305 030 14 305 030 15 305 030 16
Gases	Chromatographic-grade purity	

Table 24-1. PID-PTV Parts Referenced (Continued)

Part	Description	Part Number
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for PTV Splitless Injection

Table 24-2. PID-PTV Analytical Conditions

Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Make-up: Nitrogen = 7 ml/min Sheath Gas: Nitrogen = 40 ml/min
Oven Program	Initial Temperature = 70 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 230 °C Final Time = 1 minute
Injector	Operating Mode = PTV splitless Splitless Time = 0.8 minutes Split Flow = 50 ml/min Constant Septum Purge = Yes Inject Temp = 50 °C Inject Time = 0.1 minute Transfer ramp = 10 °C/sec Transfer Temperature = 260 °C Transfer time = 1 minute
Detector	Base Temperature = 230 °C High Current = No Detector Signal Range = 10 ⁰
Injected Volume	1 µl of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Replace the liner.
The liner currently installed in your injector should be carefully removed and replaced with the 2 mm ID Silcosteel liner, as required for the checkout, with the appropriate liner seal.
2. Replace the septum
A new septum should be installed properly in your injector.
3. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
4. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
5. Perform Column Evaluation and Leak Test.
6. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

PID-PTV Checkout in PTV Splitless Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Right inlet	PTV
Right carrier	He (helium)
Right detector	PID

1. Use **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to `Flow mode`, press **MODE/TYPE** to access the selection menu, then select `con pres`. Scroll to `Pressure` and set the pressure value to have the required carrier gas flow rate `Col.flow`.

RIGHT CARRIER ¹		
Pressure	30.0	30.0
Col.flow	3.00	
Lin. veloc.		(60.9)<

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		230
Final time 1		1.00<
Ramp 2		Off

3. Use **RIGHT INLET** to display the appropriate Programmable Temperature Vaporizing Injector Control Table. Set the required temperature setpoint

Temp. Verify to operate in **PTV splitless** mode. If not, scroll to Mode, press **MODE/TYPE** to access the selection menu, then select PTV splitless. Scroll to Splitless time to set the required setpoint.

RIGHT INLET (PTV)		
Temp	70	70
Pressure	30.0	30.0
Mode:	PTV Splitless	
Total flow	(53.0)	
Split Flow	50.0	50.0
Splitless time	0.80	0.80
Constant sept purge?		Y
Inject phase menu:		Y<

4. Scroll to Inject phase menu. Press **MODE/TYPE** to enter the **PTV Phase Menu**.

PTV PHASE MENU	
Ramped pressure?	N
Inject temp	50
Inject time	0.1
Transfer ramp	10
Transfer temp	260
Transfer time	1.00<

5. Select Ramped pressure? **NO**. Set the required Inject temp and *Inject time* setpoints as required. Then, set the Transfer ramp, the Transfer temp and the Transfer time required setpoints.
6. Use **RIGHT DETECTOR** to display the appropriate PID Detector Control Table. Set the required temperature Base Temp and the detector gases Mkup and Sheath Gas required setpoints.

RIGHT DETECTOR (PID)		
Lamp		On
Base temp	230	230
Lamp current		low
Signal pA		(15.4)
Mkup N2	7.0	7.0
Sheath Gas	40	40

7. Verify that High current is set to **NO**.
8. Scroll to Lamp and press **ON**. This start the ignition sequence. The baseline level Signal pA, will suddenly increase meaning that the lamp is lit inside the detector. A short period of conditioning is required in order to obtain a stable baseline.
9. Use **RIGHT SIGNAL** to display the appropriate PID Detector Signal Control Table. Scroll to *Range* and set the electrometer amplifier input range required.

RIGHT SIGNAL (PID)	
Output	(1000)
Offset	100
Autozero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Analog filter	Off

10. Activate your Data System and set the parameters required for the checkout.
11. In the PID Detector Signal Control Table scroll to Auto zero? and turn it **YES**.
12. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

**NOTE****Baseline Acquisition and Analysis**

Refer to the Acceptance Values reported in the Table 24-3 according to the data handling in use.

13. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
14. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
15. Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatograms should look like the one shown in *Figure 24.1, 24.2, 24.3 or 24.4*.

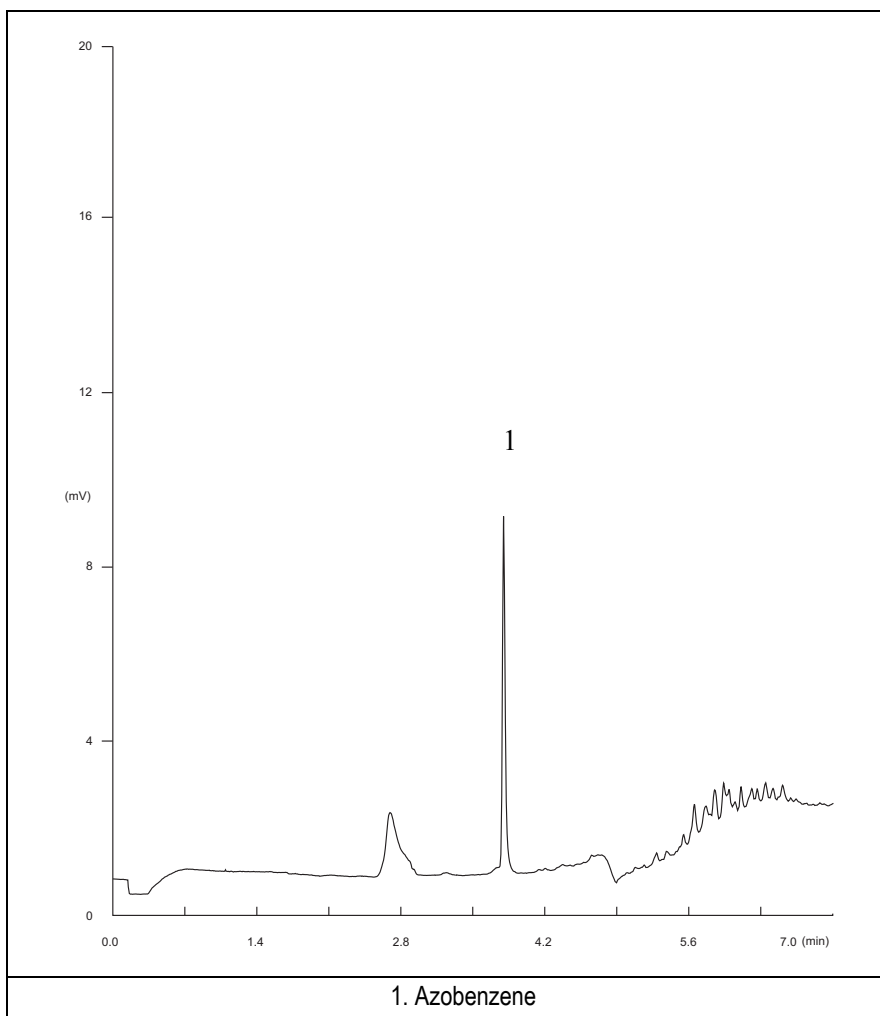


Figure 24-1. PID-PTV Injection with 8.4 eV UV Lamp

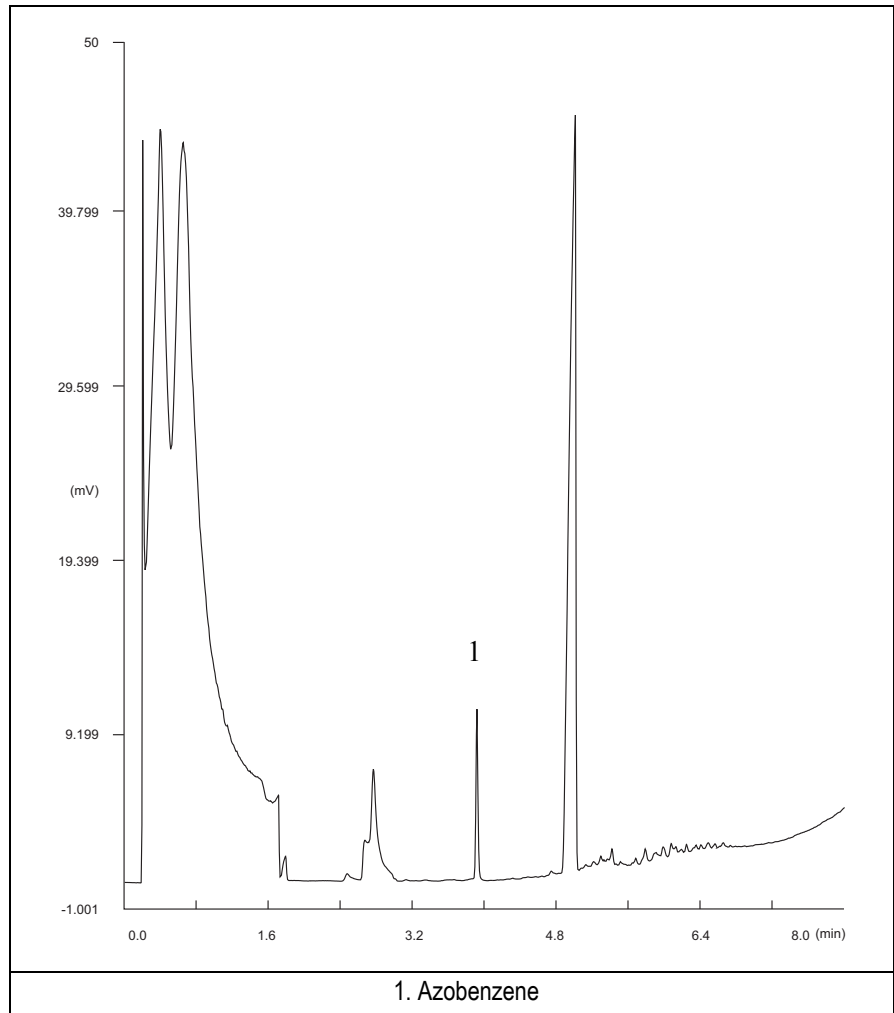


Figure 24-2. PID-PTV Injection with 9.6 eV UV Lamp

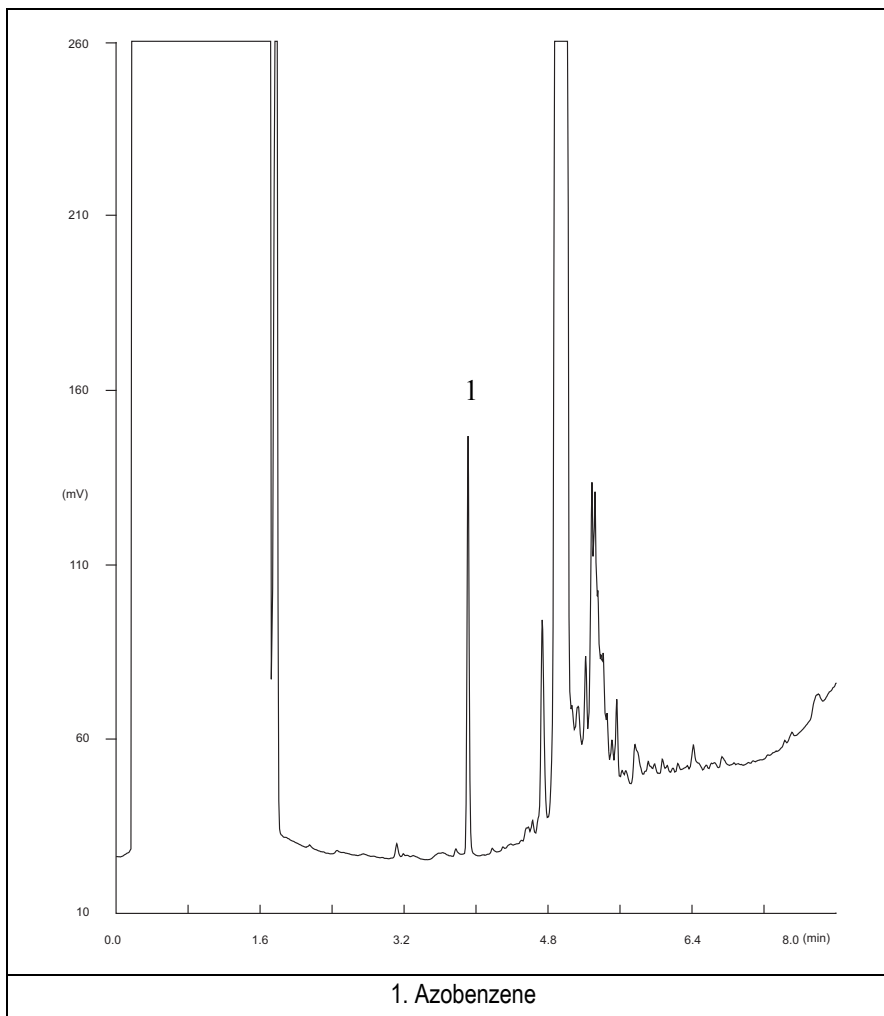


Figure 24-3. PID-PTV Injection with 10.6 eV UV Lamp

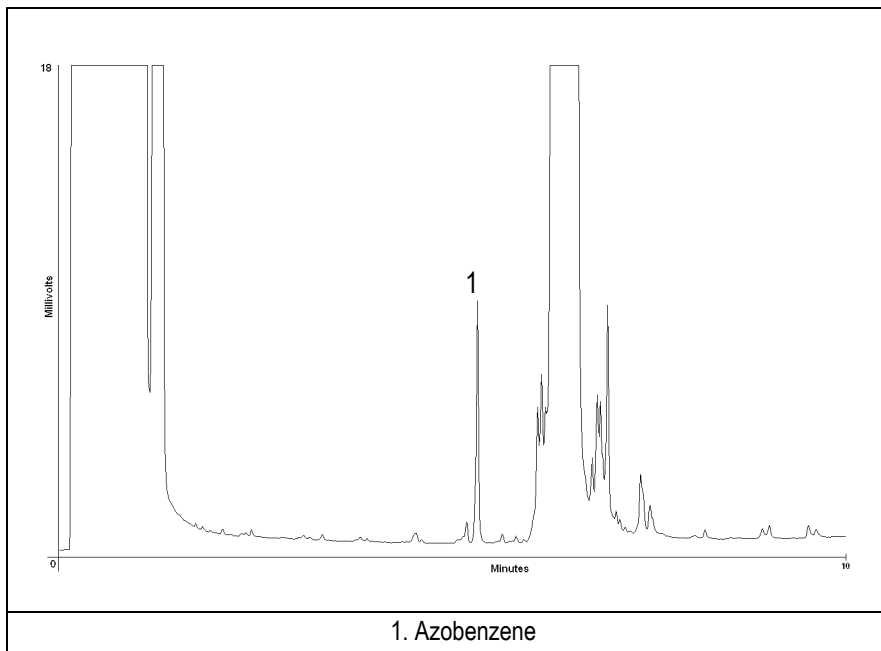





Figure 24-4. PID-PTV Injection with 11.8 eV UV Lamp


- 16. The following criteria indicate successful completion of PID-PTV checkout.
- 17. If these criteria are not met, repeat the test.


Table 24-3. PID-PTV Acceptance Criteria

CHROM-CARD									
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)				Digital (10V Full Scale)			
	Lamp (eV)	8.4	9.6	10.6	11.8	8.4	9.6	10.6	11.8
	Noise (μV)	< 100	< 50	< 150	< 150	< 1 000	< 500	< 1 500	< 1 500
	Wander (μV)	< 150	< 100	< 100	< 100	< 1 500	< 1 000	< 1 000	< 1 000
	Drift (μV/h)	< 100	< 100	< 100	< 100	< 1 000	< 1 000	< 1 000	< 1 000
	Analytical Results	Analog (1V Full Scale) Area Counts (0.1 μVs)				Digital (10V Full Scale) Area Counts (0.1 μVs)			
	Lamp (eV)	8.4	9.6	10.6	11.8	8.4	9.6	10.6	11.8
	Azobenzene	> 60*	> 110*	> 1 300*	> 24*	> 600*	> 1 100*	> 13 000*	> 240*
									

Computing-integrator (e.g. ChromJet)	
	

CHROMQUEST					
Acceptance Values	Baseline Parameters (1V Full Scale)				
	Lamp (eV)	8.4	9.6	10.6	11.8
	Noise (μV)	< 100	< 50	< 150	< 150
	Wander (μV)	< 150	< 100	< 100	< 100
	Drift (μV/h)	< 100	< 100	< 100	< 100
	Analytical Results (1V Full Scale) - Area Counts (0.01 μVs)				
	Lamp (eV)	8.4	9.6	10.6	11.8
	Azobenzene	> 600 000	> 1 100 000	> 13 000 000	> 240 000
					

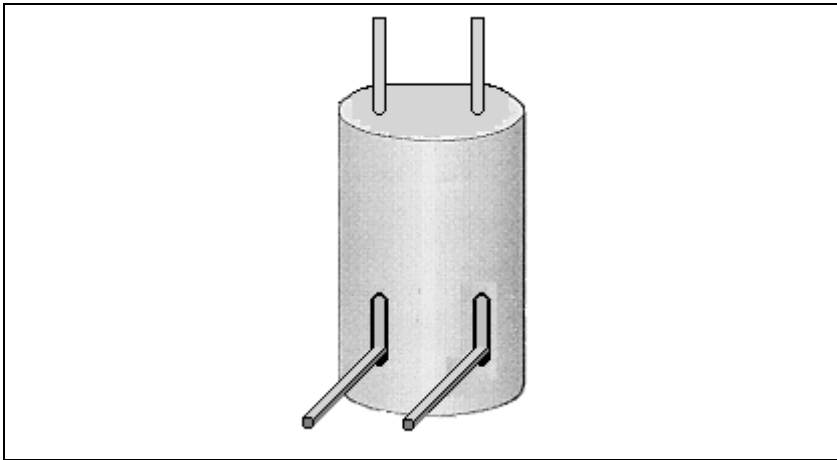
ATLAS				
Baseline Parameters (10V Full Scale)				
Lamp (eV)	8.4	9.6	10.6	11.8
Noise (μV)	< 1 000	< 500	< 1 500	< 1 500
Wander (μV)	< 1 500	< 1 000	< 1 000	< 1 000
Drift ($\mu\text{V/h}$)	< 1 000	< 1 000	< 1 000	< 1 000
Analytical Results (10V Full Scale) - Area Counts (μVs)				
Lamp (eV)	8.4	9.6	10.6	11.8
Azobenzene	> 60 000	> 110 000	> 1 300 000	> 24 000
				

XCALIBUR				
Baseline Parameters (Acquisition Frequency = 10 Hz)				
Lamp (eV)	8.4	9.6	10.6	11.8
Noise (Counts)	< 10 000	< 5 000	< 15 000	< 15 000
Wander Counts)	< 15 000	< 10 000	< 10 000	< 10 000
Drift (Counts/h)	< 10 000	< 10 000	< 10 000	< 10 000
Analytical Results Area Counts (Cts*s)				
Lamp (eV)	8.4	9.6	10.6	11.8
Azobenzene	> 600 000	> 1 100 000	> 13 000 000	> 240 000
				

SECTION

VII

SOPs Using TCD



The *SOPs Using TCD* section, contains the procedures to test the TRACE GC Ultra with the Thermal Conductivity Detector (TCD) using different injectors.

Chapter 25, *Checkout Using TCD with S/SL Injector*.

Chapter 26, *Checkout Using TCD with PKD Injector*.

Chapter 27, *Checkout Using TCD with PPKD Injector*.

Checkout Using TCD with S/SL Injector

Chapter at a Glance...

SOP Number: P0326/11/E - 01 September 2009	340
Scope	340
Parts Referenced	340
Analytical Conditions Required for Splitless Injection	341
Recommended Initial Operations	342

Operating Procedures

TCD-S/SL Checkout in Splitless Mode	343
---	-----

SOP Number: P0326/11/E - 01 September 2009

Scope

Use the following procedure to verify proper TCD operation with the Split/Splitless Injector.

Parts Referenced

Table 25-1. TCD-S/SL Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long; 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	3 mm ID for splitless injection	453 200 32
Liner Seal	Graphite seal for glass liner	290 334 06
Retaining Nut	M4 capillary column retaining nut	350 324 23
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Septum	Standard septum for S/SL injector	313 032 11
Syringe	10 µl size; 70 mm needle length	365 001 03
Test Mixture	Three components in n-Hexane: <i>Component</i> <i>Concentration</i> Dodecane 200 µg/ml Tetradecane 200 µg/ml Hexadecane 200 µg/ml	338 190 16
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for Splitless Injection

Table 25-2. TCD-S/SL Analytical Conditions

Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Make-up: Helium = 27.5 ml/min Reference Gas: Helium = 30 ml/min
Oven Program	Initial Temperature = 50 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 190 °C Final Time = 1 minute
Injector	Operating Mode = Splitless Temperature = 200 °C Splitless Time = 0.8 minutes Split Flow = 60 ml/min Constant Septum Purge = Yes
Detector	Block Temperature = 200 °C Transfer Temperature = 190 °C Constant Filament Temperature = No Filament Voltage = 10V Filament Temperature limit = 350 °C (*) Gain = x 10 Negative Polarity = No
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

(*) In case of TCD with the polyimide coated filaments, set the filament temperature limit to 320 °C.

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Replace the glass liner.
The glass liner currently installed in your injector should be carefully removed and replaced with the 3 mm ID glass liner for splitless application, as required for the checkout, with the appropriate graphite seal.
2. Replace the septum
A new septum should be installed properly in your injector.
3. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
4. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
5. Perform Column Evaluation and Leak Test
6. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

TCD-S/SL Checkout in Splitless Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	S/SL
Left carrier or Right carrier	He (helium)
Left detector or Right detector	TCD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to Pressure and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	50.0	50.0
Initial Time		1.00
Ramp 1		20.0
Final temp		190
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate Split/Splitless Injector Control Table and set the required temperature setpoint *Temp*. Verify

to operate in **Splitless** mode. If not, scroll to Mode, press **MODE/TYPE** to access the selection menu, then select Splitless. Scroll to Splitless time to set the required setpoint.

LEFT INLET (S/SL) ¹		
Temp	200	200
Pressure	30.0	30.0
Mode:	Splitless	
Total flow	(63.0)	
Split Flow	60.0	60.0
Splitless time	0.80	0.80
Constant sept purge?	Y<	

1. These settings could also be for a right inlet.

- Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate TCD Detector Control Table. Sequentially scroll to Block Temp, Transf Temp, Ref flow and mkup flow and set the required values.

LEFT DETECTOR (TCD) ¹		
Filament power	Off	
Fil status	(not rdy)	
Block temp	200	200
Transf temp	190	190
Const fil temp	Y/N	
Fil volts (CV)	10	
Fil temp limit ²	350	
Ref flow	30.0	30.0
Mkup flow	27.5	27.5<

1. These settings could also be for a right detector.

2. In case of TCD with the polyimide coated filaments, the temperature limit is 320 °C

- Scroll to Const fil temp and select it **NO**.
- Scroll to Fil volts (CV) and set the filament voltage. Scroll to Fil temp limit and set the required limit temperature setpoint.

7. Scroll to `Filament power` and turn it **ON**. After a few seconds the `ready` is displayed on `Fil status` line.
8. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate TCD Detector Signal Control Table. Scroll to `Gain` and set the desired value.

LEFT SIGNAL (TCD) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Gain (x1..x10)	10<
Neg polarity?	N
Baseline comp	Off

1. These settings could also be for a right signal.

9. Activate your Data System and set the parameters required for the checkout.
10. In the TCD Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
11. Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 25-3 according to the data handling in use.

12. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
13. After the baseline the baseline evaluation has been completed, set-up the data system to acquire a single run.
14. Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 25.1*.

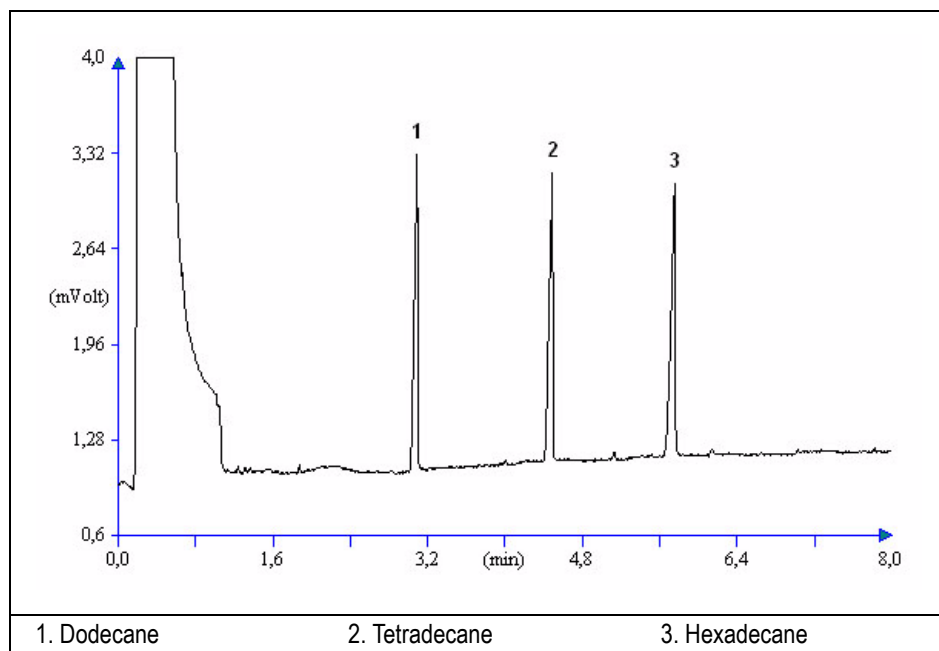





Figure 25-1. TCD-Splitless Injection


15. The following criteria indicate successful completion of TCD-S/SL checkout.
16. If these criteria are not met, repeat the test.


Table 25-3. TCD-S/SL Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 30	< 300
	Wander (μV)	< 140	< 1 400
	Drift ($\mu\text{V/h}$)	< 200	< 2 000
	Analytical Results	Analog (1V Full Scale) Area Counts (0.1 μVs)	Digital (10V Full Scale) Area Counts (0.1 μVs)
	Components	> 33 000 for each component	> 330 000 for each component
			

Computing-integrator (e.g. ChromJet)	
	

CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 30
	Wander (μV)	< 140
	Drift ($\mu\text{V/h}$)	< 200
	Analytical Results (1V Full Scale) - Area Counts (0.01 μVs)	
	Components	> 330 000 for each component
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 300
Wander (μV)	< 1 400
Drift ($\mu\text{V}/\text{h}$)	< 2 000
Analytical Results (10V Full Scale) - Area Counts (μVs)	
Components	> 33 000 for each component
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 3 000
Wander Counts)	< 14 000
Drift (Counts/h)	< 20 000
Analytical Results Area Counts (Cts*s)	
Components	> 330 000 for each component
	

Analytical Acceptance Comments	
1	In case of TCD equipped with the polyimide coated filaments, the acceptance values of the Components Area will result to be 4 times lower than the values reported in Table 25-3.

Checkout Using TCD with PKD Injector

Chapter at a Glance...

SOP Number: P0328/11/E - 01 September 2009	350
Scope	350
Parts Referenced	350
Analytical Conditions Required for Packed Injector.....	352
Recommended Initial Operations	353

Operating Procedures

TCD-PKD Checkout	354
------------------------	-----

SOP Number: P0328/11/E - 01 September 2009

Scope

Use the following procedure to verify proper TCD operation with Packed Injector.

Parts Referenced

Table 26-1. TCD-PKD Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	2 mm ID tapered	453 220 50
Liner Seal	Graphite seal for glass liner	290 334 05
Pre-column	Fused Silica Capillary Column 0.5 mt long 0.53 mm ID	260 603 75
Retaining Nut	M4 capillary column retaining nut	350 324 23
Press-fit connections	For columns 0.53/0.32 mm ID	350 438 16
Injection side adapter	For Wide bore column	347 003 03
Retaining Nut	For Injection side adapter	350 024 04
Detector side adapter	For Wide bore column	347 103 04
Ferrule	6 mm ID double brass ferrule	290 341 37
Nut	1/4" G-6 mm ID nut	350 201 18
Graphite Ferrule	Graphite ferrule for 0.53 mm ID Column Graphite ferrule for 0.32 mm ID Column	290 134 86 290 134 87
Septum	Standard septum for Packed Injector	313 032 26
Syringe	10 µl size; 50 mm needle length	365 005 25

Table 26-1. TCD-PKD Parts Referenced (Continued)

Part	Description	Part Number
Test Mixture	Three components in n-Hexane: <i>Component</i> <i>Concentration</i> Dodecane 200 µg/ml Tetradecane 200 µg/ml Hexadecane 200 µg/ml	338 190 16
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for Packed Injector

Table 26-2. TCD-PKD Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Make-up: Helium = 27.5 ml/min Reference Gas: Helium = 30 ml/min
Oven Program	Iso Temperature = 50 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 190 °C Final Time = 1 minute
Injector	Operating Mode = Packed Temperature = 180 °C
Detector	Block Temperature = 200 °C Transfer Temperature = 190 °C Constant Filament Temperature = No Filament Voltage = 10V Filament Temperature limit = 350 °C (*) Gain = x 10 Negative Polarity = No
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

(*) In case of TCD with the polyimide coated filaments, set the filament temperature limit to 320 °C.

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Insert the glass liner.
Remove the column and insert the 2 mm ID tapered glass liner, as required for the checkout, from the bottom of the injector with the appropriate liner seal. Fix the liner by using the appropriate adapter for capillary column. Refer to the TRACE GC Ultra Maintenance and Troubleshooting Manual.
2. Mount the adapter for capillary column on the detector base body.
3. Replace the septum
A new septum should be installed properly in your injector.
4. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
5. Install the precolumn.
Connect the precolumn to the injector.
6. Install the test column
Connect the test column to the precolumn by using the press fit connections provided.
7. Perform the manual leak check following the Operating Procedure “*Performing a Leak Check*” in Chapter 14 of the TRACE GC Ultra Operating Manual.
8. Connect the other end of the test column to the detector base body.
9. Connect your data handling.
Verify that your data handling is properly connected to your GC system

OPERATING PROCEDURE

TCD-PKD Checkout

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	PKD
Left carrier or Right carrier	He (helium)
Left detector or Right detector	TCD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select con pres. Scrool to Pressure and set the pressure value to have the required carrier gas flow rate Col.flow.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col.flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	50.0	50.0
Initial Time		1.00
Ramp 1		20.0
Final temp		190
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate PKD Injector Control Table. Set the required temperature setpoint Temp. Verify to operate in

Packed mode. If not, scroll to Mode, press **MODE/TYPE** to access the selection menu, then select Packed.

LEFT INLET (PKD) ¹		
Temp	180	180
Pressure	30.0	30.0
Mode:	Packed	

1. These settings could also be for a right inlet.

- Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate TCD Detector Control Table. Sequentially scroll to Block Temp, Transf Temp, Ref flow and mkup flow and set the required values.

LEFT DETECTOR (TCD) ¹		
Filament power	Off	
Fil status	(not rdy)	
Block temp	200	200
Transf temp	190	190
Const fil temp	Y/N	
Fil volts (CV)	10	
Fil temp limit ²	350	
Ref flow	30.0	30.0
Mkup flow	27.5	27.5<

1. These settings could also be for a right detector.

2. In case of TCD with the polyimide coated filaments, the temperature limit is 320 °C

- Scroll to Const fil temp and select it **NO**.
- Scroll to Fil volts (CV) and set the filament voltage. Scroll to Fil temp limit and set the required limit temperature setpoint.
- Scroll to Filament power and turn it **ON**. After a few seconds the ready is displayed on Fil status line.

- Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate TCD Detector Signal Control Table. Scroll to `Gain` and set the desired value.

LEFT SIGNAL (TCD) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Gain (x1..x10)	10<
Neg polarity?	N
Baseline comp	Off

1. These settings could also be for a right signal.

- Activate your Data System and set the parameters required for the checkout.
- In the TCD Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
- Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 26-3 according to the data handling in use.

- With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
- After the baseline evaluation has been completed, set-up the data system to acquire a single run.
- Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 26.1*

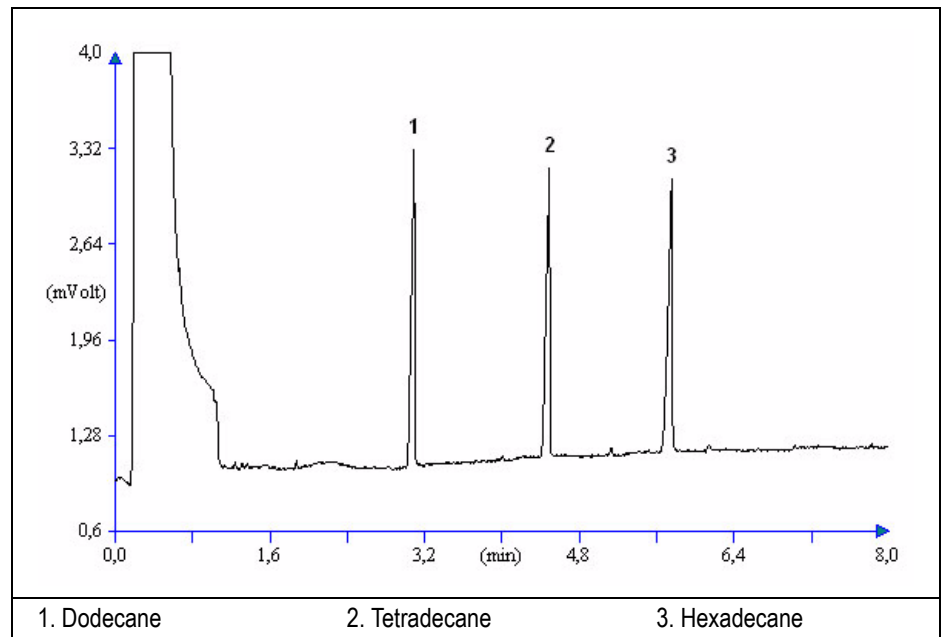





Figure 26-1. TCD-Packed Injection


15. The following criteria indicate successful completion of TCD-PKD checkout.
16. If these criteria are not met, repeat the test.


Table 26-3. TCD-PKD Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 30	< 300
	Wander (μV)	< 140	< 1 400
	Drift ($\mu\text{V/h}$)	< 200	< 2 000
	Analytical Results	Analog (1V Full Scale) Area Counts (0.1 μVs)	Digital (10V Full Scale) Area Counts (0.1 μVs)
	Components	> 30 000 for each component	> 300 000 for each component
			

Computing-integrator (e.g. ChromJet)	
	

CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 30
	Wander (μV)	< 140
	Drift ($\mu\text{V/h}$)	< 200
	Analytical Results (1V Full Scale) - Area Counts (0.01 μVs)	
	Components	> 300 000 for each component
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 300
Wander (μV)	< 1 400
Drift ($\mu\text{V}/\text{h}$)	< 2 000
Analytical Results (10V Full Scale) - Area Counts (μVs)	
Components	> 30 000 for each component
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 3 000
Wander Counts)	< 14 000
Drift (Counts/h)	< 20 000
Analytical Results Area Counts (Cts*s)	
Components	> 300 000 for each component
	

Analytical Acceptance Comments	
1	In case of TCD equipped with the polyimide coated filaments, the acceptance values of the Components Area will result to be 4 times lower than the values reported in Table 26-3.

Checkout Using TCD with PPKD Injector

Chapter at a Glance...

SOP Number: P0329/11/E - 01 September 2009	362
Scope	362
Parts Referenced.....	362
Analytical Conditions Required for Purged Packed Injector	364
Recommended Initial Operations	365

Operating Procedures

TCD-PPKD Checkout	366
-------------------------	-----

SOP Number: P0329/11/E - 01 September 2009

Scope

Use the following procedure to verify proper TCD operation with Purged Packed Injector.

Parts Referenced

Table 27-1. TCD-PPKD Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	2 mm ID tapered	453 220 50
Liner Seal	Graphite seal for glass liner	290 334 05
Pre-column	Fused Silica Capillary Column 0.5 mt long 0.53 mm ID	260 603 75
Retaining Nut	M4 capillary column retaining nut	350 324 23
Press-fit connections	For columns 0.53/0.32 mm ID	350 438 16
Injection side adapter	For Wide bore column	347 003 03
Retaining Nut	For Injection side adapter	350 024 04
Detector side adapter	For Wide bore column	347 103 04
Ferrule	6 mm ID double brass ferrule	290 341 37
Nut	1/4" G-6 mm ID nut	350 201 18
Graphite Ferrule	Graphite ferrule for 0.53 mm ID Column Graphite ferrule for 0.32 mm ID Column	290 134 86 290 134 87
Septum	Standard septum for Purged Packed Injector	313 032 26
Syringe	10 µl size; 50 mm needle length	365 005 25

Table 27-1. TCD-PPKD Parts Referenced (Continued)

Part	Description	Part Number
Test Mixture	Three components in n-Hexane: <i>Component</i> <i>Concentration</i> Dodecane 200 µg/ml Tetradecane 200 µg/ml Hexadecane 200 µg/ml	338 190 16
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for Purged Packed Injector

Table 27-2. TCD-PPKD Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Make-up: Helium = 27.5 ml/min Reference Gas: Helium = 30 ml/min
Oven Program	Initial Temperature = 50 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 190 °C Final Time = 1 minute
Injector	Operating Mode = Wide bore Septum Purge = Yes Temperature = 180 °C
Detector	Block Temperature = 200 °C Transfer Temperature = 190 °C Constant Filament Temperature = No Filament Voltage = 10V Filament Temperature limit = 350 °C (*) Gain = x 10 Negative Polarity = No
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

(*) In case of TCD with the polyimide coated filaments, set the filament temperature limit to 320 °C.

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Insert the glass liner.
Remove the column and insert the 2 mm ID tapered glass liner, as required for the checkout, from the bottom of the injector with the appropriate liner seal. Fix the liner by using the appropriate adapter for capillary column. Refer to the TRACE GC Ultra Maintenance and Troubleshooting Manual.
2. Mount the adapter for capillary column on the detector base body.
3. Replace the septum
A new septum should be installed properly in your injector.
4. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
5. Install the precolumn.
Connect the precolumn to the injector.
6. Install the test column
Connect the test column to the precolumn by using the press fit connections provided.
- 7.
8. Connect the other end of the test column to the detector base body.
9. Perform Column Evaluation.
10. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

TCD-PPKD Checkout

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	PPKD
Left carrier or Right carrier	He (helium)
Left detector or Right detector	TCD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to Pressure and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	50.0	50.0
Initial Time		1.00
Ramp 1		20.0
Final temp		200
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate PPKD Injector Control Table. Set the required temperature setpoint **Temp**. Verify to operate in

Wide bore mode. If not, scroll to Mode, press **MODE/TYPE** to access the selection menu, then select Wide bore.

LEFT INLET (PPKD) ¹		
Temp	180	180
Pressure	30.0	30.0
Mode:	Wide bore<	
Constant sept purge?	Y<	

1. These settings could also be for a right inlet.

- Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate TCD Detector Control Table. Sequentially scroll to Block Temp, Transf Temp, Ref flow and mkup flow and set the required values.

LEFT DETECTOR (TCD) ¹		
Filament power	Off	
Fil status	(not rdy)	
Block temp	200	200
Transf temp	190	190
Const fil temp	Y/N	
Fil volts (CV)	10	
Fil temp limit ¹	350	
Ref flow	30.0	30.0
Mkup flow	27.5	27.5<

1. These settings could also be

2. In case of TCD with the polyimide coated filaments, the temperature limit is 320 °C

- Scroll to Const fil temp and select it **NO**.
- Scroll to Fil volts (CV) and set the filament voltage. Scroll to Fil temp limit and set the required limit temperature setpoint.
- Scroll to Filament power and turn it **ON**. After a few seconds the ready is displayed on Fil status line.

- Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate TCD Detector Signal Control Table. Scroll to `Gain` and set the desired value.

LEFT SIGNAL (TCD) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Gain (x1..x10)	10<
Neg polarity?	N
Baseline comp	Off

1. These settings could also be for a right signal.

- Activate your Data System and set the parameters required for the checkout.
- In the TCD Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
- Perform a blank analysis injecting pure iso-octane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 27-3 according to the data handling in use.

- With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
- After the baseline evaluation has been completed, set-up the data system to acquire a single run.
- Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 27.1*

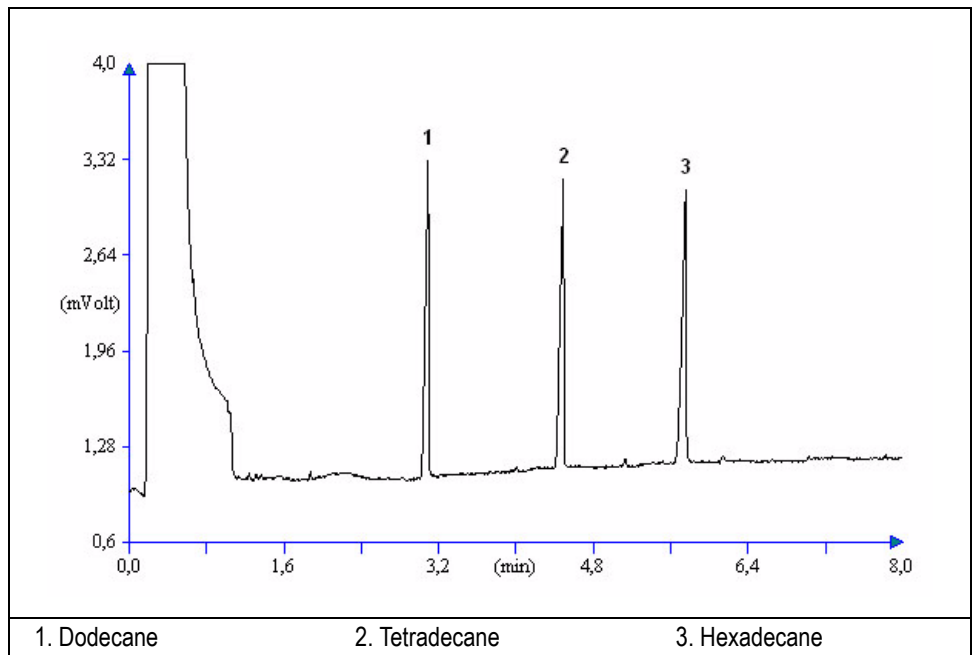





Figure 27-1. TCD-PPKD Injection


15. The following criteria indicate successful completion of TCD-PPKD checkout.
16. If these criteria are not met, repeat the test.


Table 27-3. TCD-PPKD Acceptance Criteria

Acceptance Values	CHROM-CARD		
	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 30	< 300
	Wander (μV)	< 140	< 1 400
	Drift ($\mu\text{V/h}$)	< 200	< 2 000
	Analytical Results	Analog (1V Full Scale) Area Counts (0.1 μVs)	Digital (10V Full Scale) Area Counts (0.1 μVs)
Components	> 30 000 for each component	> 300 000 for each component	
			

Computing-integrator (e.g. ChromJet)	
	

Acceptance Values	CHROMQUEST	
	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 30
	Wander (μV)	< 140
	Drift ($\mu\text{V/h}$)	< 200
	Analytical Results (1V Full Scale) - Area Counts (0.01 μVs)	
Components	> 300 000 for each component	
		

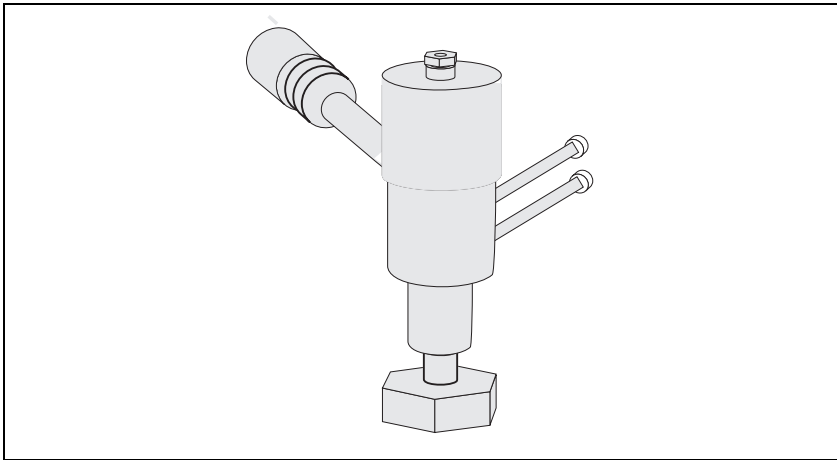
ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 300
Wander (μV)	< 1 400
Drift ($\mu\text{V}/\text{h}$)	< 2 000
Analytical Results (10V Full Scale) - Area Counts (μVs)	
Components	> 30 000 for each component
	

XCALIBUR		
Acceptance Values	Baseline Parameters (Acquisition Frequency = 10 Hz)	
	Noise (Counts)	< 3 000
	Wander Counts)	< 14 000
	Drift (Counts/h)	< 20 000
	Analytical Results Area Counts (Cts*s)	
	Components	> 300 000 for each component
		

Analytical Acceptance Comments	
1	In case of TCD equipped with the polyimide coated filaments, the acceptance values of the Components Area will result to be 4 times lower than the values reported in Table 27-3.

SECTION VIII

SOPs Using PDD



The *SOPs Using PDD* section, contains the procedures to test the TRACE GC Ultra with the Pulsed Discharge Detector (PDD) using different injectors.

Chapter 28, *Checkout Using PDD with S/SL Injector*.

Chapter 29, *Checkout Using PDD with OC Injector*.

Checkout Using PDD with S/SL Injector

Chapter at a Glance...

SOP Number: P0381/05/E - 01 September 2009	376
Scope	376
Parts Referenced	376
Analytical Conditions Required for Splitless Injection	377
Recommended Initial Operations	378

Operating Procedures

PDD-S/SL Checkout in Splitless Mode	379
---	-----

SOP Number: P0381/05/E - 01 September 2009

Scope

Use the following procedure to verify proper PDD operation with the Split/Splitless Injector.

Parts Referenced

Table 28-1. PDD-S/SL Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness.	260 800 01
Glass Liner	3 mm ID for splitless injection	453 200 32
Liner Seal	Graphite seal for glass liner	290 334 06
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
Septum	Standard septum for S/SL injector	313 032 11
Syringe	10 µl size; 70 mm needle length	365 001 03
Test Mixture	Three components in n-Hexane: <i>Component</i> <i>Concentration</i> Dodecane 1 µg/ml Tetradecane 1 µg/ml Hexadecane 1 µg/ml	338 190 32
Gases	Helium Chromatographic high grade purity (99.999%)	
Helium Purifier	Helium Purifier (VALCO)	432 100 76
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for Splitless Injection

Table 28-2. PDD-S/SL Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Discharge Gas: Helium = 30 ml/min (fixed value)
Oven Program	Initial Temperature = 60 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 160 °C Final Time = 1 minute
Injector	Operating Mode = Splitless Temperature = 250 °C Splitless Time = 0.5 minutes Split Flow = 60 ml/min Constant Septum Purge = Yes
Detector	Base Temperature = 280 °C Detector Signal Range = 10 ⁰
Injected Volume	1 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Replace the glass liner.
The glass liner currently installed in your injector should be carefully removed and replaced with the 3 mm ID glass liner for splitless application, as required for the checkout, with the appropriate liner seal.
2. Replace the septum
A new septum should be installed properly in your injector.
3. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
4. Verify that the helium purifier has been properly installed and purged.
5. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
6. Perform Column Evaluation and Leak Test.
7. Connect your data handling.
Verify that your data handling is properly connected to your GC system.



CAUTION

When leak check has been successfully carried out, power the helium purifier On. Before starting checkout, wait about 2.5 hours to reach the complete activation of the helium purifier.

OPERATING PROCEDURE

PDD-S/SL Checkout in Splitless Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	S/SL
Left carrier or Right carrier	He (helium)
Left detector or Right detector	PDD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to Pressure and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	60.0	60.0
Initial Time		1.00
Ramp 1		20.0
Final temp		160
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate Split/Splitless Injector Control Table. Set the required temperature *Temp* setpoint. Verify to

operate in **Splitless** mode. If not, scroll to *Mode*, press **MODE/TYPE** to access the selection menu, then select *Splitless*. Scroll to *Splitless time* to set the required setpoint.

LEFT INLET (S/SL) ¹		
Temp	250	250
Pressure	30.0	30.0
Mode:	Splitless	
Total flow	(63.0)	
Split Flow	60.0	60.0
Splitless time	0.50	0.50
Constant sept purge?		Y<

1. These settings could also be for a right inlet.

4. Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate PDD Detector Control Table. Set the required temperature *Base Temp* then turn the *Pulse generator* **ON**.

LEFT DETECTOR (PDD) ¹		
Pulse generator		On
Base temp	250	250
Signal pA	(1900.0)	

1. These settings could also be for a right detector.

5. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate PDD Detector Signal Control Table. Scroll to *Range* and set the electrometer amplifier input range required.

LEFT SIGNAL (PDD) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline Comp	Off

1. These settings could also be for a right signal.

6. Activate your Data System and set the parameters required for the checkout.
7. In the PDD Detector Signal Control Table, scroll to `Auto zero?` and turn it **YES**.
8. Perform a blank analysis injecting pure hexane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 28-3 according to the data handling in use.

9. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
10. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
11. Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 28.1*.

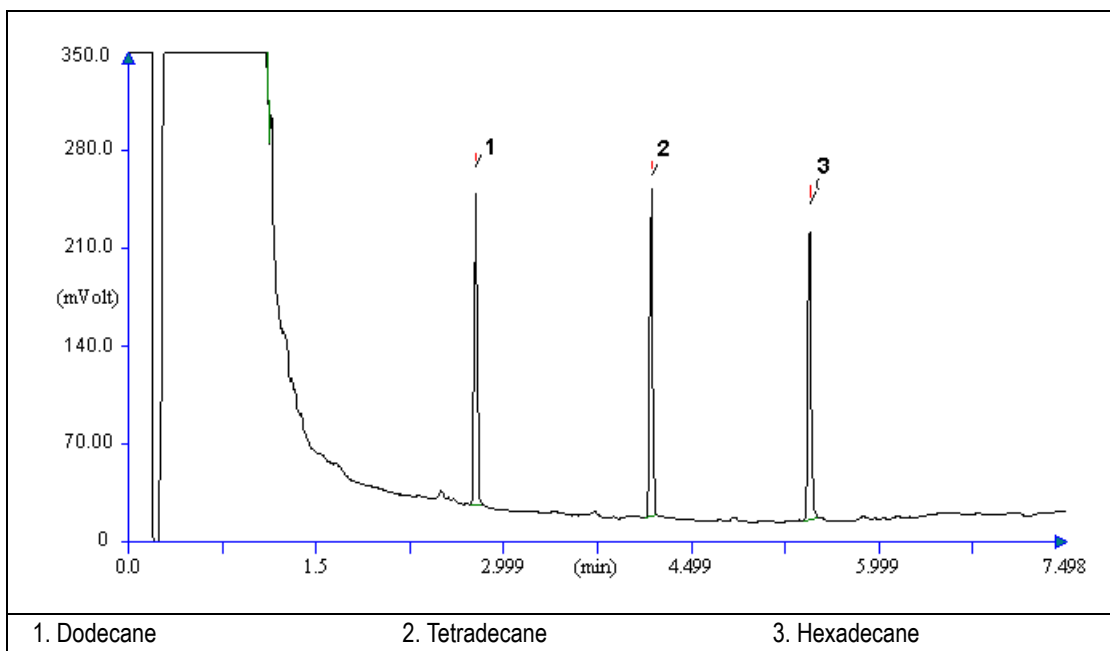





Figure 28-1. PDD-Splitless Injection


12. The following criteria indicate successful completion of PDD-S/SL checkout.
13. If these criteria are not met, repeat the test.


Table 28-3. PDD-S/SL Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 50	< 500
	Wander (μV)	< 500	< 5 000
	Drift ($\mu\text{V}/\text{h}$)	< 500	< 5 000
	Analytical Results	Analog (1V Full Scale) Area Counts (0.1 μVs)	Digital (10V Full Scale) Area Counts (0.1 μVs)
	Components	> 2 000 000 for each component	> 20 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1	1 ± 0.1
			

Computing-integrator (e.g. ChromJet)	
	

CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 50
	Wander (μV)	< 500
	Drift ($\mu\text{V}/\text{h}$)	< 500
	Analytical Results (1V Full Scale) - Area Counts (0.01 μVs)	
	Components	> 20 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 500
Wander (μV)	< 5 000
Drift ($\mu\text{V/h}$)	< 5 000
Analytical Results (10V Full Scale) - Area Counts (μVs)	
Components	> 2 000 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

XCALIBUR		
Acceptance Values	Baseline Parameters (Acquisition Frequency = 10 Hz)	
	Noise (Counts)	< 5 000
	Wander Counts)	< 50 000
	Drift (Counts/h)	< 50 000
	Analytical Results Area Counts (Cts*s)	
	Components	> 20 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1
		

Checkout Using PDD with OC Injector

Chapter at a Glance...

SOP Number: P0382/06/E - 01 September 2009	386
Scope	386
Parts Referenced	386
Analytical Conditions Required for On-Column Injection	387
Recommended Initial Operations	388

Operating Procedures

PDD-OCI Checkout in On-Column Mode	389
--	-----

SOP Number: P0382/06/E - 01 September 2009

Scope

Use the following procedure to verify proper PDD operation with the Cold On-Column Injector.

Parts Referenced

Table 29-1. PDD-OCI Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film.thickness.	260 800 01
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
Vespel Ferrule	Vespel Ferrule for 0.32 mm ID Column	290 134 60
Syringe	10 µl size; 75 mm needle length	365 020 07
Test Mixture	Three components in n-Hexane: <i>Component</i> <i>Concentration</i> Dodecane 1 µg/ml Tetradecane 1 µg/ml Hexadecane 1 µg/ml	338 190 32
Gases	Helium Chromatographic high grade purity (99.999%)	
Helium Purifier	Helium Purifier (VALCO)	432 100 76
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	
Syringe	10 µl size; 80 mm needle length	365 020 19
Pre-column	2 m long; 0.53 mm ID	260 603 75
Press-fit set	Set of Press-fir connectors for TRACE OC	350 038 45

Analytical Conditions Required for On-Column Injection

Table 29-2. PDD-OCI Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Discharge Gas: Helium = 30 ml/min (fixed value)
Oven Program	Initial Temperature = 70 °C Initial Time = 1 minute Ramp 1 = 20 °C/minute Final Temperature = 160 °C Final Time = 1 minute
Injector	Secondary Cooling = 0.2 minutes
Detector	Base Temperature = 280 °C Detector Signal Range = 10 ⁰
Injected Volume	1 µl of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
2. Install the test column.
The column currently installed should be carefully removed and replaced with the required test column.
In case of automatic On-column for TriPlus sampler, install the pre-column and connect it to the test column by press-fit connector.
3. Install and connect the TriPlus sampler and its components.
4. Perform Column Evaluation and Leak Test.
5. Connect your data handling.
Verify that your data handling is properly connected to your GC system.
6. Verify the opening/closing of the OC injector actuator by using the proper commands.
7. Verify the alignment of the syringe on the OC injector.



CAUTION

When leak check has been successfully carried out, power the helium purifier On. Before starting checkout, wait about 2.5 hours to reach the complete activation of the helium purifier.

OPERATING PROCEDURE

PDD-OCI Checkout in On-Column Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	OCI
Left carrier or Right carrier	He (helium)
Left detector or Right detector	PDD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to Pressure and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature *Temp* and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		160
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate Cold On-Column Injector Control Table. Scroll to `Sec. cool time` and set the required secondary cooling time.

LEFT INLET (OCI) ¹		
Pressure	30.0	30.0
Sec. Cool Time		0.2<

1. These settings could also be for a right inlet.

4. Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate PDD Detector Control Table. Set the required temperature `Base Temp` then turn the `Pulse generator` **ON**.

LEFT DETECTOR (PDD) ¹		
Pulse generator		On
Base temp	250	250
Signal pA		(1900.0)

1. These settings could also be for a right detector.

5. Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate PDD Detector Signal Control Table. Scroll to `Range` and set the electrometer amplifier input range required.

LEFT SIGNAL (PDD) ¹		
Output		(1000)
Offset		100
Auto zero?		Y/N
Range 10^(0..3)		0<
Analog filter		Off
Baseline comp		Off

1. These settings could also be for a right signal.

6. Activate your Data System and set the parameters required for the checkout.

7. In the PDD Detector Signal Control Table scroll to `Auto zero?` and turn it **YES**.
8. Perform a blank analysis injecting pure hexane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



Refer to the Acceptance Values reported in the Table 29-3 according to the data handling in use.

9. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
10. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
11. Perform the analysis.

Manual injection

- Inject the test mixture and press **START** on the GC to begin the checkout run.

Automatic injection with TriPlus sampler

- Fill a vial with the standard mix and place that vial in the sample tray.
- Load the method for OC and perform the sampling.

The resulting chromatogram should look like the one shown in *Figure 29.1*.

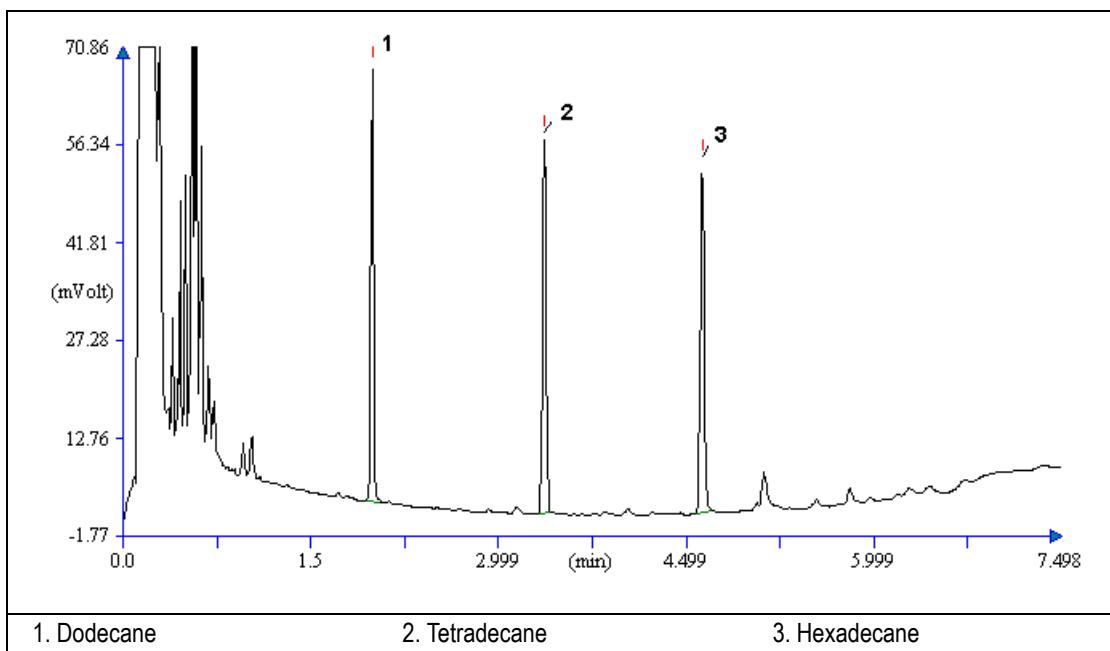





Figure 29-1. PDD-On-Column Injection


12. The following criteria indicate successful completion of PDD-OCI checkout.
13. If these criteria are not met, repeat the test.


Table 29-3. PDD-OCI Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 50	< 500
	Wander (μV)	< 500	< 5 000
	Drift ($\mu\text{V/h}$)	< 500	< 5 000
	Analytical Results	Analog (1V Full Scale) Area Counts (0.1 μVs)	Digital (10V Full Scale) Area Counts (0.1 μVs)
	Components	> 1 000 000 for each component	> 10 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1	1 ± 0.1
			

Computing-integrator (e.g. ChromJet)	
	

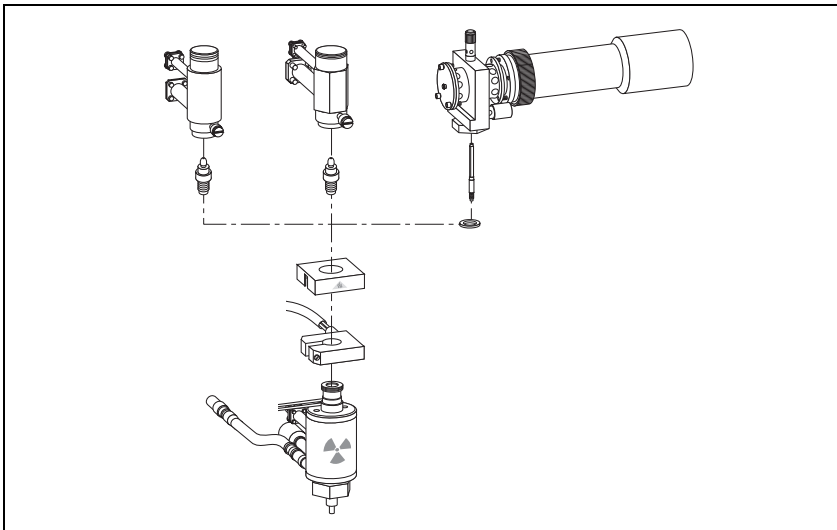
CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 50
	Wander (μV)	< 500
	Drift ($\mu\text{V/h}$)	< 500
	Analytical Results (1V Full Scale) - Area Counts (0.01 μVs)	
	Components	> 10 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 500
Wander (μV)	< 5 000
Drift ($\mu\text{V}/\text{h}$)	< 5 000
Analytical Results (10V Full Scale) - Area Counts (μVs)	
Components	> 1 000 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 5 000
Wander Counts)	< 50 000
Drift (Counts/h)	< 50 000
Analytical Results Area Counts (Cts*s)	
Components	> 10 000 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

SECTION IX

SOPs Using FID-NPD-FPD in Stacked Configuration



The *SOPs Using FID-NPD-FPD in Stacked Configuration* section, contains the procedures to test the TRACE GC Ultra with the Flame Ionization Detector (FID), Nitrogen-Phosphorus Detector (NPD) or Flame Photometric Detector (FPD) in series (stacked configuration) with the Electron Capture Detector ECD using different injectors.

Chapter 30, *Checkout Using Tandem FID*.

Chapter 31, *Checkout Using Tandem NPD*.

Chapter 32, *Checkout Using Tandem FPD*.

Checkout Using Tandem FID

Chapter at a Glance...

SOP Number: P0383/05/E - 01 September 2009	398
Scope	398
Checkout Overview	399
Important Considerations	400

Operating Procedures

Example of FID Tandem Checkout	401
--------------------------------------	-----

SOP Number: P0383/05/E - 01 September 2009

Scope

Use the following procedure to verify proper Flame Ionization Detector (FID) installed in series (stacked configuration, see Figure 30-1) with the non-destructive Electron Capture Detector ECD using different injectors.

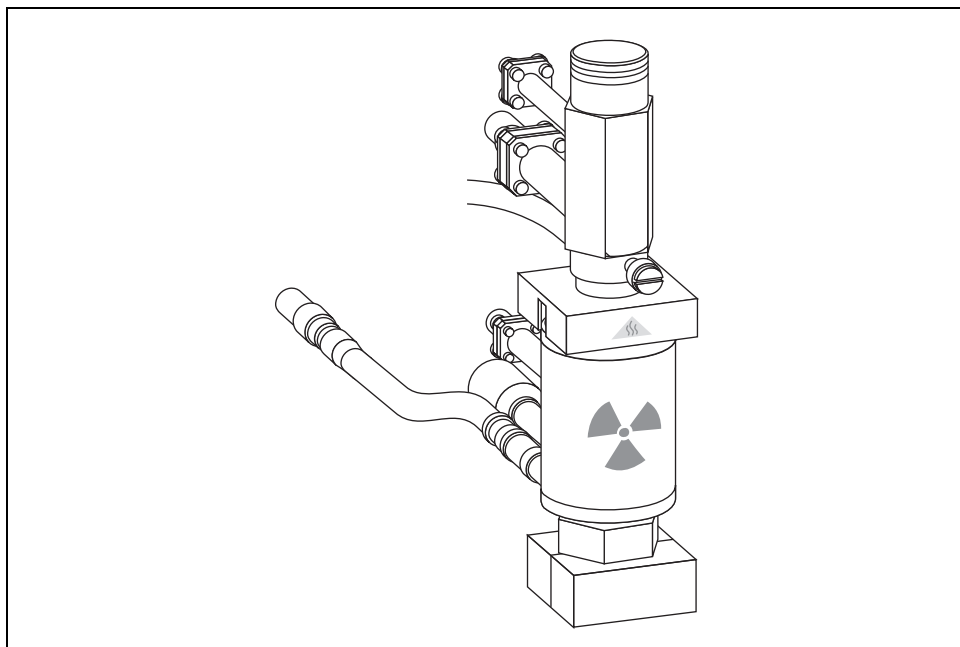


Figure 30-1. FID Tandem

Checkout Overview

The checkout must be carried out for each single ECD and FID detector, according to the injector used, referring to the relevant SOPs as reported in the following table.

Table 30-1. FID-ECD SOPs Reference

Detector	Refer To:
ECD	<i>Checkout Using ECD with S/SL Injector</i> on page 95.
	<i>Checkout Using ECD with OC Injector</i> on page 107.
	<i>Checkout Using ECD with PKD Injector</i> on page 119.
	<i>Checkout Using ECD with PPKD Injector</i> on page 131.
	<i>Checkout Using ECD with PTV Injector</i> on page 143.
FID	<i>Checkout Using FID with S/SL Injector</i> on page 37.
	<i>Checkout Using FID with OC Injector</i> on page 47.
	<i>Checkout Using FID with PKD Injector</i> on page 57.
	<i>Checkout Using FID with PPKD Injector</i> on page 69.
	<i>Checkout Using FID with PTV Injector</i> on page 81.



WARNING! To perform ECD checkout, refer to the relevant operating procedures reporting in Section III of this manual.

Before starting FID checkout procedures, it is strongly recommended to read the paragraph *Important Considerations* on page 400.

Important Considerations

This paragraph details the differences for FID checkout respect to the standard one reported in Section II of this manual.

FID Gas Required

In FID Tandem (stacked) configuration the FID only requires air and hydrogen as fuel gas to supply the flame.

The make-up gas is supplied by ECD.

Column Installation

When performing the checkout of the FID in stacked configuration it is not necessary any adjustment of the test column insertion depth. The test column remains connected to the ECD with the column insertion depth defined for this detector (109 mm measured from the bottom of the ferrule).

FID Detector and Signal Menus

When in stacked configuration, the FID is configured as **Auxiliary Detector**, then **AUX DETECTOR** and **AUX SIGNAL** instead of **LEFT/RIGHT DETECTOR** and **LEFT/RIGHT SIGNAL** must be pressed to access the relevant detector and signal menus.

The *Example of FID Tandem Checkout* operating procedure, on page 401, details the different procedure points respect to the standard FID checkout procedures reported in Section II of this manual.

OPERATING PROCEDURE

Example of FID Tandem Checkout

This procedure reports the different sequence points respect to the standard operating procedures reported in Section II. In the example, the S/SL injector is considered.

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	According to injector in use
Left carrier or Right carrier	He (helium)
Aux detector	FID

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scrool to **Pressure** and set the pressure value to have the required carrier gas flow rate **Col. flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	50.0	50.0
Initial Time		1.00
Ramp 1		20.0
Final temp		200
Final time 1		1.00 <
Ramp 2		Off

- Use **LEFT INLET** or **RIGHT INLET** to display the appropriate Split/Splitless Injector Control Table. Set the required temperature *Temp* setpoint. Verify to operate in **Splitless** mode. If not, scroll to *Mode*, press **MODE/TYPE** to access the selection menu, then select *Splitless*. Scroll to *Splitless time* to set the required setpoint.

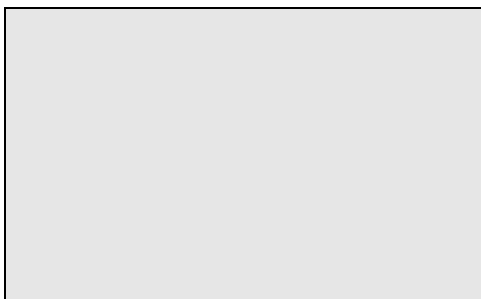
LEFT INLET (S/SL) ¹		
Temp	230	230
Pressure	30.0	30.0
Mode:	Splitless	
Total flow	(63.0)	
Split Flow	60.0	60.0
Splitless time	0.80	0.80
Constant sept purge?		Y<

1. These settings could also be for a right inlet.

- Use **AUX DETECTOR** to display the appropriate FID Detector Control Table. Set the required temperature **Base Temp** and the detector gases **H2** and **Air** required setpoints.

AUX DETECTOR (FID)		
Flame		Off
Base temp	250	250
Signal pA		(5.5)
Ign. thresh		2.0
Flameout retry		Off
H2	35	35
Air	350	350<

- Since the make-up gas is supplied by ECD detector, set the value into the ECD detector Control Table.



1. These settings could also be for a right detector.

6. Ignite the FID flame scrolling to `Flame` and pressing **ON**.
7. Use **AUX SIGNAL** to display the appropriate FID Detector Signal Control Table. Observe the FID flame signal at the display. This is the flame-on background offset. Scroll to `Range` and set the electrometer amplifier input range required.

AUX SIGNAL (FID)	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline Comp	Off

8. Activate your Data System and set the parameters required for the checkout.
9. In the Aux Detector Signal Control Table, scroll to `Auto zero?` and turn it **YES**.
10. Perform a blank analysis injecting pure hexane and press **START** on the GC to begin the checkout run.

Continue with Baseline Acquisition and Analysis.

Checkout Using Tandem NPD

Chapter at a Glance...

SOP Number: P0384/06/E - 01 September 2009	406
Scope	406
Checkout Overview	407
Important Considerations	408

Operating Procedures

Example of NPD Tandem Checkout	409
--------------------------------------	-----

SOP Number: P0384/06/E - 01 September 2009

Scope

Use the following procedure to verify proper Nitrogen Phosphorus Detector (NPD) installed in series (stacked configuration, see Figure 31-1) with the non-destructive Electron Capture Detector ECD using different injectors.

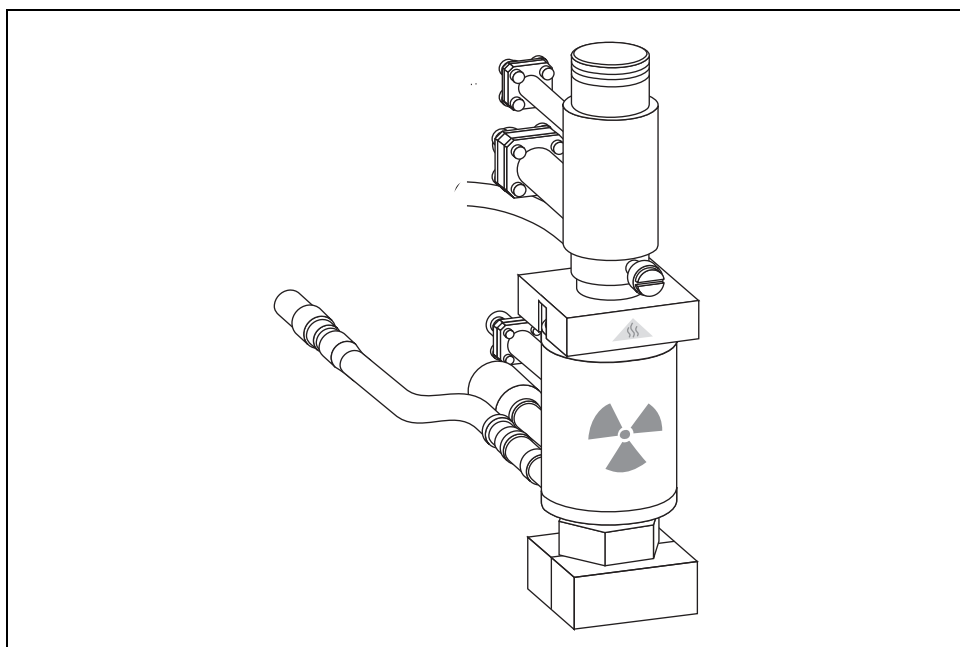


Figure 31-1. NPD Tandem

Checkout Overview

The checkout must be carried out for each single ECD and NPD detector, according to the injector used, referring to the relevant SOPs as reported in the following table.

Table 31-1. NPD-ECD SOPs Reference

Detector	Refer To:
ECD	<i>Checkout Using ECD with S/SL Injector</i> on page 95.
	<i>Checkout Using ECD with OC Injector</i> on page 107.
	<i>Checkout Using ECD with PKD Injector</i> on page 119.
	<i>Checkout Using ECD with PPKD Injector</i> on page 131.
	<i>Checkout Using ECD with PTV Injector</i> on page 143.
NPD	<i>Checkout Using NPD with S/SL Injector</i> on page 157.
	<i>Checkout Using NPD with OC Injector</i> on page 169.
	<i>Checkout Using NPD with PKD Injector</i> on page 181.
	<i>Checkout Using NPD with PPKD Injector</i> on page 195.
	<i>Checkout Using NPD with PTV Injector</i> on page 209.



WARNING! To perform ECD checkout, refer to the relevant operating procedures reporting in Section III of this manual.

Before starting NPD checkout procedures, it is strongly recommended to read the paragraph *Important Considerations* on page 408.

Important Considerations

This paragraph details the differences for NPD checkout respect to the standard one reported in Section IV of this manual.

NPD Gas Required

In NPD Tandem (stacked) configuration the NPD only requires air and hydrogen as fuel gas to supply the flame.

The make-up gas is supplied by ECD.

Column Installation

When performing the checkout of the NPD in stacked configuration it is not necessary any adjustment of the test column insertion depth. The test column remains connected to the ECD with the column insertion depth defined for this detector (109 mm measured from the bottom of the ferrule). For that reason, it is strongly recommended the use of the silcostealed jet instead of the standard one.

NPD Detector and Signal Menus

When in stacked configuration, the NPD is configured as **Auxiliary Detector**, then **AUX DETECTOR** and **AUX SIGNAL** instead of **LEFT/RIGHT DETECTOR** and **LEFT/RIGHT SIGNAL** must be pressed to access the relevant detector and signal menus.

The *Example of NPD Tandem Checkout* operating procedure, on page 409, details the different procedure points respect to the standard NPD checkout procedures reported in Section IV of this manual.

OPERATING PROCEDURE

Example of NPD Tandem Checkout

This procedure reports the different sequence points respect to the standard operating procedures reported in Section IV. In the example, the S/SL injector is considered.

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	According to injector in use
Left carrier or Right carrier	He (helium)
Aux detector	NPD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scroll to **Pressure** and set the pressure value to have the required carrier gas flow rate **Col.flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col.flow	3.00	
Lin. veloc.		(60.9)<

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		230
Final time 1		1.00<
Ramp 2		Off

- Use **LEFT INLET** or **RIGHT INLET** to display the appropriate Split/Splitless Injector Control Table. Set the required temperature *Temp* setpoint. Verify to operate in **Splitless** mode. If not, scroll to *Mode*, press **MODE/TYPE** to access the selection menu, then select *Splitless*. Scroll to *Splitless time* to set the required setpoint.

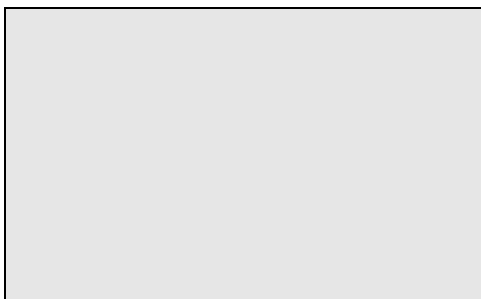
LEFT INLET (S/SL) ¹		
Temp	230	230
Pressure	30.0	30.0
Mode:	Splitless	
Total flow	(63.0)	
Split Flow	60.0	60.0
Splitless time	0.80	0.80
Constant sept purge?		Y<

1. These settings could also be for a right inlet.

- Use **AUX DETECTOR** to display the appropriate NPD Detector Control Table. Set the required temperature **Base Temp** and the detector gases **H2** and **Air** required setpoints.

AUX DETECTOR (NPD)		
Source cur,A		Off
Base temp	300	300
Signal pA		(10.4)
Target curr. pA		(X.XX)
Auto adjust		No
Polarizer V		3.5
H2 delay time		Off
H2	2.3	2.3
Air	60	60

- Since the make-up gas is supplied by ECD detector, set the value into the ECD detector Control Table.



1. These settings could also be for a right detector.
6. Scroll to `Polarizer V` and set the required setpoint (3.5 V).
7. Scroll to `Source cur,A` and turn on the source as described in the relevant *Checkout Using NPD* procedure.
8. Use AUX Signal to display the appropriate NPD Detector Signal Control Table. Scroll to `Range` and set the electrometer amplifier input range required.

AUX SIGNAL (NPD)	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline Comp	Off

9. Activate your Data System and set the parameters required for the checkout.
10. In the Aux Detector Signal Control Table, scroll to `Auto zero?` and turn it **YES**.
11. Perform a blank analysis injecting pure hexane and press **START** on the GC to begin the checkout run.

Continue with Baseline Acquisition and Analysis.

Checkout Using Tandem FPD

Chapter at a Glance...

SOP Number: P0385/06/E - 01 September 2009	414
Scope	414
Checkout Overview	415
Important Considerations	416

Operating Procedures

Example of FPD Tandem Checkout	417
Example of FPD Tandem Checkout	417

SOP Number: P0385/06/E - 01 September 2009

Scope

Use the following procedure to verify proper Flame Photometric Detector (FPD) installed in series (stacked configuration, see Figure 32-1) with the non-destructive Electron Capture Detector ECD using different injectors.

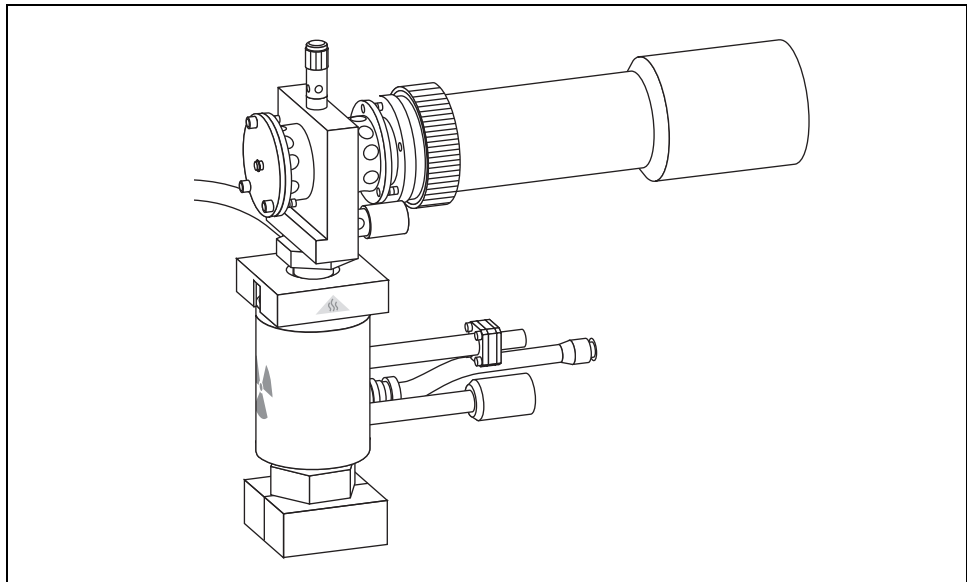


Figure 32-1. FPD Tandem

Checkout Overview

The checkout must be carried out for each single ECD and FPD detector, according to the injector used, referring to the relevant SOPs as reported in the following table.

Table 32-1. FPD-ECD SOPs Reference

Detector	Refer To:
ECD	<i>Checkout Using ECD with S/SL Injector</i> on page 95.
	<i>Checkout Using ECD with OC Injector</i> on page 107.
	<i>Checkout Using ECD with PKD Injector</i> on page 119.
	<i>Checkout Using ECD with PPKD Injector</i> on page 131.
	<i>Checkout Using ECD with PTV Injector</i> on page 143.
FPD	<i>Checkout Using FPD with S/SL Injector</i> on page 225.
	<i>Checkout Using FPD with OC Injector</i> on page 237.
	<i>Checkout Using FPD with PKD Injector</i> on page 249.
	<i>Checkout Using FPD with PPKD Injector</i> on page 263.
	<i>Checkout Using FPD with PTV Injector</i> on page 277.



WARNING! To perform ECD checkout, refer to the relevant operating procedures reporting in Section III of this manual.

Before starting FPD checkout procedures, it is strongly recommended to read the paragraph *Important Considerations* on page 416.

Important Considerations

This paragraph details the differences for FPD checkout respect to the standard one reported in Section V of this manual.

FPD Gas Required

In FPD Tandem (stacked) configuration the FPD only requires air and hydrogen as fuel gas to supply the flame.

The make-up gas supplied by ECD has to be maintained.

Column Installation

When performing the checkout of the FPD in stacked configuration it is not necessary any adjustment of the test column insertion depth. The test column remains connected to the ECD with the column insertion depth defined for this detector (109 mm measured from the bottom of the ferrule). For that reason, it is strongly recommended the use of the silcosteeled jet instead of the standard one.

FPD Detector and Signal Menus

When in stacked configuration, the FPD is configured as **Auxiliary Detector**, then **AUX DETECTOR** and **AUX SIGNAL** instead of **LEFT/RIGHT DETECTOR** and **LEFT/RIGHT SIGNAL** must be pressed to access the relevant detector and signal menus.

The *Example of FPD Tandem Checkout* operating procedure, on page 417, details the different procedure points respect to the standard FPD checkout procedures reported in Section V of this manual.

OPERATING PROCEDURE

Example of FPD Tandem Checkout

This procedure reports the different sequence points respect to the standard operating procedures reported in Section V. In the example, the S/SL injector is considered.

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	According to injector in use
Left carrier or Right carrier	He (helium)
Aux detector	FPD

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select **con pres.** Scrool to **Pressure** and set the pressure value to have the required carrier gas flow rate **Col.flow**.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col.flow	3.00	
Lin. veloc.		(60.9)<

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	70.0	70.0
Initial Time		1.00
Ramp 1		20.0
Final temp		230
Final time 1		1.00<
Ramp 2		Off

- Use **LEFT INLET** or **RIGHT INLET** to display the appropriate Split/Splitless Injector Control Table. Set the required temperature *Temp* setpoint. Verify to operate in **Splitless** mode. If not, scroll to *Mode*, press **MODE/TYPE** to access the selection menu, then select *Splitless*. Scroll to *Splitless time* to set the required setpoint.

LEFT INLET (S/SL) ¹		
Temp	230	230
Pressure	30.0	30.0
Mode:	Splitless	
Total flow	(63.0)	
Split Flow	60.0	60.0
Splitless time	0.80	0.80
Constant sept purge?		Y<

1. These settings could also be for a right inlet.

- Use **AUX DETECTOR** to display the appropriate FPD Detector Control Table. Set the required temperature **FPD Temp** and the detector gases **H2** and **Air** required setpoints.

AUX DETECTOR (FPD)		
Flame		Off
FPD temp	150	150
Signal pA	(1.4)	
High voltage mode?		N
H2	90	90
Air	115	115

- Since the make-up gas is supplied by ECD detector, set the value into the ECD detector Control Table.

Base temp	250	250
ECD Temp	300	300
Ref current nA		1.0
Freq kHz		(2.20)
Pulse amp V		50
Pulse width μ s		1.0
Mkup (N2)	30	30<

1. These settings could also be for a right detector.

6. Verify that High voltage mode is set to **NO**.
7. Scroll to Flame and press **ON**. This start the ignition sequence. When ignition is confirmed, the photomultiplier tube is energized. The baseline level Signal pA, will suddenly increase meaning that the flame is lit inside the detector. After a few seconds, the baseline should stabilize to the standing current of the system.
8. Use AUX Signal to display the appropriate FPD Detector Signal Control Table. Scroll to Range and set the electrometer amplifier input range required.

AUX SIGNAL (FPD)	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Baseline Comp	Off

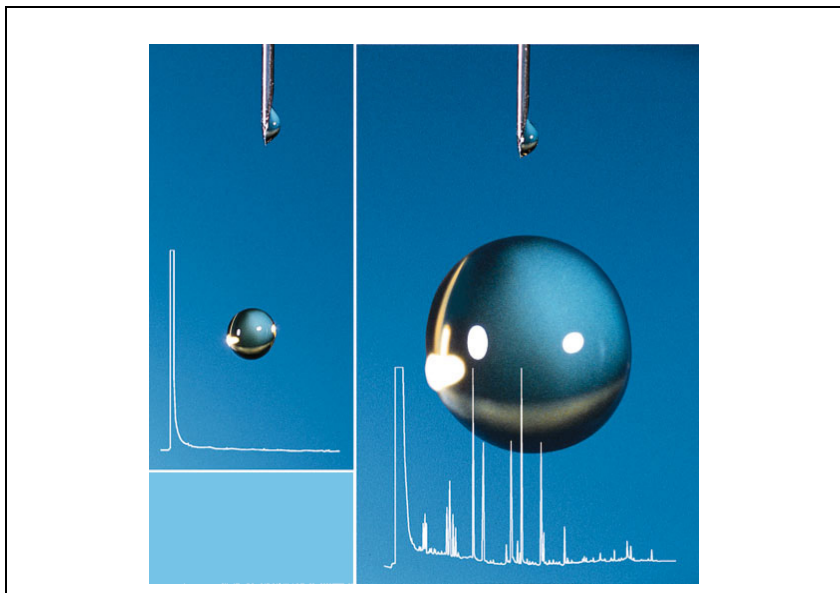
9. Activate your Data System and set the parameters required for the checkout.
10. In the Aux Detector Signal Control Table, scroll to Auto zero? and turn it **YES**.
11. Perform a blank analysis injecting pure hexane and press **START** on the GC to begin the checkout run.

Continue with Baseline Acquisition and Analysis.

SECTION

X

SOPs for Large Volume Applications



The *SOPs for Large Volume Applications* section, contains the procedures to test the TRACE GC Ultra for large volume application by using different injectors.

Chapter 33, *Checkout Using FID with LVSL Injector*

Chapter 34, *Checkout Using FID with LVOC Injector*

Checkout Using FID with LVSL Injector

Chapter at a Glance...

SOP Number: P0430/04/E - 01 September 2009	424
Scope	424
Parts Referenced	424
Analytical Conditions Required for LVSL Injection.....	426
Recommended Initial Operations	427

Operating Procedures

FID-LVSL Checkout in Splitless Mode	428
---	-----

SOP Number: P0430/04/E - 01 September 2009

Scope

Use the following procedure to verify proper FID operation with the LVSL Injector.

Parts Referenced

Table 33-1. FID-LVSL Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness	260 800 01
Pre-column	Retention gap DPTMDS deactivated; 5 m long 0.32 mm ID	260 800 10
Press fit	Deactivated universal press-tight connector (set of 5)	350 038 50
Glass Liner	5 mm ID for LVSL injection	453 020 65
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
Retaining Nut	M4 capillary column retaining nut	350 324 23
LVSL Adapter	LVSL adapter for 9 mm septa	347 093 46
Vespel seal	Vespel seal for LVSL adapter	356 034 50
Septum Cap	Cap for LVSL adapter	350 010 55
Septum holder	Metallic holder for 9 mm septa	233 030 20
Septum	9 mm septa for LVSL injector (set of 10)	313 032 41
Syringe	50 µl size; 50 mm needle length, 0.63C, conic tip	365 030 15

Table 33-1. FID-LVSL Parts Referenced

Part	Description	Part Number
Test Mixture	Three components in n-Hexane: <i>Component</i> <i>Concentration</i> Dodecane 1 µg/ml Tetradecane 1 µg/ml Hexadecane 1 µg/ml	338 190 32
Gases	Chromatographic-grade purity	
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur, Computing-integrator	

Analytical Conditions Required for LVSL Injection

Table 33-2. FID-LVSL Analytical Conditions

Parameters Setting	
Gases	Carrier Gas: Helium = 30 kPa Constant Pressure Hydrogen = 35 ml/min Air = 350 ml/min Make-up Gas: Nitrogen = 30 ml/min
Oven Program	Initial Temperature = 55 °C Initial Time = 2 minutes Ramp 1 = 25 °C/minute Final Temperature = 200 °C Final Time = 1 minute
Injector	Operating Mode = Splitless Temperature = 230 °C Splitless Time = 0.8 minutes Split Flow = 60 ml/min Constant Septum Purge = Off Stop Purge for = 0.8 minutes
Detector	Base Temperature = 250 °C Detector Signal Range = 10 ⁰
Injected Volume	20 µl + needle of Test Mixture
Analog Signal Output	Chrom-Card Acquisition Frequency = Medium
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Verify a “DPFC 2001” pneumatic module is installed.
2. Verify the LVSL kit is correctly installed (LVSL adapter, LV liner packed with glass wool).
3. Replace the 9-mm septum.
A new septum should be installed properly in your injector.
4. Connect the required gas lines
Verify the required gas supplies are properly connected to your GC.
5. Install the pre-column and connect the test column by means of the press fit.
The column currently installed should be carefully removed and replaced with the required test column.
6. Perform Column Evaluation and Leak Test.
7. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

FID-LVSL Checkout in Splitless Mode

Before beginning, press **CONFIG** to verify the GC configuration:

Left inlet or Right inlet	S/SL
Left carrier or Right carrier	He (helium)
Left detector or Right detector	FID

1. Use **LEFT CARRIER** or **RIGHT CARRIER** to display the appropriate Carrier Gas Control Table. Verify to operate in constant pressure mode. If not, scroll to Flow mode, press **MODE/TYPE** to access the selection menu, then select con pres. Scrol to Pressure and set the pressure value to have the required carrier gas flow rate Col. flow.

LEFT CARRIER ¹		
Pressure	30.0	30.0
Col. flow	3.00	
Lin. veloc.		(60.9) <

1. These settings could also be for a right carrier.

2. Use **OVEN** to display the Column Oven Control Table. Set the oven temperature and the Oven Program required.

OVEN		
Temp	55.0	55.0
Initial Time		2.00
Ramp 1		25.0
Final temp		200
Final time 1		1.00 <
Ramp 2		Off

3. Use **LEFT INLET** or **RIGHT INLET** to display the appropriate Split/Splitless Injector Control Table. Set the required temperature *Temp* setpoint. Verify to

operate in **Splitless** mode. If not, scroll to **Mode**, press **MODE/TYPE** to access the selection menu, then select **Splitless**. Scroll to **Splitless time** to set the required setpoint.

LEFT INLET (S/SL) ¹		
Temp	230	230
Pressure	30.0	30.0
Mode:	Splitless	
Total flow	(63.0)	
Split Flow	60.0	60.0
Splitless time	0.80	0.80
Constant sept purge?	N	
Stop purge for	0.8	

1. These settings could also be for a right inlet.

- Use **LEFT DETECTOR** or **RIGHT DETECTOR** to display the appropriate FID Detector Control Table. Set the required temperature **Base Temp** and the detector gases **H2**, **Air** and **Mkup** required setpoints.

LEFT DETECTOR (FID) ¹		
Flame	Off	
Base temp	250	250
Signal pA	(5.5)	
Ign.thresh	2.0	
Flameout retry	Off	
H2	35	35
Air	350	350
Mkup N2	30	30<

1. These settings could also be for a right detector.

- Ignite the FID flame scrolling to **Flame** and pressing **ON**.
- Use **LEFT SIGNAL** or **RIGHT SIGNAL** to display the appropriate FID Detector Signal Control Table. Observe the FID flame signal at the display. This is the

flame-on background offset. Scroll to `Range` and set the electrometer amplifier input range required.

LEFT SIGNAL (FID) ¹	
Output	(1000)
Offset	100
Auto zero?	Y/N
Range 10^(0..3)	0<
Analog filter	Off
Baseline Comp	Off

1. These settings could also be for a right signal.

7. Activate your Data System and set the parameters required for the checkout.
8. In the FID Detector Signal Control Table, scroll to `Auto zero?` and turn it **YES**.
9. Perform a blank analysis injecting pure hexane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 33-3 according to the data handling in use.

10. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
11. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
12. Inject the test mixture and press **START** on the GC to begin the checkout run. The resulting chromatogram should look like the one shown in *Figure 2.1*.

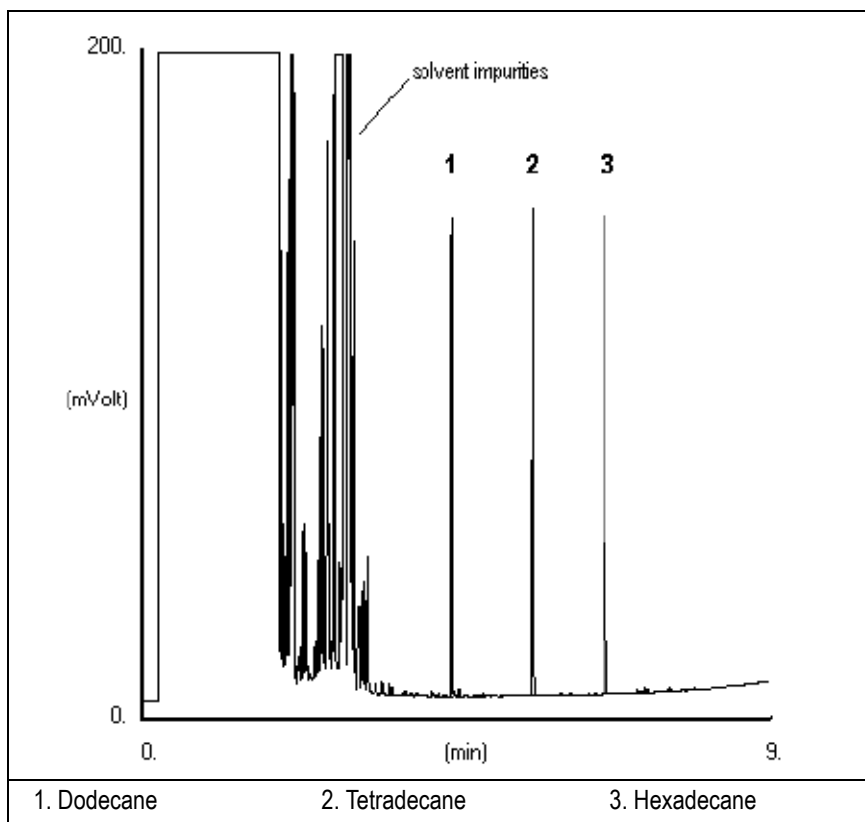





Figure 33-1. FID-LV Splitless Injection


13. The following criteria indicate successful completion of FID-LVSL checkout.
14. If these criteria are not met, repeat the test.

Table 33-3. FID-LVSL Acceptance Criteria

CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 30	< 300
	Wander (μV)	< 50	< 500
	Drift (μV/h)	< 100	< 1 000
	Analytical Results	Analog (1V Full Scale) Area Counts (0.1 μVs)	Digital (10V Full Scale) Area Counts (0.1 μVs)
	Components	> 1 500 000 for each component	> 15 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1	1 ± 0.1
			

CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 30
	Wander (μV)	< 50
	Drift (μV/h)	< 100
	Analytical Results (1V Full Scale) - Area Counts (0.01 μVs)	
	Components	> 15 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 300
Wander (μV)	< 500
Drift ($\mu\text{V}/\text{h}$)	< 1 000
Analytical Results (10V Full Scale) - Area Counts (μVs)	
Components	> 1 500 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 3 000
Wander Counts)	< 5 000
Drift (Counts/h)	< 10 000
Analytical Results Area Counts (Cts*s)	
Components	> 15 000 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

Analytical Acceptance Comments	
1	When the make-up gas is not used, the acceptance values will result to be 2.5 times lower than the values reported in Table 33-3.
2	When helium is used as make-up gas, the acceptance values will result to be 10 times lower than the values reported in Table 33-3.

Checkout Using FID with LVOC Injector

Chapter at a Glance...

SOP Number: TE P0609/02/E-01 September 2009	436
Scope	436
Parts Referenced	436
Recommended Initial Operations	440
Analytical Conditions Required for LVOC Injection	438

Operating Procedures

FID-LVOC Checkout in LVOC Mode	441
--------------------------------------	-----

SOP Number: TE P0609/02/E-01 September 2009

Scope

Use the following procedure to verify proper FID operation with the LVOC Injector.

Parts Referenced

Table 34-1. FID-LVOC Parts Referenced

Part	Description	Part Number
Test Column	Fused Silica Capillary Column TR-5; 7 m long 0.32 mm ID; 0.25 µm film thickness	260 800 01
Pre-column	UNCORET column, 15 m, 3 m coated with SE 54, 0.25 µm f.t.	260 604 27
Restrictor	Restrictor for SVE valve, 25 cm x 25µm ID	260 603 88
Connector	T-shaped connector	347 084 48
Vespel Ferrule	Vespel ferrule 0.9 mm hole (for connecting T-piece to SVE)	290 334 98
	Vespel ferrule for for 0.53 mm ID Column	290 134 71
Graphite Ferrule	Graphite ferrule for 0.32 mm ID Column	290 134 87
	Graphite ferrule for 0.53 mm ID Column	290 134 86
Retaining Nut	M4 capillary column retaining nut	350 324 23
	Column retaining nut and backwasher for OC	452 100 01
Seal	O-ring	290 113 02
Syringe	250 µl size; removable needle	365 004 90
Needle	(0,47/80 cone) for syringe p/n 365 004 90	365 664 80
Vial	2 ml screw-top vials	240 140 21
Cap	Screw-top cap for 2 ml vials p/n 240 140 21	386 060 92

Table 34-1. FID-LVOC Parts Referenced (Continued)

Part	Description	Part Number	
Test Mixture	Three components in n-Hexane:	338 190 32	
	<i>Component</i> <i>Concentration</i>		
	Dodecane 1 µg/ml		
	Tetradecane 1 µg/ml		
Hexadecane 1 µg/ml			
Sampler	TriPlus autosampler		
Gases	Chromatographic-grade purity		
Data Acquisition	Chrom-Card, ChromQuest, Atlas, Xcalibur		

Analytical Conditions Required for LVOC Injection

Analytical Condition for the GC

Table 34-2. FID-LVOC Analytical Conditions for the GC

GC Parameters Setting	
Gases	Carrier Gas: Helium = 60 kPa Constant Pressure SVE flow = > 10 ml/min Sweep flow = $\geq 0,05$ ml/min Hydrogen = 35 ml/min Air = 350 ml/min Make-up Gas: Nitrogen = 30 ml/min
Oven Program	Initial Temperature = 72 °C Initial Time = 2.5 minutes Ramp 1 = 50 °C/minute Final Temperature = 220 °C Final Time = 1 minute
Injector	Operating Mode = LVOC Secondary cooling = 6 s SVE Temperature = 150 °C SVE Duration = 12 s
Detector	Base Temperature = 250 °C Detector Signal Range = 10^0
Injected Volume	100 μ l of Test Mixture
Digital Signal Output	Chrom-Card, ChromQuest, Atlas, Xcalibur Acquisition Frequency = 10 Hz

Analytical Condition for the TriPlus Autosampler

Table 34-3. FID-LVOC Analytical Conditions for the TriPlus Autosampler

TriPlus Parameters Setting	
Sampling	Sampling vol (μL) = 100 Plunger strokes = 0 Air volume (μL) = 10 Filling volume (μL) = 110
Injection depth mode	Pre-injection dwell time (s) = 0,2 Post-injection dwell time (s) = 0,3 Injection depth (mm) = 71 Injection speed ($\mu\text{L/s}$) = 10
Sampling depth in vial	Sampling vial depth % = 100
Sample viscosity	Sample pull up speed ($\mu\text{L/s}$) = 5 Delay after bubble elimination (s) = 1 Viscosity delay (s) = 0.3
GC syncro start	Syncro type = Delayed
Advanced parameters	Wash solvent depth % = 96 Waste depth % = 20 Needle speed into vial (mm/s) = 100 Solvent filling pull-up speed ($\mu\text{L/s}$) = 10 Bubble elimination pull-up speed ($\mu\text{L/s}$) = 10 Delay between strokes (s) = 0.1

Recommended Initial Operations

Before starting the checkout, the following operations should be carried out:

1. Connect the required gas lines.
Verify the required gas supplies are properly connected to your GC.
2. Verify a “DPFC 2001” pneumatic module is installed.
3. Verify the LVOC actuator for TriPlus is correctly installed.
4. Replace the O-ring.
A new O-ring should be installed properly in your injector
5. Installation of the pre-column and test column
The column currently installed should be carefully removed and replaced with the required test column.
Install the pre-column and connect it to the test column and the SVE by using the T-shaped connector. Insert the column into the precolumn for about two cm.
6. Install and connect the TriPlus sampler and its components.
7. Verify the opening/closing of the OC injector actuator by using the proper commands.
8. Verify the alignment of the syringe on the OC injector.
9. Perform Column Evaluation and Leak Test.
10. Check the sweep flow at the outlet of the restrictor.
It should be ≥ 0.05 ml/min.
11. Check the SVE flow. It should be > 10 ml/min.
12. Connect your data handling.
Verify that your data handling is properly connected to your GC system.

OPERATING PROCEDURE

FID-LVOC Checkout in LVOC Mode

1. Set the GC parameters required to perform a LVOC injection listed in Table 34-2 on page 438.
2. Set the TriPlus parameters required to perform a LVOC injection listed in Table 34-3 on page 439
3. In the FID Detector Signal Control Table, scroll to `Auto zero?` and turn it **YES**.
4. Perform a blank analysis injecting pure hexane and press **START** on the GC to begin the checkout run.

Baseline Acquisition and Analysis



NOTE

Refer to the Acceptance Values reported in the Table 34-4 according to the data handling in use.

5. With the GC in Stand-by/Prep Run condition, activate the data system for 10 minutes to evaluate your baseline in isothermal condition.
6. After the baseline evaluation has been completed, set-up the data system to acquire a single run.
7. Fill one 2 ml vial with the sample and cap. Install the vial into the autosampler tray position 1.
8. Set up in the data system a sequence of three runs.
9. Perform the sequence. The resulting chromatogram should look like the one shown in *Figure 34-1*.

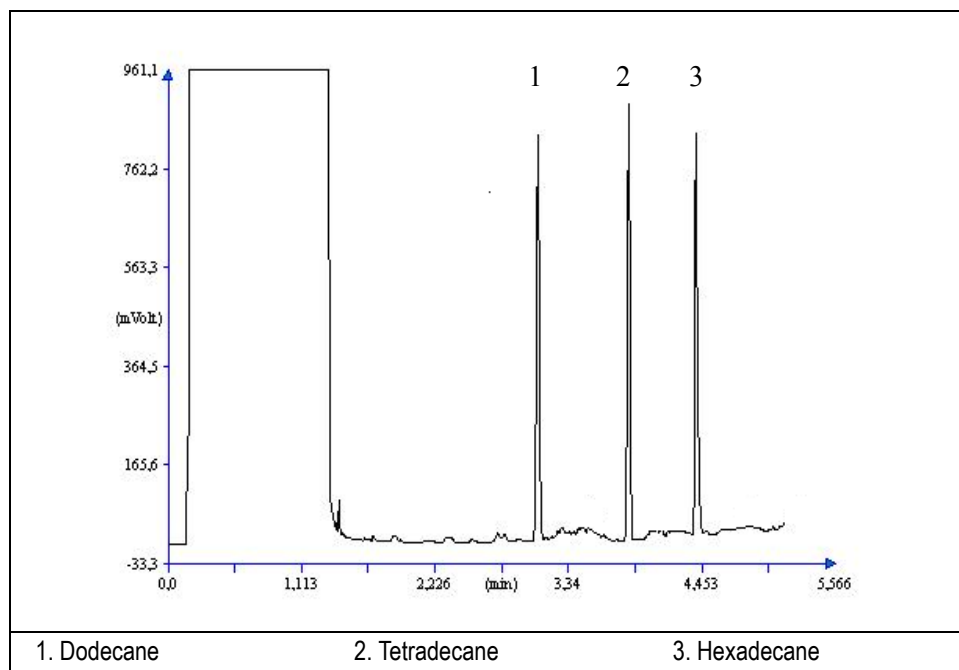





Figure 34-1. FID-LVOC Injection


10. The following criteria indicate successful completion of FID-LVOC checkout.
11. If these criteria are not met, repeat the test.

Table 34-4. FID-LVOC Acceptance Criteria

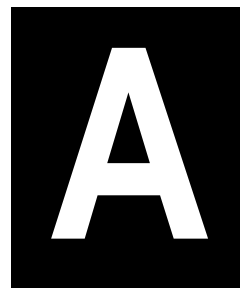
CHROM-CARD			
Acceptance Values	Baseline Parameters	Analog (1V Full Scale)	Digital (10V Full Scale)
	Noise (μV)	< 30	< 300
	Wander (μV)	< 50	< 500
	Drift ($\mu\text{V/h}$)	< 100	< 1 000
	Analytical Results	Analog (1V Full Scale) Area Counts (0.1 μVs)	Digital (10V Full Scale) Area Counts (0.1 μVs)
	Components	> 10 000 000 for each component	> 100 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1	1 ± 0.1
			

CHROMQUEST		
Acceptance Values	Baseline Parameters (1V Full Scale)	
	Noise (μV)	< 30
	Wander (μV)	< 50
	Drift ($\mu\text{V/h}$)	< 100
	Analytical Results (1V Full Scale) - Area Counts (0.01 μVs)	
	Components	> 100 000 000 for each component
	Area Count Ratio Calculated as C12/C16	1 ± 0.1
		

ATLAS	
Baseline Parameters (10V Full Scale)	
Noise (μV)	< 300
Wander (μV)	< 500
Drift ($\mu\text{V}/\text{h}$)	< 1 000
Analytical Results (10V Full Scale) - Area Counts (μVs)	
Components	> 10 000 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

XCALIBUR	
Baseline Parameters (Acquisition Frequency = 10 Hz)	
Noise (Counts)	< 3 000
Wander Counts)	< 5 000
Drift (Counts/h)	< 10 000
Analytical Results Area Counts (Cts*s)	
Components	> 100 000 000 for each component
Area Count Ratio Calculated as C12/C16	1 ± 0.1
	

Analytical Acceptance Comments	
1	When the make-up gas is not used, the acceptance values will result to be 2.5 times lower than the values reported in Table 34-4.
2	When helium is used as make-up gas, the acceptance values will result to be 10 times lower than the values reported in Table 34-4.



Customer Communication

Thermo Fisher Scientific provides comprehensive technical assistance worldwide and is dedicated to the quality of our customer relationships and services.

This appendix also contains a one-page *Reader Survey*. Use this survey to give us feedback on this manual and help us improve the quality of our documentation

How To Contact Us

Use http://www.thermo.com/com/cda/resources/resource_detail/1,,12512,00.html address for products information.

Use <http://www.gc-gcms-customersupport.com/WebPage/Share/Default.aspx> address to contact your local Thermo Fisher Scientific office or affiliate GC-GC/MS Customer Support.

Reader Survey

Product: TRACE GC Ultra
Manual: Standard Operating Procedures Manual
Part No.: 317 092 00

**Please help us improve the quality of our documentation by completing and returning this survey.
 Circle one number for each of the statements below.**

	Strongly Agree	Agree	Neutral	Disagree	Strongly Disagree
The manual is well organized.	1	2	3	4	5
The manual is clearly written.	1	2	3	4	5
The manual contains all the information I need.	1	2	3	4	5
The instructions are easy to follow.	1	2	3	4	5
The instructions are complete.	1	2	3	4	5
The technical information is easy to understand.	1	2	3	4	5
Examples of operation are clear and useful.	1	2	3	4	5
The figures are helpful.	1	2	3	4	5
I was able to install the system using this manual.	1	2	3	4	5

If you would like to make additional comments, please do. (Attach additional sheets if necessary.)

Fax or mail this form to:

Thermo Fisher Scientific S.p.A.
 Strada Rivoltana km 4
 20090 Rodano (MI)
 ITALY
 Fax: 39 02 95059388

This section contains an alphabetical list and descriptions of terms used in this guide and the help diskette. This also includes abbreviations, acronyms, metric prefixes, and symbols.

A

A	ampere
ac	alternating current
ADC	analog-to-digital converter

B

b	bit
B	byte (8 b)
baud rate	data transmission speed in events per second

C

°C	Celsius
CIP	Carriage and Insurance Paid To
cm	centimeter
COC	Cold On-Column Injector
CPU	central processing unit (of a computer)
CSE	Customer Service Engineer
<Ctrl>	control key on the terminal keyboard

D

d	depth
DAC	digital-to-analog converter
dc	direct current
DS	data system

E

ECD	Electron Capture Detector
EMC	electromagnetic compatibility
ESD	electrostatic discharge

F

°F	Fahrenheit
FID	Flame Ionization Detector
FOB	Free on Board
FPD	Flame Photometric Detector
ft	foot

G

g	gram
GC	gas chromatograph
GND	electrical ground

H

<i>h</i>	height
h	hour
harmonic distortion	A high-frequency disturbance that appears as distortion of the fundamental sine wave.
HV	high voltage
Hz	hertz (cycles per second)

I

IEC	International Electrotechnical Commission
-----	---

impulse	See <i>transient</i>
in.	inch
I/O	input/output
K	
k	kilo (10^3 or 1024)
K	Kelvin
kg	kilogram
kPa	kilopascal
L	
l	length
l	liter
LAN	Local Area Network
lb	pound
LED	light-emitting diode
LVOCI	Large Volume On-Column Injector
LVSL	Large Volume Splitless
M	
m	meter (or milli [10^{-3}])
M	mega (10^6)
μ	micro (10^{-6})
min	minute
mL	milliliter
mm	millimeter

Glossary

m/z	mass-to-charge ratio
N	
n	nano (10^{-9})
NPD	Nitrogen Phosphorous Detector
O	
Ω	ohm
P	
p	pico (10^{-12})
Pa	pascal
PCB	printed circuit board
PDD	Pulsed Discharge Detector
PID	Photo Ionization Detector
PN	part number
psi	pounds per square inch
R	
RAM	random access memory
<Return>	<Return> key on the keyboard
RF	radio frequency
ROM	read-only memory
RS-232	industry standard for serial communications
S	
s	second

sag	See <i>surge</i>
slow average	A gradual, long-term change in average RMS voltage level, with typical durations greater than 2 s.
SOP	Standard Operating Procedure
surge	A sudden change in average RMS voltage level, with typical duration between 50 μ s and 2 s.

T

TCD	Thermal Conductivity Detector
transient	A brief voltage surge of up to several thousand volts, with a duration of less than 50 μ s.

U

UFM	Ultra Fast Module
-----	-------------------

V

V	volt
V ac	volts, alternating current
V dc	volts, direct current
VGA	Video Graphics Array

W

<i>w</i>	Width
W	Watt

NOTE The symbol for a compound unit that is a quotient (for example, degrees Celsius per minute or grams per liter) is written with a negative exponent with the denominator. For example:
 $^{\circ}\text{C min}^{-1}$ instead of $^{\circ}\text{C/min}$
 g L^{-1} instead of g/L

A

Acceptance Criteria

- ATLAS 168, 180, 193, 207, 221, 236, 248, 261, 275, 289, 306, 319, 335, 348, 359, 371, 384, 394, 433, 444
- CHROM-CARD 55, 66, 78, 90, 104, 116, 129, 141, 153, 167, 179, 192, 206, 220, 235, 247, 260, 274, 288, 358, 370, 383, 393, 432, 443
- CHROMQUEST 55, 78, 104, 116, 129, 153, 167, 179, 192, 206, 220, 235, 247, 260, 274, 288, 305, 318, 334, 347, 358, 370, 383, 393, 432, 443
- Computing-integrator (e.g. ChromJet) 55, 66, 78, 90, 167, 179, 192, 206, 220, 305, 318, 334, 347, 358, 370, 383, 393
- XCALIBUR 105, 117, 130, 142, 154, 168, 180, 193, 207, 221, 236, 248, 261, 275, 289, 306, 319, 335, 348, 359, 371, 384, 394, 433, 444
- adapter for capillary column 61, 73, 123, 135, 185, 199, 253, 267, 353, 365
- Air 27, 39, 49, 60, 72, 83, 159, 171, 184, 198, 211, 227, 239, 252, 266, 279, 426, 438
- Aldrin 96, 108, 121, 133, 144
- Analytical Acceptance Comments 46, 56, 67, 79, 91, 105, 117, 130, 142, 154, 236, 248, 261, 275, 289, 348, 359, 371, 433, 444
- Analytical Column 38, 58, 70, 82, 120, 132, 144, 182, 196, 210, 250, 264, 278, 322, 350, 362, 376, 424, 436
- Analytical Conditions Required for LVSL Injection 426, 438
- Analytical Conditions Required for On-Column Analysis 171, 239, 309
- Analytical Conditions Required for On-Column Injection 49, 109, 387
- Analytical Conditions Required for PTV Splitless Injection 83, 145, 211, 279, 324
- Analytical Conditions Required for Splitless Injection 39, 60, 72, 83, 97, 122, 134, 145, 159, 184, 198, 211, 227, 252, 266, 279, 341, 352, 364, 377, 425, 437
- ATLAS 433, 444
- Atlas 27, 38, 48, 59, 71, 82, 96, 108, 121, 133, 144, 158, 170, 183, 197, 210, 226, 238, 251, 265, 278, 295, 308, 323, 340, 351, 363, 376, 386, 425, 437
- Azobenzene 158, 165, 170, 177, 183, 190, 197, 204, 210, 218, 226, 238, 251, 265, 278, 294, 301, 302, 303,

304, 308, 314, 315, 316, 317, 322, 330, 331, 332, 333

B

- Base Temperature 39, 49, 60, 72, 83, 97, 109, 122, 134, 145, 159, 171, 184, 198, 211, 227, 239, 252, 266, 279, 296, 309, 324, 341, 352, 364, 377, 387, 426, 438

C

- Carrier Gas 27, 31, 39, 49, 60, 72, 83, 97, 109, 122, 134, 145, 159, 171, 184, 198, 211, 227, 239, 252, 266, 279, 296, 309, 324, 341, 352, 364, 377, 387, 426, 438
- Checkout Using Electron Capture Detector with Cold On- Column Injector 107
- Checkout Using Electron Capture Detector with Split Splitless Injector 95
- Checkout Using FID Tandem 397
- Checkout Using Flame Ionization Detector with Cold On-Column Injector 47, 385, 397, 405, 413
- Checkout Using Flame Ionization Detector with Split Splitless Injector 37, 57, 81, 119, 143, 181, 209, 249, 277, 321, 349, 375, 423, 435
- Checkout Using FPD Tandem 413
- Checkout Using Nitrogen Phosphorus Detector with Cold On-Column Injector 169, 237, 307
- Checkout Using Nitrogen Phosphorus Detector with Split Splitless Injector 157, 225, 293, 339
- Checkout Using NPD Tandem 405
- Checkout Using PDD with OC Injector 385, 397, 405, 413
- Checkout Using PDD with S/SL Injector 375
- ChromQuest 27, 38, 48, 59, 71, 82, 96, 108, 121, 133, 144, 158, 170, 183, 197, 210, 226, 238, 251, 265, 278, 295, 308, 323, 340, 351, 363, 376, 386, 425, 437
- Column Conditioning 29
- Column Conditioning Parameters 31
- Column Evaluation 29
- Column Installation 28
- Column Leak Check 29
- Computing -integrator 38, 48, 59, 71, 82, 96, 108, 121, 133, 144, 158, 170, 183, 197, 210, 226, 238, 251, 265, 278, 295, 308, 323, 340, 351, 363, 376, 386, 425, 437

Index

- Computing-integrator 27
 - Connect the required gas lines 40, 50, 61, 73, 84, 98, 110, 123, 135, 146, 160, 172, 185, 199, 212, 228, 240, 253, 267, 280, 297, 310, 325, 342, 353, 365, 378, 388, 427, 440
 - Connect your data handling 40, 50, 61, 73, 84, 98, 110, 123, 135, 146, 160, 172, 185, 199, 212, 228, 240, 253, 267, 280, 297, 310, 325, 342, 353, 365, 378, 388, 427, 440
 - Constant Filament Temperature 341, 352, 364
 - Constant Septum Purge 39, 83, 97, 145, 159, 211, 227, 279, 296, 324, 341, 377, 426
- ## D
- Data Acquisition 27, 38, 48, 59, 71, 82, 96, 108, 121, 133, 144, 158, 170, 183, 197, 210, 226, 238, 251, 265, 278, 295, 308, 323, 340, 351, 363, 376, 386, 425, 437
 - Data Handling Connections 28
 - Detector 39, 49, 60, 72, 83, 97, 109, 122, 134, 145, 159, 184, 211, 227, 239, 252, 266, 279, 296, 309, 324, 341, 352, 364, 377, 387, 426, 438
 - Detector Connections 29
 - Detector side adapter 58, 120, 182, 250, 350, 362
 - Detector Signal Range 39, 49, 60, 72, 83, 159, 171, 184, 198, 211, 227, 239, 252, 266, 279, 296, 309, 324, 377, 387, 426, 438
 - Detector UV Lamp 294, 308, 322
 - Digital Chrom-Card 27, 38, 48, 59, 71, 82, 96, 108, 133, 144, 158, 170, 183, 197, 238, 251, 265, 278, 295, 308, 323, 340, 363, 376, 386, 425, 437
 - Dodecane 38, 48, 59, 71, 82, 340, 351, 363, 376, 386, 392, 425, 437
 - Drift 45, 46, 55, 56, 66, 67, 78, 79, 90, 91, 104, 105, 116, 117, 129, 130, 141, 142, 153, 154, 167, 168, 179, 180, 192, 193, 206, 207, 220, 221, 235, 236, 247, 248, 260, 261, 274, 275, 288, 289, 305, 306, 318, 319, 334, 335, 347, 348, 358, 359, 370, 371, 383, 384, 393, 394, 432, 433, 443, 444
- ## E
- ECD 127, 139, 151
 - ECD temperature 97, 109, 122, 134, 145
 - ECD-COC Checkout in On-Column Mode 111
 - ECD-COC Parts Referenced 108
 - ECD-OCI Analytical Conditions 109
 - ECD-OCI Checkout in On-Column Mode 111
 - ECD-OCI Parts Referenced 108
 - ECD-On-Column Injection 114
 - ECD-PTV Checkout in PTV Splitless Mode 147
 - ECD-S/SL Analytical Conditions 97
 - ECD-S/SL Checkout in Splitless Mode 99
 - ECD-S/SL Parts Referenced 96
 - ECD-Splitless Injection 102
 - ECD-SSL Analytical Conditions 97
 - ECD-SSL Checkout in Splitless Mode 99
 - ECD-SSL Injection 102
 - ECD-SSL Parts Referenced 96
- ## F
- FID 65, 77, 89
 - FID Tandem
 - Checkout Overview 399
 - Connecting Test Column 401
 - Important Considerations 400
 - FID-ECD SOPs Reference 399
 - FID-LVSL Analytical Conditions 426, 438
 - FID-LVSL Checkout in Splitless Mode 428, 441
 - FID-LVSL Parts Referenced 424, 436
 - FID-OC Parts Referenced 48, 386
 - FID-OCI Analytical Conditions 49
 - FID-OCI Checkout in On-Column Mode 51, 389
 - FID-OCI Checkout in On-Column Mode 327 385, 397, 413
 - FID-OCI Parts Referenced 48, 386
 - FID-On-Column Injection 54, 392
 - FID-PTV Checkout in PTV Splitless Mode 85
 - FID-S/SL
 - Analytical Acceptance Criteria 45, 55, 66, 78, 90, 104, 116, 129, 141, 153, 167, 179, 192, 206, 220, 235, 247, 260, 274, 288, 305, 318, 334, 347, 358, 370, 383, 393, 432, 443
 - FID-S/SL Acceptance Criteria
 - ATLAS 46, 56, 67, 79, 91
 - CHROM-CARD 45
 - CHROMQUEST 45
 - Computing-integrator (e.g. ChromJet) 45
 - XCALIBUR 46, 56, 67, 79, 91
 - FID-S/SL Analytical Conditions 39, 60, 72, 83, 122, 134, 145, 184, 198, 211, 252, 266, 279, 324, 352, 364, 377, 426, 438

- FID-S/SL Checkout in Splitless Mode 41, 85, 123, 147, 185, 213, 253, 281, 326, 353, 379, 428, 441
- FID-S/SL Parts Referenced 38, 58, 70, 82, 120, 132, 144, 182, 196, 210, 250, 264, 278, 322, 350, 362, 424, 436
- FID-Splitless Injection 44, 65, 77, 89, 127, 139, 151, 190, 204, 218, 357, 369, 431, 442
- FID-SSL Analytical Conditions 39, 60, 72, 83, 122, 134, 145, 184, 198, 211, 252, 266, 279, 324, 352, 364, 377, 426, 438
- FID-SSL Checkout in Splitless Mode 41, 123, 147, 185, 213, 253, 281, 326, 353, 379, 428, 441
- Filament Temperature limit 341, 352, 364
- Filament Voltage 341, 352, 364
- Final Temperature 31, 39, 49, 60, 72, 83, 97, 109, 134, 145, 159, 171, 184, 198, 211, 227, 239, 252, 266, 279, 296, 309, 324, 341, 352, 364, 377, 387, 426, 438
- Final Time 31, 39, 49, 60, 72, 83, 97, 109, 134, 145, 159, 171, 184, 198, 211, 227, 239, 252, 266, 279, 296, 309, 324, 341, 352, 364, 377, 387, 426, 438
- Flame Ionization Detector 35, 155
- FPD Tandem
 - Checkout Overview 415
 - Connecting Test Column 417
 - Important Considerations 416
- FPD Temperature 227, 239, 252, 266, 279
- FPD-ECD SOPs Reference 415
- FPD-On-Column Injection with 394 nm 245
- FPD-On-Column Injection with 526 nm filter 244
- FPD-PKD Injection with 394 nm 258
- FPD-PKD Injection with 526 nm filter 257
- FPD-PPKD Injection with 394 nm 272
- FPD-PPKD Injection with 526 nm filter 271
- FPD-PTV Checkout in PTV Splitless Mode 281
- FPD-PTV Injection with 394 nm filter 286
- FPD-PTV Injection with 526 nm filter 285
- FPD-Splitless Injection with 394 nm 233, 245, 258
- FPD-Splitless Injection with 526 nm filte 232, 244, 257, 271
- Fuel Gases 27
- G**
- Gain 341, 352, 364
- Gas Supply Connections 28
- Gases 27, 38, 39, 48, 49, 59, 60, 71, 72, 82, 83, 96, 97, 108, 109, 121, 122, 133, 134, 144, 145, 158, 159, 170, 171, 183, 184, 197, 198, 210, 211, 226, 227, 238, 239, 251, 252, 265, 266, 278, 279, 294, 296, 308, 309, 322, 324, 340, 341, 351, 352, 363, 364, 376, 377, 386, 387, 425, 426, 437, 438
- General Overview 25
- Getting Started 28
- Glass Liner 38, 58, 70, 82, 96, 120, 132, 144, 158, 182, 196, 210, 226, 250, 264, 294, 340, 350, 362, 376, 424
- Glass Liner Installation 28
- Graphite Ferrule 48, 96, 108, 158, 170, 226, 238, 294, 308, 340, 386
- Graphite ferrule for 0.32 mm ID Column 38, 58, 70, 82, 120, 132, 144, 182, 196, 210, 250, 264, 278, 322, 350, 362, 376, 424, 436
- Graphite seal for glass liner 38, 58, 70, 82, 120, 132, 144, 158, 182, 196, 210, 226, 250, 264, 278, 294, 322, 340, 350, 362, 376
- Graphite seal glass liner 96
- H**
- Helium 27, 39, 49, 60, 72, 83, 97, 109, 122, 134, 145, 159, 171, 184, 198, 211, 227, 239, 252, 266, 279, 296, 309, 324, 341, 352, 364, 377, 387, 426, 438
- Helium Purifier 376, 386
- helium purifier
 - Verify 378
- Hexadecane 38, 48, 59, 71, 82, 340, 351, 363, 376, 386, 392, 425, 437
- High Current 296, 309, 324
- High voltage mode 252, 266, 279
- Hydrogen 27, 39, 49, 60, 72, 83, 159, 171, 184, 198, 211, 227, 239, 252, 266, 279, 426, 438
- I**
- Initial Temperature 31, 39, 49, 60, 72, 83, 97, 109, 122, 134, 145, 159, 171, 184, 198, 211, 227, 239, 252, 266, 279, 296, 309, 324, 341, 352, 364, 377, 387, 426, 438
- Initial Time 31, 39, 49, 60, 72, 83, 97, 109, 122, 134, 145, 159, 171, 184, 198, 211, 227, 239, 252, 266, 279, 296, 309, 324, 341, 352, 364, 377, 387, 426, 438
- Inject Temp 83, 145, 211, 279, 324
- Inject Time 83, 145, 211, 279, 324

Index

- Injected Volume 39, 49, 60, 72, 83, 97, 109, 122, 134, 145, 159, 171, 184, 198, 211, 227, 239, 252, 266, 279, 296, 309, 324, 341, 352, 364, 377, 387, 426, 438
- Injection side adapte 182
- Injection side adapter 58, 70, 120, 132, 196, 250, 350, 362
- Injector 31, 39, 49, 60, 72, 83, 97, 109, 122, 134, 145, 159, 171, 184, 198, 211, 227, 239, 252, 266, 279, 296, 309, 324, 341, 352, 364, 377, 387, 426, 438
- Install the analytical column 40, 61, 73, 84, 98, 123, 135, 146, 185, 199, 212, 253, 267, 280, 325, 353, 365, 378, 427, 440
- Install the test column 50, 61, 73, 110, 123, 135, 160, 172, 185, 199, 228, 240, 253, 297, 310, 342, 353, 365, 388
- Interferential Filter 226, 238, 250, 264, 278
- L**
- Lamp Current 296, 309, 324, 341, 352, 364
- Leak Test 30, 98, 160, 212, 228, 280, 297, 325, 342
- Lindane 96, 108, 121, 133, 144
- Liner Seal 38, 58, 70, 82, 96, 120, 132, 144, 158, 182, 196, 210, 226, 250, 264, 278, 294, 322, 340, 350, 362, 376
- M**
- Make-up 97, 109, 159, 171, 296, 309, 324, 341, 352, 364
- Make-up Gas 27, 39, 49, 60, 72, 83, 122, 134, 145, 184, 198, 211, 426, 438
- manual leak check 61, 73, 123, 135, 185, 253, 267, 353, 365
- N**
- Negative Polarity 341, 352, 364
- Nitrogen 27, 39, 49, 60, 72, 83, 97, 109, 122, 134, 145, 159, 171, 184, 198, 211, 296, 309, 324, 341, 352, 364, 426, 438
- Noise 45, 46, 55, 56, 66, 67, 78, 79, 90, 91, 104, 105, 116, 117, 129, 130, 141, 142, 153, 154, 167, 168, 179, 180, 192, 193, 206, 207, 220, 221, 235, 236, 247, 248, 260, 261, 274, 275, 288, 289, 305, 306, 318, 319, 334, 335, 347, 348, 358, 359, 370, 371, 383, 384, 393, 394, 432, 433, 443, 444
- NPD 190, 204, 218
- NPD Tandem
Checkout Overview 407
- Important Considerations 408
- NPD Tandem Checkout
Example 409
- NPD-COC Analytical Conditions 171, 239, 309
- NPD-COC Checkout in On-Column Mode 173, 241, 311
- NPD-COC Parts Referenced 170, 238, 308
- NPD-ECD SOPs Reference 407
- NPD-OCI Analytical Conditions 171, 239, 309
- NPD-OCI Checkout in On-Column Mode 173, 241, 311
- NPD-OCI Parts Referenced 170, 238, 308
- NPD-On-Column Injection 177
- NPD-On-column Injection 177
- NPD-PTV Checkout in PTV Splitless Mode 213
- NPD-S/SL Analytical Conditions 159, 227, 296, 341
- NPD-S/SL Checkout in Splitless Mode 161, 229, 298, 343
- NPD-S/SL Parts Referenced 158, 226, 294, 340
- NPD-Splitless Injection 165, 232, 244, 257, 271, 301, 302, 303, 304, 314, 315, 316, 317, 330, 331, 332, 346
- NPD-SSL Analytical Conditions 159, 227, 296, 341
- NPD-SSL Checkout in Splitless Mode 161, 229, 298, 343
- NPD-SSL Injection 165, 232, 244, 257, 271, 301, 302, 303, 304, 314, 315, 316, 317, 330, 331, 332, 346
- NPD-SSL Parts Referenced 158, 226, 294, 340
- O**
- Octadecan 165, 177, 190, 204, 218, 301, 302, 303, 304, 314, 315, 316, 317, 330, 331, 332, 333
- Octadecane 158, 170, 183, 197, 210, 226, 238, 251, 265, 278, 294, 308, 322
- Operating Mode 39, 60, 72, 83, 97, 122, 134, 145, 159, 184, 198, 211, 227, 252, 266, 279, 296, 324, 341, 352, 364, 377, 426, 438
- Operating Procedure 28
- Oven Program 31, 39, 49, 60, 72, 83, 97, 109, 122, 134, 145, 159, 171, 184, 198, 211, 227, 239, 252, 266, 279, 296, 309, 324, 341, 352, 364, 377, 387, 426, 438
- P**
- Parathion methyl 158, 170, 183, 197, 210, 226, 238, 251, 265, 278, 294, 308, 322

- Parts Referenced 27, 48, 96, 108, 158, 170, 226, 238, 294, 308, 340, 386, 424, 436
- PDD-OCI Analytical Conditions 387
- PDD-OCI Checkout in On-Column Mode 389
- PDD-OCI Parts Referenced 386
- PDD-On-Column Injection 392
- PDD-S/SL
- Analytical Conditions 377
 - Checkout in Splitless Mode 379
 - Parts Referenced 376
- PDD-Splitless Injection 382
- Perform Column Evaluation 73, 135, 199, 365
- Perform Leak Test and Column Evaluation 40, 50, 84, 98, 110, 146, 160, 172, 212, 228, 240, 280, 297, 310, 325, 342, 378, 388, 427, 440
- PID-On-Column Injection with 10.6 eV or 11.8 eV UV Lamp 316, 317
- PID-On-Column Injection with 10.6 eV UV Lamp 316
- PID-On-Column Injection with 11.8 eV UV Lamp 317
- PID-On-Column Injection with 9.6 eV UV Lamp 315
- PID-PKD Injection with 10.6 eV or 11.8 eV UV Lamp 332
- PID-PKD Injection with 8.4 eV UV Lamp 330
- PID-PKD Injection with 9.6 eV UV Lamp 331
- PID-PPKD Injection with 10.6 eV or 11.8 eV UV Lamp 332
- PID-PPKD Injection with 8.4 eV UV Lamp 330
- PID-PPKD Injection with 9.6 eV UV Lamp 331
- PID-PTV Injection with 10.6 eV or 11.8 eV UV Lamp 332
- PID-PTV Injection with 10.6 eV UV Lamp 332
- PID-PTV Injection with 11.8 eV UV Lamp 333
- PID-PTV Injection with 8.4 eV UV Lamp 330
- PID-PTV Injection with 9.6 eV UV Lamp 331
- PID-Splitless Injection with 10.6 eV or 11.8 eV UV Lamp 303, 304, 316, 317, 332
- PID-Splitless Injection with 10.6 eV UV Lamp 303
- PID-Splitless Injection with 11.8 eV UV Lamp 304
- PID-Splitless Injection with 8.4 eV UV Lamp 301, 302, 303, 304, 314, 315, 316, 317, 330, 331, 332
- PID-Splitless Injection with 9.6 eV UV Lamp 302, 303, 304, 315, 316, 317, 331, 332
- Polarizer voltage 159, 171, 184, 198, 211, 239, 252
- Press 70
- Press-fit cconnections 70
- Press-fit connections 58, 70, 120, 132, 182, 196, 250, 350, 362
- Pulse Amplitude 97, 109, 122, 134, 145
- Pulse generator 380, 390
- Pulse Width 97, 109, 122, 134, 145
- R**
- Ramp 1 31, 39, 49, 60, 72, 83, 97, 109, 134, 145, 159, 171, 184, 198, 211, 227, 239, 252, 266, 279, 296, 309, 324, 341, 352, 364, 377, 387, 426, 438
- Recommended 40, 73, 135, 146, 267, 280, 325, 365, 378, 427, 440
- Recommended Initial Operations 40, 50, 73, 83, 98, 110, 122, 135, 146, 160, 172, 184, 198, 212, 228, 240, 252, 267, 280, 310, 325, 342, 365, 378, 388, 427, 440
- Reference Current 97, 109, 122, 134, 145
- Reference Gas 341, 352, 364
- Registration and Revision Number 26
- Replace the glass liner 98, 160, 228, 297, 342
- Replace the glass liner. 40, 61, 73, 84, 123, 135, 146, 185, 199, 212, 253, 267, 280, 325, 353, 365, 378
- Replace the septum 40, 61, 73, 84, 98, 123, 135, 146, 160, 185, 199, 212, 228, 253, 267, 280, 297, 325, 342, 353, 365, 378, 427
- S**
- Scope 26, 38, 48, 58, 70, 82, 96, 108, 120, 132, 144, 158, 170, 182, 196, 210, 226, 238, 250, 264, 278, 294, 308, 322, 340, 350, 362, 376, 386, 398, 406, 414, 424, 436
- Secondary Cooling 49, 109, 387
- Secondary cooling 171, 239, 309
- Septum 38, 58, 70, 82, 96, 120, 132, 144, 158, 182, 196, 210, 226, 250, 264, 278, 294, 322, 340, 350, 362, 376, 424, 436
- Septum Installation 28
- Shealth Gas 296, 309, 324, 341, 352, 364
- SOPs Parts Referenced 27
- Source Current 159, 171, 184, 198, 211, 227, 239, 252, 266, 279
- Split Flow 39, 83, 97, 145, 159, 211, 227, 279, 296, 324, 341, 377, 426

Index

- Splitless Time 39, 83, 97, 145, 159, 211, 227, 279, 296,
324, 341, 377, 426
- Standard 26
- Standard Operating Procedures 26
- Syringe 27, 38, 48, 58, 70, 82, 96, 108, 120, 132, 144, 158,
170, 182, 196, 210, 226, 238, 250, 264, 278, 294,
308, 322, 340, 350, 362, 376, 386, 424, 436
- T**
- TCD 357, 369
- Temperature 39, 60, 72, 97, 122, 134, 159, 184, 198, 227,
252, 266, 296, 341, 352, 364, 377, 426
- Test Column 27, 48, 96, 108, 158, 170, 226, 238, 294,
308, 340, 386
- Test Mixture 38, 48, 59, 71, 82, 96, 108, 121, 133, 144,
158, 170, 183, 197, 210, 226, 238, 251, 265, 278,
294, 308, 322, 340, 351, 363, 376, 386, 425, 437
- Test Mixtures 27
- Tetradecane 38, 48, 59, 71, 82, 340, 351, 363, 376, 386,
392, 425, 437
- Transfer ram 83, 145, 279, 324
- Transfer ramp 83, 145, 211, 279, 324
- Transfer Temperature 83, 145, 211, 279, 324
- Transfer time 83, 145, 211, 279, 324
- W**
- Wander 45, 46, 55, 56, 66, 67, 78, 79, 90, 91, 104, 105,
116, 117, 129, 130, 141, 142, 153, 154, 167, 168,
179, 180, 192, 193, 206, 207, 220, 221, 235, 236,
247, 248, 260, 261, 274, 275, 288, 289, 305, 306,
318, 319, 334, 335, 347, 348, 358, 359, 370, 371,
383, 384, 393, 394, 432, 433, 443, 444
- X**
- Xcalibur 27, 38, 48, 59, 71, 82, 96, 108, 121, 133, 144,
158, 170, 183, 197, 210, 226, 238, 251, 265, 278,
295, 308, 323, 340, 351, 363, 376, 386, 425, 437

Operating Procedures

Test Column Conditioning	31
FID-S/SL Checkout in Splitless Mode	41
FID-OCI Checkout in On-Column Mode	51
FID-PKD Checkout	62
FID-PPKD Checkout	74
FID-PTV Checkout in PTV Splitless Mode	85
ECD-S/SL Checkout in Splitless Mode	99
ECD-OCI Checkout in On-Column Mode	111
ECD-PKD Checkout	124
ECD-PPKD Checkout	136
ECD-PTV Checkout in PTV Splitless Mode	147
NPD-S/SL Checkout in Splitless Mode	161
NPD-OCI Checkout in On-Column Mode	173
NPD-PKD Checkout	186
NPD-PPKD Checkout	200
NPD-PTV Checkout in PTV Splitless Mode	213
FPD-S/SL Checkout in Splitless Mode	229
FPD-OCI Checkout in On-Column Mode	241
FPD-PKD Checkout	254
FPD-PPKD Checkout	268
FPD-PTV Checkout in PTV Splitless Mode	281
PID-S/SL Checkout in Splitless Mode	298
PID-OCI Checkout in On-Column Mode	311
PID-PTV Checkout in PTV Splitless Mode	326
TCD-S/SL Checkout in Splitless Mode	343
TCD-PKD Checkout	354
TCD-PPKD Checkout	366
PDD-S/SL Checkout in Splitless Mode	379
PDD-OCI Checkout in On-Column Mode	389
Example of FID Tandem Checkout	401
Example of NPD Tandem Checkout	409
Example of FPD Tandem Checkout	417
FID-LVSL Checkout in Splitless Mode	428
FID-LVOC Checkout in LVOC Mode	441

Figures

Figure 2-1.	FID-Splitless Injection.....	44
Figure 3-1.	FID-On-Column Injection.....	54
Figure 4-1.	FID-Packed Injection.....	65
Figure 5-1.	FID-PPKD Injection.....	77
Figure 6-1.	FID-PTV Injection.....	89
Figure 7-1.	ECD-Splitless Injection.....	102
Figure 8-1.	ECD-On-Column Injection.....	114
Figure 9-1.	ECD-PKD Injection.....	127
Figure 10-1.	ECD-PPKD Injection.....	139
Figure 11-1.	ECD-PTV Injection.....	151
Figure 12-1.	NPD-Splitless Injection.....	165
Figure 13-1.	NPD-On-Column Injection.....	177
Figure 14-1.	NPD-Packed Injection.....	190
Figure 15-1.	NPD-PPKD Injection.....	204
Figure 16-1.	NPD-PTV Injection.....	218
Figure 17-1.	FPD-Splitless Injection with 526 nm filter.....	232
Figure 17-2.	FPD-Splitless Injection with 394 nm.....	233
Figure 18-1.	FPD-On-Column Injection with 526 nm filter.....	244
Figure 18-2.	FPD-On-Column Injection with 394 nm.....	245
Figure 19-1.	FPD-PKD Injection with 526 nm filter.....	257
Figure 19-2.	FPD-PKD Injection with 394 nm.....	258
Figure 20-1.	FPD-PPKD Injection with 526 nm filter.....	271
Figure 20-2.	FPD-PPKD Injection with 394 nm.....	272
Figure 21-1.	FPD-PTV Injection with 526 nm filter.....	285
Figure 21-2.	FPD-PTV Injection with 394 nm filter.....	286
Figure 22-1.	PID-Splitless Injection with 8.4 eV UV Lamp.....	301
Figure 22-2.	PID-Splitless Injection with 9.6 eV UV Lamp.....	302
Figure 22-3.	PID-Splitless Injection with 10.6 eV UV Lamp.....	303
Figure 22-4.	PID-Splitless Injection with 11.8 eV UV Lamp.....	304
Figure 23-1.	PID-On-Column Injection with 8.4 eV UV Lamp.....	314
Figure 23-2.	PID-On-Column Injection with 9.6 eV UV Lamp.....	315
Figure 23-3.	PID-On-Column Injection with 10.6 eV UV Lamp.....	316
Figure 23-4.	PID-On-Column Injection with 11.8 eV UV Lamp.....	317
Figure 24-1.	PID-PTV Injection with 8.4 eV UV Lamp.....	330
Figure 24-2.	PID-PTV Injection with 9.6 eV UV Lamp.....	331
Figure 24-3.	PID-PTV Injection with 10.6 eV UV Lamp.....	332
Figure 24-4.	PID-PTV Injection with 11.8 eV UV Lamp.....	333

Figure 25-1. TCD-Splitless Injection	346
Figure 26-1. TCD-Packed Injection	357
Figure 27-1. TCD-PPKD Injection	369
Figure 28-1. PDD-Splitless Injection	382
Figure 29-1. PDD-On-Column Injection	392
Figure 30-1. FID Tandem.....	398
Figure 31-1. NPD Tandem	406
Figure 32-1. FPD Tandem.....	414
Figure 33-1. FID-LV Splitless Injection	431
Figure 34-1. FID-LVOC Injection	442



Tables

Table 1-1.	SOPs Parts Referenced	27
Table 1-2.	Column Conditioning Parameters	31
Table 2-1.	FID-S/SL Parts Referenced	38
Table 2-2.	FID-S/SL Analytical Conditions	39
Table 2-3.	FID-S/SL Acceptance Criteria	45
Table 3-1.	FID-OCI Parts Referenced	48
Table 3-2.	FID-OCI Analytical Conditions	49
Table 3-3.	FID-OCI Acceptance Criteria	55
Table 4-1.	FID-PKD Parts Referenced	58
Table 4-2.	FID-PKD Analytical Conditions	60
Table 4-3.	FID-PKD Acceptance Criteria	66
Table 5-1.	FID-PPKD Parts Referenced	70
Table 5-2.	FID-PPKD Analytical Conditions	72
Table 5-3.	FID-PPKD Acceptance Criteria	78
Table 6-1.	FID-PTV Parts Referenced	82
Table 6-2.	FID-PTV Analytical Conditions	83
Table 6-3.	FID-PPKD Acceptance Criteria	90
Table 7-1.	ECD-S/SL Parts Referenced	96
Table 7-2.	ECD-S/SL Analytical Conditions	97
Table 7-3.	ECD-S/SL Acceptance Criteria	104
Table 8-1.	ECD-OCI Parts Referenced	108
Table 8-2.	ECD-OCI Analytical Conditions	109
Table 8-3.	ECD-OCI Acceptance Criteria	116
Table 9-1.	ECD-PKD Parts Referenced	120
Table 9-2.	ECD-PKD Analytical Conditions	122
Table 9-3.	ECD-PKD Acceptance Criteria	129
Table 10-1.	ECD-PPKD Parts Referenced	132
Table 10-2.	ECD-PPKD Analytical Conditions	134
Table 10-3.	ECD-PPKD Acceptance Criteria	141
Table 11-1.	ECD-PTV Parts Referenced	144
Table 11-2.	ECD-PTV Analytical Conditions	145
Table 11-3.	ECD-PTV Acceptance Criteria	153
Table 12-1.	NPD-S/SL Parts Referenced	158
Table 12-2.	NPD-S/SL Analytical Conditions	159
Table 12-3.	NPD-S/SL Acceptance Criteria	167
Table 13-1.	NPD-OCI Parts Referenced	170
Table 13-2.	NPD-OCI Analytical Conditions	171

Table 13-3.	NPD-OCI Acceptance Criteria	179
Table 14-1.	NPD-PKD Parts Referenced.....	182
Table 14-2.	NPD-PKD Analytical Conditions.....	184
Table 14-3.	NPD-PKD Acceptance Criteria.....	192
Table 15-1.	NPD-PPKD Parts Referenced	196
Table 15-2.	NPD-PPKD Analytical Conditions	198
Table 15-3.	NPD-PPKD Acceptance Criteria.....	206
Table 16-1.	NPD-PTV Parts Referenced.....	210
Table 16-2.	NPD-PTV Analytical Conditions.....	211
Table 16-3.	NPD-PTV Acceptance Criteria	220
Table 17-1.	FPD-S/SL Parts Referenced	226
Table 17-2.	FPD-S/SL Analytical Conditions	227
Table 17-3.	FPD-S/SL Acceptance Criteria.....	235
Table 18-1.	FPD-OCI Parts Referenced	238
Table 18-2.	FPD-OCI Analytical Conditions	239
Table 18-3.	FPD-OCI Acceptance Criteria.....	247
Table 19-1.	FPD-PKD Parts Referenced	250
Table 19-2.	FPD-PKD Analytical Conditions	252
Table 19-3.	FPD-PKD Acceptance Criteria.....	260
Table 20-1.	FPD-PPKD Parts Referenced	264
Table 20-2.	FPD-PPKD Analytical Conditions.....	266
Table 20-3.	FPD-PPKD Acceptance Criteria	274
Table 21-1.	FPD-PTV Parts Referenced.....	278
Table 21-2.	FPD-PTV Analytical Conditions.....	279
Table 21-3.	FPD-PTV Acceptance Criteria	288
Table 22-1.	PID-S/SL Parts Referenced	294
Table 22-2.	PID-S/SL Analytical Conditions	296
Table 22-3.	PID-S/SL Acceptance Criteria	305
Table 23-1.	PID-OCI Parts Referenced	308
Table 23-2.	PID-OCI Analytical Conditions	309
Table 23-3.	PID-OCI Acceptance Criteria.....	318
Table 24-1.	PID-PTV Parts Referenced.....	322
Table 24-2.	PID-PTV Analytical Conditions.....	324
Table 24-3.	PID-PTV Acceptance Criteria.....	334
Table 25-1.	TCD-S/SL Parts Referenced.....	340
Table 25-2.	TCD-S/SL Analytical Conditions.....	341
Table 25-3.	TCD-S/SL Acceptance Criteria.....	347
Table 26-1.	TCD-PKD Parts Referenced.....	350

Table 26-2.	TCD-PKD Analytical Conditions.....	352
Table 26-3.	TCD-PKD Acceptance Criteria.....	358
Table 27-1.	TCD-PPKD Parts Referenced.....	362
Table 27-2.	TCD-PPKD Analytical Conditions.....	364
Table 27-3.	TCD-PPKD Acceptance Criteria.....	370
Table 28-1.	PDD-S/SL Parts Referenced.....	376
Table 28-2.	PDD-S/SL Analytical Conditions.....	377
Table 28-3.	PDD-S/SL Acceptance Criteria.....	383
Table 29-1.	PDD-OCI Parts Referenced.....	386
Table 29-2.	PDD-OCI Analytical Conditions.....	387
Table 29-3.	PDD-OCI Acceptance Criteria.....	393
Table 30-1.	FID-ECD SOPs Reference.....	399
Table 31-1.	NPD-ECD SOPs Reference.....	407
Table 32-1.	FPD-ECD SOPs Reference.....	415
Table 33-1.	FID-LVSL Parts Referenced.....	424
Table 33-2.	FID-LVSL Analytical Conditions.....	426
Table 33-3.	FID-LVSL Acceptance Criteria.....	432
Table 34-1.	FID-LVOC Parts Referenced.....	436
Table 34-2.	FID-LVOC Analytical Conditions for the GC.....	438
Table 34-3.	FID-LVOC Analytical Conditions for the TriPlus Autosampler.....	439
Table 34-4.	FID-LVOC Acceptance Criteria.....	443

