



PRODUCT MANUAL

IonPac[®] TRACE METAL CONCENTRATOR (TMC-1) COLUMN

Now sold under the
Thermo Scientific brand

Thermo
SCIENTIFIC

 **DIONEX**

IC | HPLC | MS | EXTRACTION | PROCESS | AUTOMATION

PRODUCT MANUAL

for the

IONPAC[®] TRACE METAL CONCENTRATOR COLUMN (TMC-1)

(P/N 049000)

© DIONEX Corporation, 2004

Document No. 034255
Revision 06
18 November 2004

TABLE OF CONTENTS

SECTION 1 - INTRODUCTION 3

SECTION 2 - OPERATION 4

2.1 Chemicals Required 4

2.2 Solutions Required 4

2.3 Concentration Methods 4

SECTION 3 - TROUBLESHOOTING GUIDE 5

3.1 High Back Pressure from a Contaminated Inlet Bed Support 5

3.2 High Background, or Noise 5

3.3 Poor Peak Shape 6

SECTION 1 - INTRODUCTION

The Trace Metal Concentrator Column (TMC-1) is used in Chelation Ion Chromatography. This manual describes the operation of the TMC-1. The following DIONEX Technical Note describes an important application which requires the TMC-1.

Technical

Note # Description

25 Determination of Transition Metals in Complex Matrices by Chelation Ion Chromatography™

The TMC-1 contains a 17- μm , fully sulfonated, high capacity cation-exchange resin and has an ion-exchange capacity of 0.3 μeq /column. The TMC-1 has high affinity for alkaline-earth, transition and lanthanide metals. In normal operation, however, Group I and II sample cations are removed in the MetPac CC-1 concentrator column. The TMC-1 is specifically designed for interfacing the MetPac CC-1 Column (P/N 042156) with the IonPac CS5 Analytical Column (P/N 037028), or the IonPac CS5A Analytical Column (P/N 046100 or 052516).

The operating back pressure of the TMC-1 should be less than 300 psi under normal operating conditions. Typically, the TMC-1 is used in conjunction with the Chelation IC System where the back pressure should not exceed 1,500 psi.

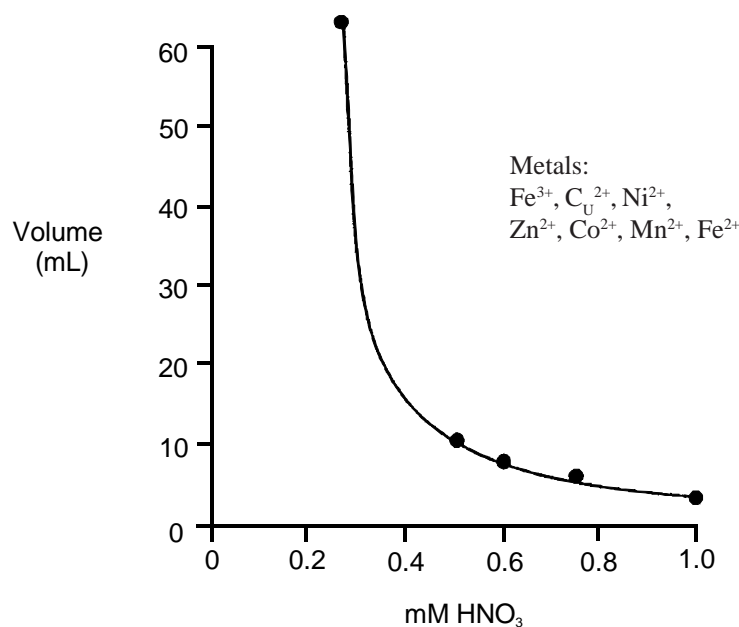


Figure 1

Retention Volume vs Concentration of HNO₃ for Transition Metals on TMC-1

SECTION 2 - OPERATION

2.1 Chemicals Required

The chemicals and water required to prepare the reagents and eluents should be of the highest purity available. Use deionized water with a specific resistance of 18.2 megohm-cm. Prepared reagents can be purchased from DIONEX. See Technical Note No. 25 for reagent ordering information and preparation procedures.

2.2 Solutions Required

Table 1
Chelation Concentration Reagents

	1 liter	6 Pack
2 M Nitric Acid	P/N 033442	P/N 033443
2 M Ammonium Acetate	P/N 033440	P/N 033441
0.1 M Ammonium Nitrate	P/N 033444	P/N 033445 (for CIC only)

2.3 Concentration Methods

The TMC-1 is used to interface the MetPac CC-1 with the IonPac CS5 or CS5A Analytical Column when performing Chelation IC (see Technical Note No. 25 for operation of the Chelation IC System).

Install the TMC-1 as indicated in Technical Note No. 25, Figure 3, "Schematic of Chelation IC System 1," or Figure 4, "Schematic of Chelation IC System 2."

The concentrated transition metals can be eluted from the TMC-1 using nitric acid. Use DIONEX 2.0 M Nitric Acid Chelation Concentration Reagent (P/N 033442 or 033443) for the most reliable results. The elution volumes of various transition metals, as a function of nitric acid concentration, is given in Figure 1, "Retention Volume vs. Concentration of HNO₃ for Transition Metals on TMC-1." The TMC-1 is compatible with most acids and bases up to a concentration of 3.0M.

Due to the relatively high capacity of the TMC-1, it should not be used in place of other cation concentrators such as the IonPac CG2 or CG3.

The recommended cleaning procedure is a 15 minute rinse with 0.02 M oxalic acid at 2.0 mL/min. This cleaning procedure effectively removes both ionic and colloidal metal contamination.

SECTION 3 - TROUBLESHOOTING GUIDE

The purpose of the Troubleshooting Guide is to help you solve operating problems that may arise while using the Trace Metal Concentrator Column (TMC-1). For more information on problems that originate with the Sample Concentration Module or the Ion Chromatograph, refer to the Troubleshooting Guide in the appropriate operator's manual. If you cannot solve the problem on your own, call the DIONEX Office nearest you (see, "DIONEX Worldwide Offices").

3.1 High Back Pressure from a Contaminated Inlet Bed Support

If the TMC-1 displays high back pressure, the bed support in the column inlet may be contaminated. Follow the instructions below to change the bed support assembly using one of the two spare bed support assemblies included in the ship kit provided with the column.

- a) Disconnect the column from the system.
- b) Carefully unscrew the inlet (top) column end fitting using two open-end wrenches.
- c) Remove the old bed support. Turn the end fitting over and tap it against a benchtop or other hard, flat surface to remove the bed support and seal assembly. If the bed support must be pried out of the end fitting, use a sharp pointed object such as a pair of tweezers, but be careful that you do not scratch the walls of the end fitting. Discard the old assembly.
- d) Place a new bed support assembly in the end fitting. Use the end of the column to carefully start the bed support assembly into the end fitting.

Bed Support Assembly	P/N042955
Seal Washer	P/N042956
Bed Support	P/N053889
End Fitting	P/N042367

- e) Screw the end fitting back onto the column. Tighten it fingertight and then using two open-end wrenches, tighten it an additional 1/4 turn (25 in/lb). Tighten further only if leaks are observed.

NOTE: *If any of the column packing becomes lodged between the end of the column and the bed support washer assembly, no amount of tightening will seal the column. Make sure that the washer and the end of the column are clean before screwing the end fitting back onto the column.*

- f) Reconnect the column to the system and resume operation.

3.2 High Background, or Noise

Normally, problems such as high background, noise or baseline instability will not be attributable to the TMC-1. These problems usually originate in either the analytical column or the post-column detection chemistry. Before checking the TMC-1 as the source of system background noise, consult the appropriate troubleshooting sections in the analytical column Product Manual, the Ion Chromatograph Operator's Manual and the detector manual.

If the source of the high background noise is isolated to the TMC-1, then proceed with the following steps:

- a) Make sure that the eluents and regenerant are correctly formulated.
 - b) Make sure that the eluents are made from chemicals with the recommended purity (see Section 2, "Operation").
 - c) Make sure that deionized water used to prepare the reagents has a specific resistance of 18.2 megohm-cm.
-

3.3 Poor Peak Shape

In some instances, poor peak shape in Chelation IC may be caused by a contaminated TMC-1. To clean the TMC-1, wash with 2.0M HNO₃ for 10 minutes at 3.0 mL/min. Following the HNO₃ wash, rinse the column with deionized water for 3 min at 3.0 mL/min. Replace the cleaned column in the Chelation IC System and run through the Chelation IC program once before doing an analytical run.

Corporate Headquarters

Dionex Corporation
1228 Titan Way
P.O. Box 3603
Sunnyvale, CA 94088-3603
Phone: 408 737-0700
Fax: 408 730-9403

**Salt Lake City
Technical Center**

1515 West 2200 South, Suite A
Salt Lake City, UT 84119-1484
Phone: 801 972-9292
Fax: 801 972-9291

LC Packings Netherlands

LC Packings
Abberdaan 114
1046 AA Amsterdam
The Netherlands
Phone: 31-20-683 9768
Fax: 31-20-685 3452

Dionex Softron GmbH

Dornierstrasse 4
D-82110 Germering
Germany
Phone: 49 8989 468 0
Fax: 49 8989 468 200

US. Regional Offices**Western Region**

Dionex Corporation
445 Lakeside Drive
Sunnyvale, CA 94085
Phone: 408 737-8522
Fax: 408 733-8748

Southwestern Region

Dionex Corporation
340 N. Sam Houston Pkwy. East
Suite 199
Houston, TX 77060
Phone: 281 847-5652
Fax: 281 847-2203

Central Region

Dionex Corporation
1010 Executive Court
Suite 200
Westmont, IL 60559
Phone: 630 789-3660
Fax: 630 789-3702

Eastern Region

Dionex Corporation
14 E. Stow Road
Suite 100
Marlton, NJ 08053
Phone: 856 596-0600
Fax: 856 596-5358

Southeastern Region

Dionex Corporation
1820 Water Place
Suite 250
Atlanta, GA 30339
Phone: 770 432-8100
Fax: 770 432-8300

International Subsidiaries**Australia**

Dionex Pty Ltd
Unit 31, 2 Chaplin Drive
Lane Cove, NSW 2066
Australia
Phone: 61 2 9420 5233
Fax: 61 2 9420 5244

Austria

Dionex Austria GmbH
Laxenburger Strasse 220
A-1230 Wien
Austria
Phone: 431 616 51 25
Fax: 431 616 51 25 55

Belgium

Dionex N. V.
Draaiboornstraat 6
2160 Wommelgem
Belgium
Phone: 32 3 353 4294
Fax: 32 3 353 4293

Canada

Dionex Canada Ltd.
1540 Cornwall Road
Suite 204
Oakville, Ontario L6J 7W5
Canada
Phone: 905 844-9650
Fax: 905 844-6134

China

Dionex China Ltd.
Room 2810, Level 28
Metroplaza Tower 2
No. 223 Hing Fong Road
Kwai Chung, N.T.
Hong Kong
China
Phone: 852 2428 3282
Fax: 852 2428-7898

Denmark

Dionex Denmark A/S
Egegaardsvej 41-st
2610 Rodovre
Denmark
Phone: 45-36-36-90-90
Fax: 45-36-36-90-99

England

Dionex (UK) Ltd.
4 Albany Court
Camberley
Surrey, GU16 7QL
England
Phone: 44 1276 691722
Fax: 44 1276 691837

France

Dionex S.A.
164-166 AVENUE Joseph Kessel
78960 Voisins Le Bretonneux
France
Phone: 33 1 39 30 01 10
Fax: 33 1 39 30 01 12

Germany

Dionex GmbH
Am Woertzgarten 10
D-65510 Idstein
Germany
Phone: 49 6126 991-0
Fax: 49 6126 991277

Italy

Dionex S.r.l.
Via della Maglianella, 65R
00166 Roma
Italy
Phone: 39 06 66 51 5052
Fax: 39 06 66 51 5057

Japan

Nippon Dionex K.K.
DNX Shin-Osaka Bldg.
6-3-14 Nishi-Nakajima
Yodogawa-ku
Osaka 532-0011
Japan
Phone: 81 6 6885-1213
Fax: 81 6 6885-1215

Korea

Dionex Korea Ltd.
In Sung Building 89-106
Shinjung 2-dong
Yangcheon-gu
Seoul, 158-849
Korea
Phone: 82 2 2653-2580
Fax: 82 2 2653-2508

Switzerland

Dionex (Switzerland) AG
Solothurnerstr. 259
4600 Olten
Switzerland
Phone: 41 62 205 99 66
Fax: 41 62 205 99 60

The Netherlands

Dionex B.V.
Lange Bunder 5
4854 MB Bavel
The Netherlands
Phone: 31 161 434303
Fax: 31 161 433065

© 2004 Dionex Corporation
LPN XXXXX 10M 09/04
Printed in U.S.A.



I C | H P L C | M S | E X T R A C T I O N | P R O C E S S | A U T O M A T I O N