

Helium Carrier Gas Conserving Inlet for Gas Chromatography

*Edward B. McCauley, P. Magni, M. Santoro, S. Pelagatti
Thermo Fisher Scientific, Milan, Italy*



Overview

Purpose: We introduce a new, innovative injector, which allows conservation of helium carrier gas. While preserving the analytical GC column flow with helium, it maintains the septum purge and the split flows with another inert gas like nitrogen—even during the analytical run. The analytical performance of the injector is in line with that of the common split/splitless injector module.

Introduction

The helium supply chain crisis has negative implications on research and laboratory operations worldwide. Helium rationing, delayed deliveries and price increases still cause difficulty in production planning and uncertainty in instrument productive uptime.

Although the GC and GC-MS segment consumes less than one percent of global helium usage per year, the shortages and delivery interruptions have wide-spread consequences for many industries utilizing varied analytical techniques.

There are different ways to tackle this problem. One alternative is reducing the split ratio during the split time of the injection. Another is to switch to another carrier gas, such as hydrogen, or to switch gases only when the GC or GC/MS are not running.

All of these solutions have downsides. Reducing the split flow is not suitable for splitless injections and can affect accuracy for smaller volume injections. Hydrogen presents some safety concerns and requires translation and optimization of the methods of use. Switching gases when the instrument is not running is not applicable to routine labs and requires some time to switch back to helium before resuming operations.

The Thermo Scientific™ Instant Connect Helium Saver Injector Module offers a massive reduction in helium consumption while working with the usual split/splitless injector. The Helium Saver module benefits from the modularity concept of all of the injectors and detectors for the Thermo Scientific™ TRACE™ 1300 Series GC. It is easily installed and quickly operative, maximizing instrument uptime. The Helium Saver module can also be mounted on pre-existing TRACE 1300 Series GC systems, without modifications.

FIGURE 1. TRACE 1300 Series GC modularity with user-installable modules.

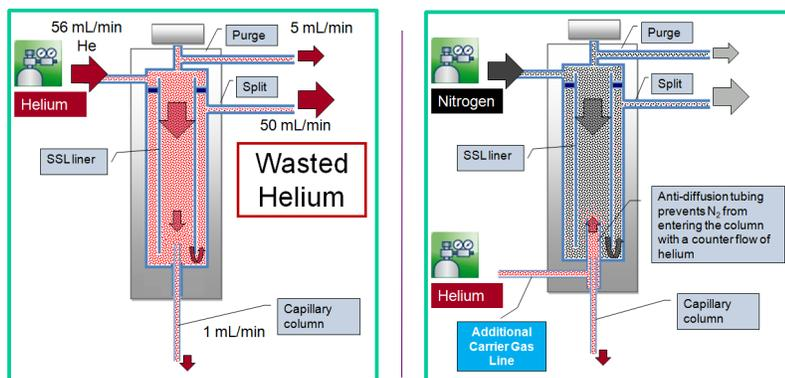


Previously, split/splitless injectors used a single gas as a carrier, for septum purge, and for sample split. Usually, only $\sim 1/10^{\text{th}}$ – $1/50^{\text{th}}$ of the total gas flow entered the column.

Purge and split flows cannot be reduced beyond a certain limit without introducing contamination into the column/detector (due to sample matrix and septa particles accumulated in the liner and lines) or air diffusing from septa seals.

The Helium Saver injector uses two different gases: nitrogen is used for septum purge and sample split flows while helium feeds only the analytical column, drastically reducing helium consumption.

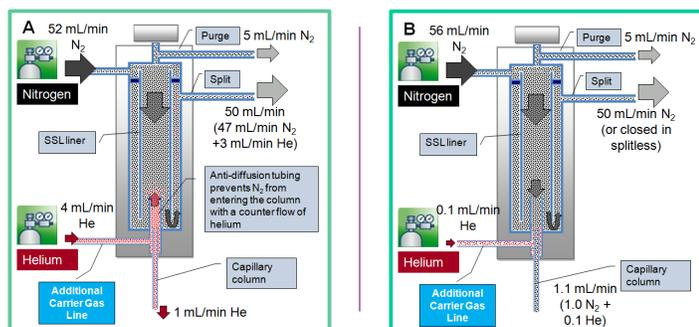
FIGURE 2. Differences between the standard SSL injector and the Helium Saver injector.



How the Helium Saver Injector Works

During all “non-injection” periods, the helium flow is slightly higher than the column flow, drastically reducing helium consumption compared to a standard split/splitless injector (Figure 3A).

FIGURE 3. Helium flow during non-injection (A) and injection (B) phases.



During the injection period, nitrogen flows into the column for sample introduction. Helium is supplied at 0.1 mL/min to keep the connection swept and to avoid dead volumes (Figure 3B).

When the GC is in stand-by, it consumes no helium. Separation is always done using helium carrier gas.

Benefits of the Helium Saver Injector

The Helium Saver injector significantly decrease helium use. A standard helium tank can last up to 3.5 years when used 24/7, 365 days of the year for GC-MS analysis and up to 14.6 years when shutting the helium off or switching to N₂ on weekends and overnight. This means that a single helium cylinder could potentially last for the lifetime of the instrument.

The Instant Connect Helium Saver module is operated exactly like an SSL injector, allowing pain-free translation of any method that involves split/splitless-type injection.

FIGURE 6. Areas and retention time comparison between SSL and Helium Saver Module.

Component	Absolute peak area (counts)			Retention times (min.)				
	Helium only	Helium Saver (1/2 + He)		Helium only	Helium Saver (1/2 + He)		Diff (min)	Diff%
		Counts	Diff%		Counts	Counts		
nC10	1509238	1643138	+8.9	6.878	6.862	-0.017	-0.24	
nC11	1501697	1631649	+8.7	8.312	8.297	-0.015	-0.18	
nC12	1525277	1634623	+7.2	9.733	9.720	-0.013	-0.14	
nC13	1539736	1643560	+6.7	11.100	11.092	-0.008	-0.08	
nC14	1528093	1626721	+6.5	12.393	12.387	-0.007	-0.05	
nC15	1535544	1631809	+6.3	13.618	13.612	-0.007	-0.05	
nC16	1555112	1644967	+5.8	14.778	14.773	-0.005	-0.03	
nC17	1564216	1652088	+5.6	15.877	15.873	-0.003	-0.02	
nC18	1566094	1653903	+5.5	16.922	16.918	-0.003	-0.02	
nC19	1573423	1658194	+5.4	17.917	17.913	-0.003	-0.02	
nC20	1588978	1677003	+5.5	18.867	18.865	-0.002	-0.01	
nC21	1584344	1669784	+5.4	19.773	19.772	-0.002	-0.01	
nC22	1585427	1668266	+5.2	20.643	20.640	-0.003	-0.02	
nC23	1582252	1661415	+5.0	21.477	21.475	-0.002	-0.01	
nC24	1588109	1665599	+4.9	22.278	22.277	-0.002	-0.01	
nC25	1568167	1640702	+4.6	23.047	23.047	0.000	0.00	

No changes on RT (few 1/1000 of a minute) nor on peak areas

Repeatability has been tested in both split and splitless mode, along with the linearity of the split ratio. The results confirm the excellent analytical performance of the module.

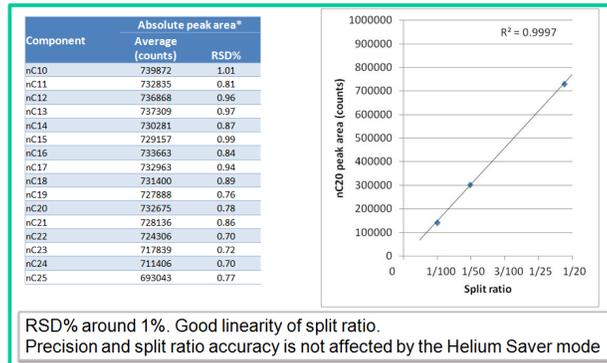
FIGURE 7. Area and retention time repeatability for splitless mode.

Component	Absolute peak area		Retention times		
	Average (counts)	RSD%	Average (min.)	SD (min.)	RSD%
nC10	1675299	0.48	6.863	0.001	0.01
nC11	1660541	0.59	8.299	0.001	0.01
nC12	1674407	0.66	9.722	0.001	0.01
nC13	1679538	0.53	11.092	0.001	0.01
nC14	1662749	0.54	12.389	0.001	0.01
nC15	1663737	0.71	13.615	0.001	0.01
nC16	1672853	0.48	14.775	0.001	0.01
nC17	1679632	0.50	15.875	0.001	0.01
nC18	1679468	0.45	16.920	0.002	0.01
nC19	1685406	0.50	17.916	0.002	0.01
nC20	1702906	0.54	18.866	0.002	0.01
nC21	1697257	0.56	19.774	0.002	0.01
nC22	1695464	0.56	20.644	0.001	0.01
nC23	1689021	0.61	21.478	0.001	0.01
nC24	1696156	0.65	22.279	0.001	0.01
nC25	1671095	0.64	23.049	0.001	0.00

RSD% below 1% on absolute peak areas and SD in the range of 1/1000 of a minute on retention time

Precision is not affected by the Helium Saver Module

FIGURE 8. Area and retention time repeatability for split mode and split ratio linearity.



The injector module has been tested for high-boiling compounds discrimination and activity towards labile compounds, always with excellent results.

FIGURE 9. Discrimination results on a C10-C40 hydrocarbon mix.

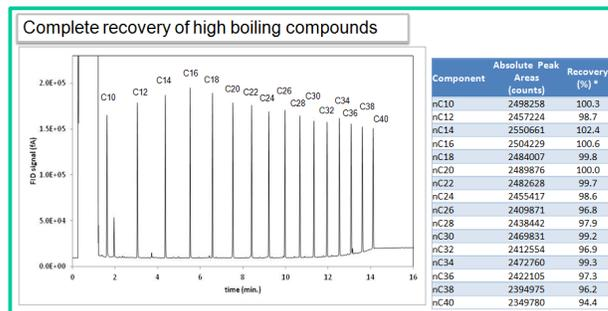
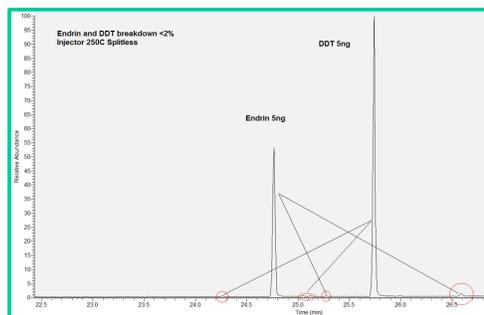


FIGURE 10. Endrin and DDT breakdown.



Conclusion

The Instant Connect Helium Saver Module is an innovative injector whose revolutionary concept benefits users with significant and continuous helium gas savings, without influencing operator routine. It guarantees analytical performance comparable to that of the most widely used split/splitless injectors. The injector is compatible with all TRACE 1300 Series GC systems and does not require the intervention of a service engineer for installation, allowing immediate savings in cost and productivity.

www.thermoscientific.com

©2014 Thermo Fisher Scientific Inc. All rights reserved. ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.



Thermo Fisher Scientific,
Austin, TX USA is
ISO 9001:2008 Certified.

Africa +43 1 333 50 34 0
Australia +61 3 9757 4300
Austria +43 810 282 206
Belgium +32 53 73 42 41
Canada +1 800 530 8447
China 800 810 5118 (free call domestic)
400 650 5118

Denmark +45 70 23 62 60
Europe-Other +43 1 333 50 34 0
Finland +358 9 3291 0200
France +33 1 60 92 48 00
Germany +49 6103 408 1014
India +91 22 6742 9494
Italy +39 02 950 591

Japan +81 45 453 9100
Latin America +1 561 688 8700
Middle East +43 1 333 50 34 0
Netherlands +31 76 579 55 55
New Zealand +64 9 980 6700
Norway +46 8 556 468 00
Russia/CIS +43 1 333 50 34 0

Singapore +65 6289 1190
Spain +34 914 845 965
Sweden +46 8 556 468 00
Switzerland +41 61 716 77 00
UK +44 1442 233555
USA +1 800 532 4752

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand