

Now sold under the  
Thermo Scientific brand

**Thermo**  
SCIENTIFIC

## Separation of Sinapine Thiocyanate in Semen Raphani Using an Acclaim Phenyl-1 Column

Semen Raphani (Figure 1), the seed of *Raphanus sativus L.*, is a Chinese medicinal plant commonly used for treatment of dyspepsia and hypertension. Its anti-hypertension properties are attributed to the presence of sinapine thiocyanate.<sup>1</sup> The Chinese Pharmacopoeia (CP) monitors the quality control of Semen Raphani with a reversed-phase (RP) high-performance liquid chromatography (HPLC) method for the determination of sinapine thiocyanate (structure shown in Figure 2).<sup>2</sup> The method specifies a stationary phase with phenyl groups bonded to silica.

This work describes a separation of sinapine thiocyanate using an Acclaim® Phenyl-1 column, which is based on covalent modification of high-purity, spherical, and porous silica particles with a specially designed silane ligand bearing a proprietary alkyl aromatic functionality.<sup>3</sup> Figure 2 shows the separation of sinapine thiocyanate extracted from Semen Raphani following the CP method. The analyte of interest eluted as a symmetrical peak with adequate retention. The UV spectra of sinapine thiocyanate collected in the standard, Semen Raphani sample, and spiked Semen Raphani sample are highly consistent. The purity of sinapine thiocyanate may be estimated using the peak purity match factor, which can be calculated by Chromeleon® Chromatography Data System (CDS) software.

The calculated peak purity match factor for sinapine thiocyanate separated from the Semen Raphani extract is 950 (the corresponding value for 100% purity is 1000). These results demonstrate that the Acclaim Phenyl-1 column provides good selectivity and is suitable for analysis of sinapine thiocyanate.



Figure 1. Semen Raphani.

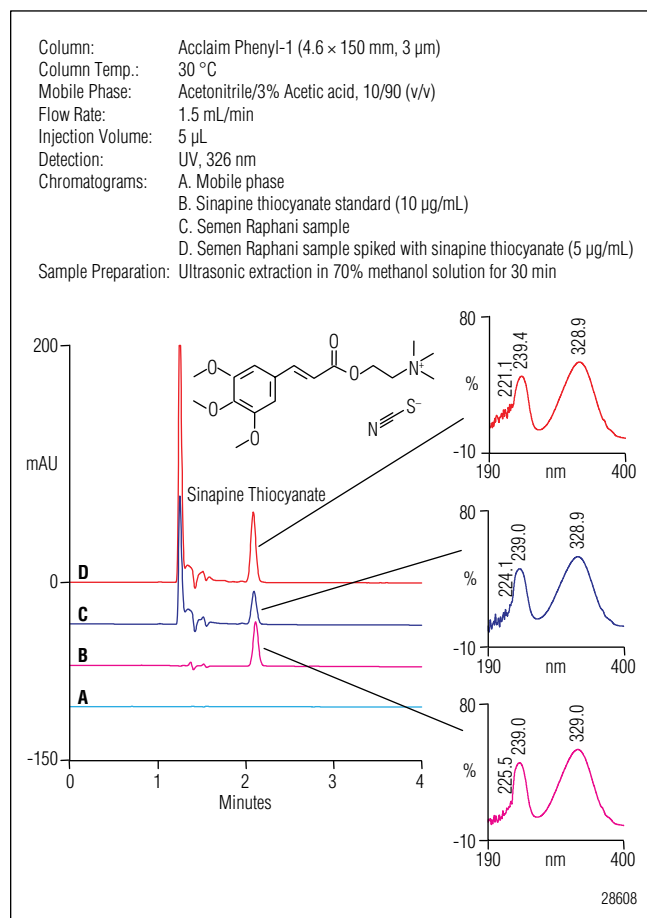


Figure 2. Chromatograms of sinapine thiocyanate on the Acclaim Phenyl-1 column following the CP method.

## EQUIPMENT

Dionex UltiMate® 3000 RSLC system, including:

LPG 3400RS pump

WPS 3000RS autosampler

TCC 3000RS thermostatted column compartment

DAD 3000RS UV-vis detector

Chromeleon 6.80 SR9 software

## REFERENCES

- Li, T.Y.; Li, T.G.; Zhang, G.X.; Guo, H.Y.; Liu, J.; Li, B.G.; Piao Z.Y.; Gai G.G. Experimental Study of Hypotensive Effect of Soluble Semen Raphani Alkaloids in Spontaneous Hypertensive Rats. *World J. Integrated Traditional and Western Med.* **2007**, 2 (1), 25.
- Chinese Pharmacopoeia (Vol. 1), 2010, p 255.
- Dionex Corporation, *Acclaim Phenyl-1 Unique Reversed-Phase Column with High Aromatic Selectivity*, Sunnyvale, CA. <http://www.dionex.com/en-us/products/columns/lc/reversed-phase/acclaim-phenyl/lp-87227.html> (accessed April 4, 2011).

Acclaim, Chromeleon, and UltiMate are registered trademarks of Dionex Corporation.

Speed • Simplicity • Solutions



### Dionex Corporation

1228 Titan Way  
P.O. Box 3603  
Sunnyvale, CA  
94088-3603  
(408) 737-0700

### North America

U.S./Canada (847) 295-7500

### South America

Brazil (55) 11 3731 5140

### Europe

Austria (43) 1 616 51 25 Benelux (31) 20 683 9768 (32) 3 353 4294  
Denmark (45) 36 36 90 90 France (33) 1 39 30 01 10 Germany (49) 6126 991 0  
Ireland (353) 1 644 0064 Italy (39) 02 51 62 1267 Sweden (46) 8 473 3380  
Switzerland (41) 62 205 9966 United Kingdom (44) 1276 691722

### Asia Pacific

Australia (61) 2 9420 5233 China (852) 2428 3282 India (91) 22 2764 2735  
Japan (81) 6 6885 1213 Korea (82) 2 2653 2580 Singapore (65) 6289 1190  
Taiwan (886) 2 8751 6655

[www.dionex.com](http://www.dionex.com)

LPN 2813 PDF 08/16  
©2016 Dionex Corporation