

Separation of Etoposide Using an Acclaim Phenyl-1 Column

INTRODUCTION

Etoposide, a semisynthetic derivative of podophyllotoxin, is used in treating patients with a variety of malignant tumors.¹ Its analysis by reversed-phase high-performance liquid chromatography (HPLC) has been published by the United States Pharmacopeia (USP) and Chinese Pharmacopoeia (CP).^{2,3}

The USP related compounds method uses a 50 min gradient on a column containing packing L11. The USP column packing L11 is defined as phenyl groups chemically bonded to porous silica particles 1.5 to 10 μm in diameter. The particle size requirement in the USP etoposide monograph for the related compounds method is less than 5 μm in diameter. The resolution (R_s) between etoposide and propylparaben required in the related compounds test should be not less than 1.1. The CP method is similar to the USP method but adds a requirement that the retention time (t_R) of etoposide be approximately 25 min.

The work shown here describes a separation of etoposide using an Acclaim[®] Phenyl-1 column. The Phenyl-1 column is based on covalent modification of high-purity, spherical, porous silica particles (3 μm), with a specially designed silane ligand-bearing proprietary alkyl aromatic functionality.⁴

Now sold under the
Thermo Scientific brand

Thermo
SCIENTIFIC

Figure 1A shows chromatograms of etoposide and propylparaben following the USP method. The R_s between etoposide and propylparaben is 8.7, much better than required in the USP method; and the t_R of etoposide is close to 25 min, which meets the requirement in the CP method.

Figure 1B shows a faster and simpler method for the separation of etoposide and propylparaben using the Phenyl-1 column with isocratic elution. The separation of etoposide is completed within 7 min with excellent resolution ($R_s = 5.2$) between etoposide and propylparaben.

EQUIPMENT

Dionex UltiMate® 3000 RSLC system including:

HPG 3400RS pump

WPS 3000RS autosampler

TCC-3000RS thermostatted column compartment

DAD-3000RS UV-vis detector

Chromleon® Chromatography Data System (CDS) software version 6.80 SR9

REFERENCES

1. Kato, Y.; Mawatari, H.; Nishimura, S.I.; Sakura, N.; Ueda, K. Determination of Etoposide Serum Concentrations in Small Pediatric Samples by an Improved Method of Reversed-Phase High-Performance Liquid Chromatography. *Acta Med. Okayama* **2003**, *57* (1), 21–24.
2. The United States Pharmacopeia, USP34–NF29, 2788.
3. Chinese Pharmacopoeia (Vol. 2), 2010, 472.
4. 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 Jan 6, 2011).

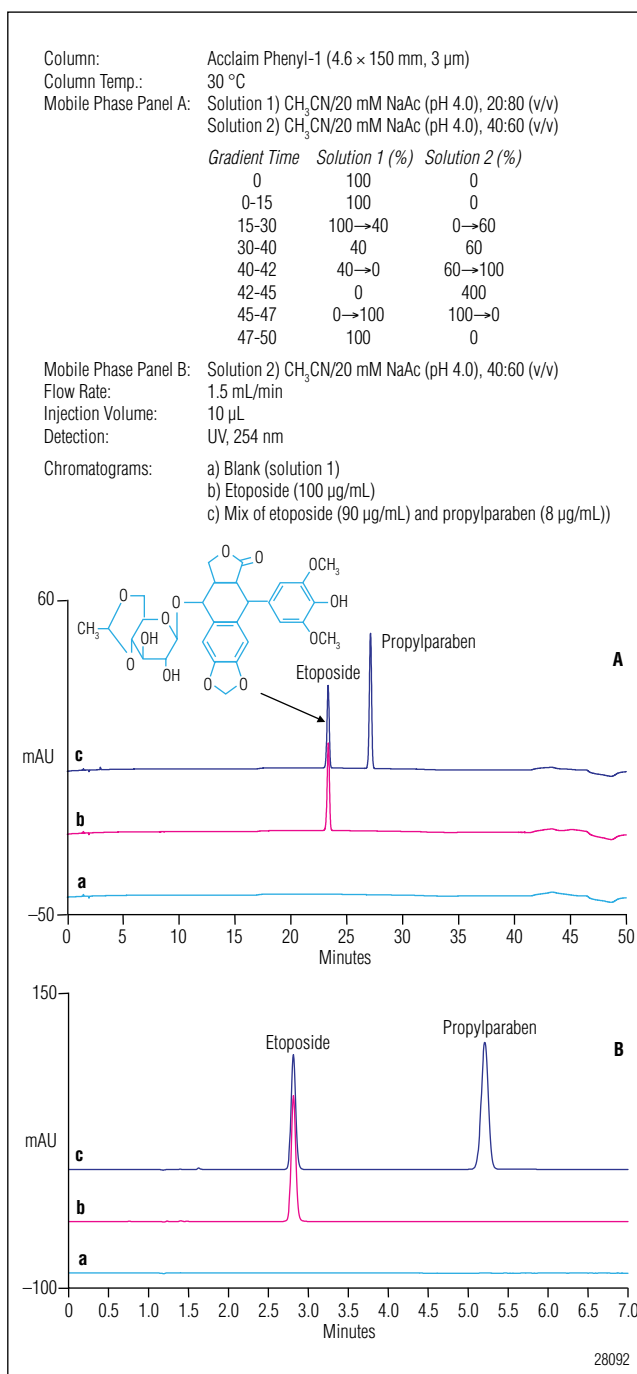


Figure 1. Chromatograms of etoposide following (A) the USP method and (B) isocratic method, using the Acclaim Phenyl-1 column.

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

Passion. Power. Productivity.



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

LPN 2730 PDF 09/16
 ©2016 Dionex Corporation