

# Experimental Protocols for Preparative Purifications Using Macrocyclic Glycopeptide Chiral Stationary Phases

---

## Introduction

---

CHIROBIOTIC phases offer unique opportunities for preparative purifications. Key factors are:

1. No solvent limitations. Halogenated solvents as well as very polar solvents are well tolerated on these CSPs. This solvent tolerance is especially useful when optimizing for sample solubility.
2. Same CSP can be run in four distinctly different mobile phase types. Use of acid/base on any one of these CSPs does not preclude their use in other mobile phases. Mobile phases listed here in the order of success:
  - a. Polar ionic mode<sup>®</sup>:  
MeOH/Acid/Base or  
MeOH/NH<sub>4</sub> salt
  - b. Reversed phase mode:  
ACN/Buffer

*Please note efficient workup procedures available for the reversed phase mode. Method outlined for analytical C18 column. See workup following Example 3: Warfarin*

- c. Polar organic mode;  
MeOH, ACN or combination
- d. Normal phase mode:  
Heptane/EtOH

*\*Note: The chiral recognition mechanism is different in each of these mobile phase types.*

3. Very long term stability with these CSPs due, in part, to the multiple linkages used in anchoring the CSP and, secondly, to the mild conditions typically required.

4. Range of capacities is compound dependent. Significantly overlaps cellulose and amylase phases based on throughput primarily because separations on these CSPs are usually very fast. Capacities on CHIROBIOTIC V2/T2 phases have been 2.5mg/gm with an alpha of 1.5. Maximum capacity achieved was 300mg "on column" using a 250x21.2mm column with an alpha of 2.0.
5. Excellent economics especially with the polar organic and polar ionic modes. Ionic interactions play a significant role in the chiral recognition mechanism on these phases. Solvents here are anhydrous and more volatile and less toxic than the typical normal phase mode.

## Optimization Studies

---

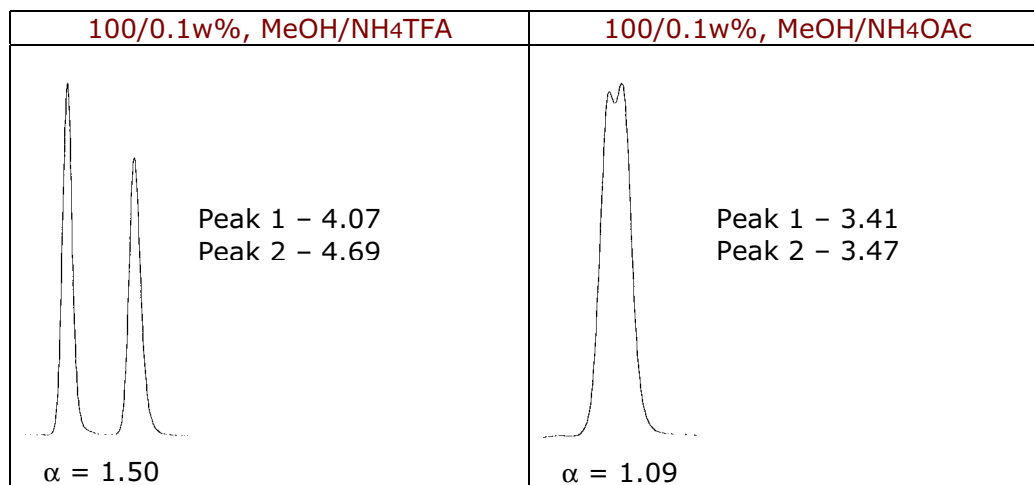
1. Prime consideration in dealing with preparative applications is the balance between selectivity and sample solubility. See Example 3.
2. Check the CHIROBIOTIC V2 and T2 if the polar organic or polar ionic modes have been chosen. These two CSPs may offer increased resolution and increased capacity in these modes only. For neutral molecules, the polar organic mode will work best eliminating the need for acid/base or volatile salt.
3. When operating in the polar ionic mode check carefully the choice of volatile salt as it may effect selectivity dramatically. Ammonium trifluoroacetate usually favors bases while ammonium acetate favors acids.

## CASE STUDY: Nicardipine

Optimization for Prep

Column: CHIROBIOTIC V

FLOW RATE: 1.0 mL/min.



## Issues on Mass Overload

1. Study injection volume and concentration effects:  
Compromise between injection volume and analyte concentration is typically required. For neutral molecules a smaller volume can be used but for separations in the Polar Ionic Mode samples have to be more dilute and larger volumes injected.
2. Salts of chiral drugs:
  - a. Increased buffer strength. When a large amount of sample is injected onto the column consideration has to be given to the quick disassociation of the sample. Usually increasing the buffer strength by a factor of 2 will accomplish this task.
  - b. Solubility. If the salt reduces the solubility of the sample, water may be added for all modes but especially in the polar ionic mode. See example 6.
3. Overload studies:  
Minimize detector overload by using a less sensitive wavelength. It is necessary to evaluate column overload

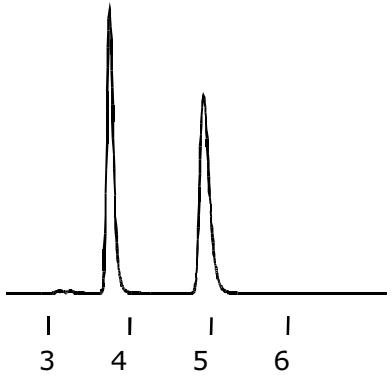
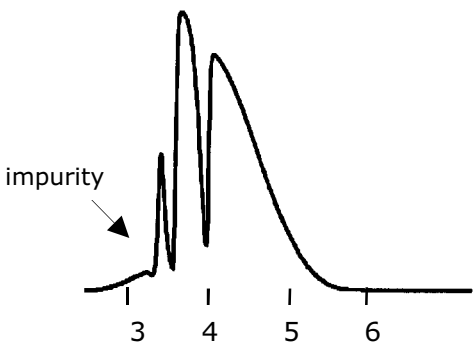
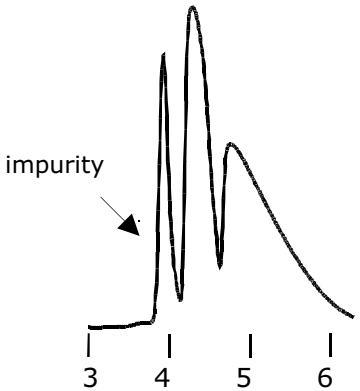
not detector overload. The analytical column is injected with increasing amounts of sample until resolution decreases to about 85%. This number can then be used for direct scale up. From a 250x4.6mm column to a 21.2mm id is a factor of 20x. From the analytical to a 50mm id column is a factor of 120x.

4. Non-linear adsorption isotherms are observed when overloading, i.e., retention decreases with increasing load and peak splitting can result at some critical point.
5. Connecting tubing. To minimize band spread keep tubing connections to 0.01" (0.25mm) or 0.02" (0.5mm) especially if recycle and shave methodologies are used.

The following examples were chosen as simple demonstrations of some of the points covered above. They do not represent the best or the worst examples of what is currently a broad base of experience. Please read CHIROBIOTIC V2/T2 literature for more detailed information about these phases.

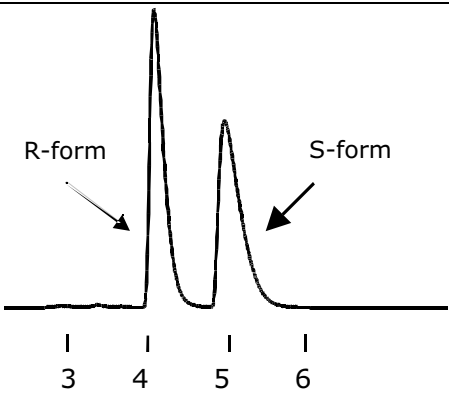
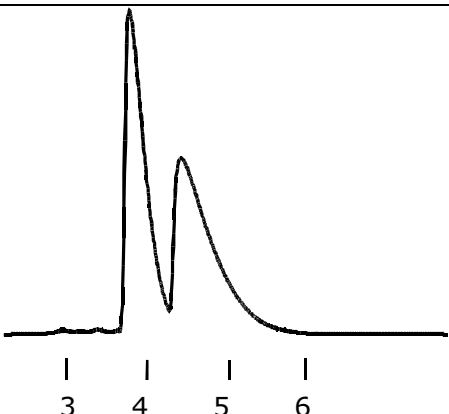
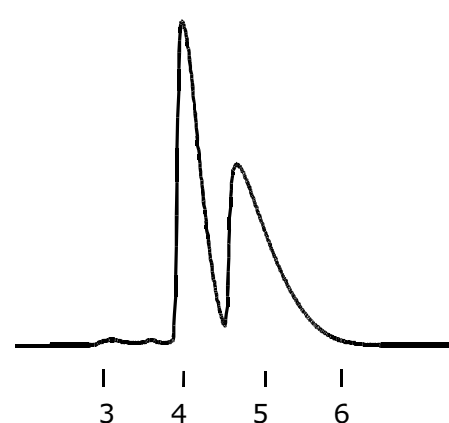
## Example 1: 5-Methyl-5-phenyl hydantoin

Solubility: >80 mg/mL in MeOH  
Column: CHIROBIOTIC T  
Mobile Phase: 100% MeOH  
Selectivity: 2.1

<p>Analytical: 250x4.6mm, 5<math>\mu</math>M Load: 2 mg/mL x 5<math>\mu</math>L (10 <math>\mu</math>g) Flow Rate: 1 mL/min UV: 220nm</p> <p>Peak 1: 3.83 min Peak 2: 5.05 min</p>	 <p>A chromatogram plot with the x-axis labeled from 3 to 6 minutes. Two sharp, well-resolved peaks are visible. The first peak is at approximately 3.83 minutes and the second is at approximately 5.05 minutes. The baseline is flat and stable.</p>
<p>Analytical: 250x4.6mm, 5<math>\mu</math>M Load: 80mg/mL x 0.1mL (8 mg) Flow Rate: 1mL/min UV: 270 nm</p>	 <p>A chromatogram plot with the x-axis labeled from 3 to 6 minutes. A large, broad peak is centered at approximately 5.05 minutes. A much smaller peak is visible at approximately 3.83 minutes, labeled with an arrow and the word "impurity".</p>
<p>Prep: 250x21.2mm, 5<math>\mu</math>M Load: 80mg/mL x 1.5mL (120 mg) Flow Rate: 18 mL/min UV: 270 nm Throughput: 20 mg/g CSP/hr</p>	 <p>A chromatogram plot with the x-axis labeled from 3 to 6 minutes. The profile is very similar to the analytical run at 270 nm, showing a large peak at approximately 5.05 minutes and a small peak at approximately 3.83 minutes labeled "impurity".</p>

## Example 2: 4-Phenyl-2 oxazolidinone

Solubility: >80mg/mL in MeOH  
Column: CHIROBIOTIC T2  
Mobile Phase: 100% MeOH  
Selectivity: 1.6

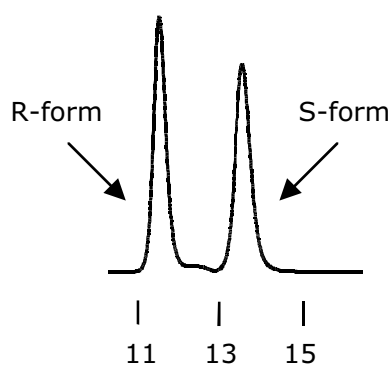
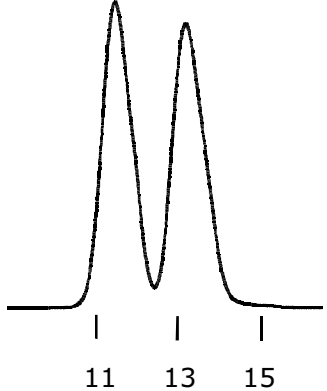
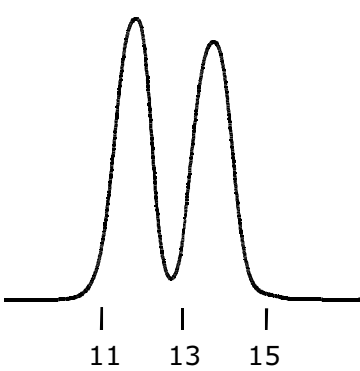
<p>Analytical: 250x4.6mm, 5<math>\mu</math> M Load: 2mg/mL x 5<math>\mu</math>L (10 <math>\mu</math>g) Flow Rate: 1 mL/min UV: 220nm</p> <p>Peak 1: 4.16 min Peak 2: 4.95 min</p>	
<p>Analytical: 250x4.6mm, 5<math>\mu</math>M Load: 80mg/mL x 0.02mL (1.6 mg) Flow Rate: 1 mL/min UV: 254 nm</p>	
<p>Prep: 250x21.2mm, 5<math>\mu</math>M Load: 80mg/mL x 0.5mL (40 mg) Flow Rate: 18 mL/min UV: 254 nm Throughput: 6 mg/g CSP/hr</p>	

## Example 3: Warfarin

Column: CHIROBIOTIC V

### Solubility vs. Selectivity

EtOH/0.1% TEAA, pH 4.1	k <sub>1</sub>	Solubility mg/mL	Selectivity
20/80	4.94	0.1	1.50
30/70	2.19	0.3	1.60
40/60	1.31	0.6	1.47
50/50	0.29	3.0	1.45
80/20	0.1	15.0	1.0

<p>Analytical: 250x 4.6 mm, 5µM  Mobile Phase:  40/60, EtOH/0.1% TEAA, pH 4.1  Selectivity: 1.47  Flow Rate: 0.5 mL/min  Load: 2mg/mL x 5µL (10 µg)  UV: 230 nm</p> <p>Peak 1: 11.33 min  Peak 2: 13.64 min</p>	
<p>Analytical: 250 x 4.6 mm, 5µM  Mobile Phase:  40/60, EtOH/0.1% TEAA, pH 4.1  Flow Rate: 0.5 mL/min  Load: 7.5mg/mL x 0.05mL (0.37 mg)  UV: 330 nm</p>	
<p>Prep: 250 x 21.2mm, 5µM  Mobile Phase:  40/60, EtOH/0.1% TEAA, pH 4.1  Flow Rate: 10 mL/min  Load: 7.5mg/mL x 1 mL (7.5 mg)  UV: 330 nm  Throughput: 0.5 mg/g CSP/hr</p>	

## Recovery From Reversed Phase Systems

The degree of hydrophobicity of the analyte is the main criterion in applying this technique. The following procedure can be used to ascertain the capacity of the C18 column and the size column required for recovery of separated analytes from a full scale prep run.

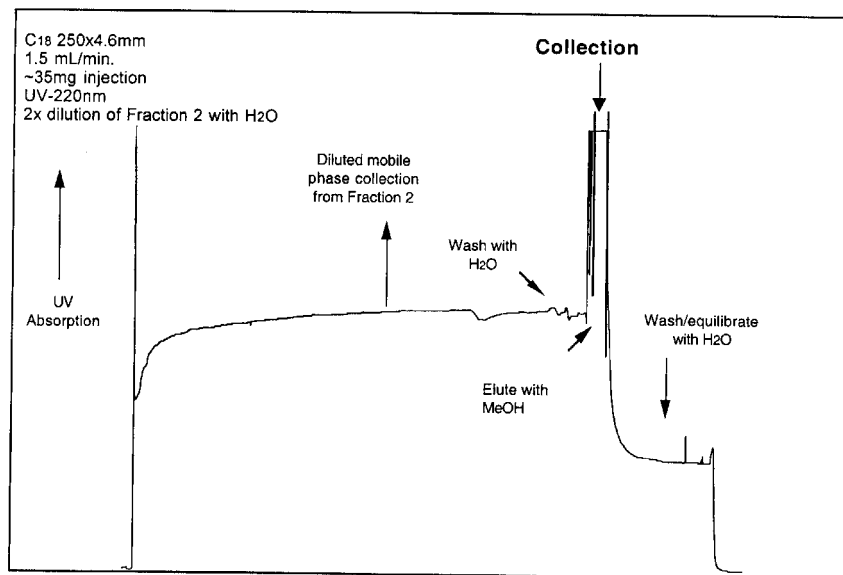
1. Equilibrate an analytical C18 column with HPLC grade water.
2. Pump the recovered chiral column eluant\* containing the enantiomers through the column until the compound breaks through.
3. Wash column with water to remove any buffer present and elute with an appropriate organic solvent. Methanol is often the best choice but ethanol or acetonitrile can be used as well. Measure the volume of collected solvent and assay for recovered analyte.

4. If recovered amount falls below anticipated capacity it is always possible to further dilute the eluant that is being charged with water. In addition, larger size C18 columns can be used.
5. After elution of the compound of interest the column is equilibrated with water for the next addition.
6. From analytical runs it is possible to calculate the size of the column required for a larger scale run.

*\*Note: Adjust pH to increase hydrophobicity and convert all cations and anions to neutral salt. Principle is to suppress the ionization of the analyte.*

Typically doubling the column diameter is equivalent to four times the volume.

### Product Recovery Case Study Warfarin



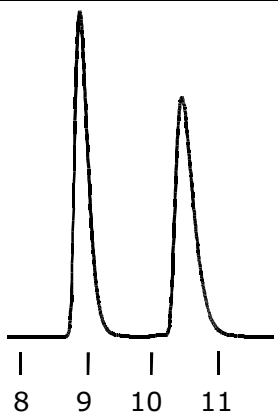
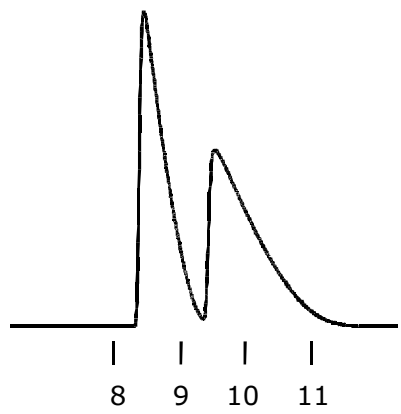
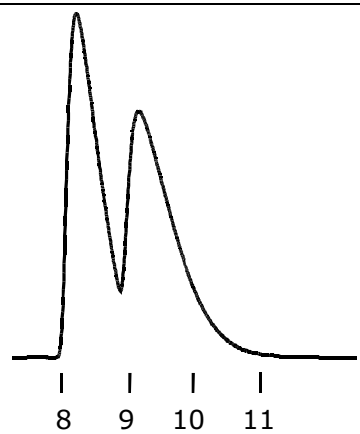
*NOTE: C18 recovery method makes for efficient use of the reversed phase mode with CHIROBIOTIC phases. Also note speed of recovery of the separated enantiomers from a C18 is faster than evaporation of an equivalent volume of heptane.*

Results:

Final Results	Peak 1	Peak 2
ee	98.5%	99%
Yield	90%	85%

## Example 4: Tolperisone

Solubility: >30mg/mL in MeOH  
Column: CHIROBIOTIC V2  
Mobile Phase: 100/0.1w%, MeOH/NH<sub>4</sub>TFA  
Selectivity: 1.33

<p>Analytical: 250x4.6mm, 5<math>\mu</math>M Load: 2mg/mL x 5<math>\mu</math>L (10 <math>\mu</math>g) Flow Rate: 1 mL/min UV: 230nm</p> <p>Peak 1: 8.80 min Peak 2: 10.69 min</p>	 <p>The chromatogram displays two well-resolved peaks. The first peak is at 8.80 minutes and the second is at 10.69 minutes. The x-axis is labeled with minutes 8, 9, 10, and 11.</p>
<p>Analytical: 250x4.6mm, 5<math>\mu</math>M Load: 30mg/mL x 0.03 mL (0.9 mg) Flow Rate: 1mL/min UV: 265 nm</p>	 <p>The chromatogram shows two peaks that are partially overlapping. The first peak is at 8.80 minutes and the second is at 10.69 minutes. The x-axis is labeled with minutes 8, 9, 10, and 11.</p>
<p>Prep: 250x21.2mm, 5<math>\mu</math>M Flow Rate: 20 mL/min Load: 30mg/mL x 0.7mL (21 mg) UV: 265 nm Throughput: 1.8 mg/g CSP/hr</p>	 <p>The chromatogram displays two overlapping peaks. The first peak is at 8.80 minutes and the second is at 10.69 minutes. The x-axis is labeled with minutes 8, 9, 10, and 11.</p>

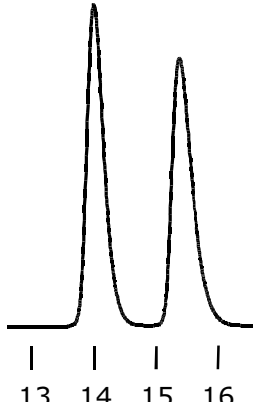
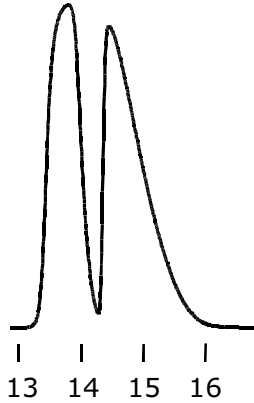
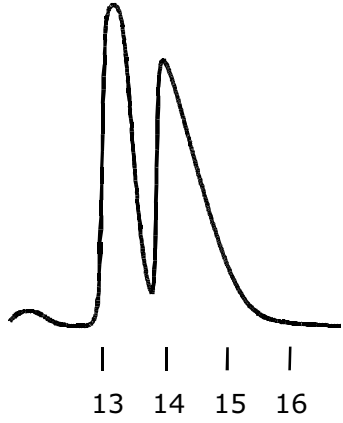
## Example 5: Propranolol

Solubility: >20mg/mL in MeOH

Column: CHIROBIOTIC T

Mobile Phase: 100/0.1w%, MeOH/NH<sub>4</sub>TFA

Selectivity: 1.14

<p>Analytical: 250x4.6mm, 5<math>\mu</math>M Load: 2mg/mL x 5<math>\mu</math>L (10 <math>\mu</math>g) Flow Rate: 1 mL/min UV: 230nm</p> <p>Peak 1: 14.03 min Peak 2: 15.62 min</p>	 <p>A chromatogram plot with the x-axis labeled from 13 to 16 minutes. Two sharp, well-resolved peaks are visible. The first peak is at 14.03 minutes and the second peak is at 15.62 minutes. The baseline is flat and stable.</p>
<p>Analytical: 250x4.6mm, 5<math>\mu</math>M Load: 15mg/mL x 0.7mL (1 mg) Flow Rate: 1mL/min UV: 270 nm</p>	 <p>A chromatogram plot with the x-axis labeled from 13 to 16 minutes. Two peaks are visible, but they are significantly overlapping. The first peak is at 14.03 minutes and the second peak is at 15.62 minutes. The baseline is relatively flat.</p>
<p>Prep: 250x21.2mm, 5<math>\mu</math>M Flow Rate: 22 mL/min Load: 15 mg/mL x 1.6mL (24 mg) UV: 270 nm Throughput: 1.6 mg/g CSP/hr</p>	 <p>A chromatogram plot with the x-axis labeled from 13 to 16 minutes. Two peaks are visible, but they are significantly overlapping. The first peak is at 14.03 minutes and the second peak is at 15.62 minutes. The baseline shows some noise and a slight upward trend towards the end of the run.</p>



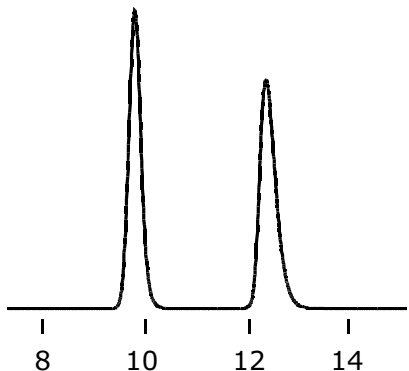
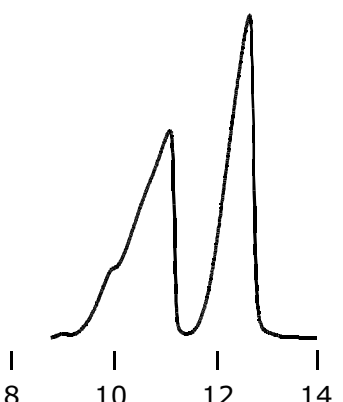
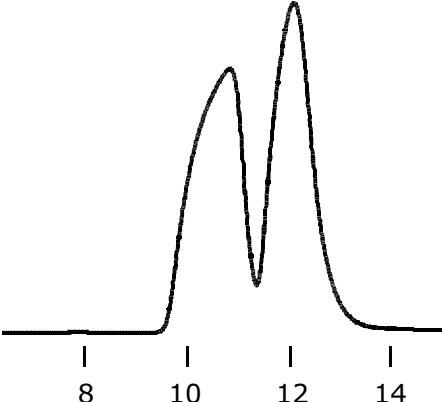
## Example 6: Terbutaline (hemisulfate salt)

Column: CHIROBIOTIC T

### Solubility vs. Selectivity

100/0.1w%, MeOH/NH <sub>4</sub> TFA	k <sub>1</sub>	Solubility mg/mL	Selectivity
No H <sub>2</sub> O*	2.50	2.2	1.37
Add 5% H <sub>2</sub> O (v/v)	2.76	12	1.29
Add 10% H <sub>2</sub> O (v/v)	3.50	21	1.24

*\*Adding water to increase sample solubility can sometimes be effective.  
This phenomena is sample dependent.*

<p>Analytical: 250x 4.6 mm, 5µM Mobile Phase: 100/0.1w%, MeOH/NH<sub>4</sub>TFA Selectivity: 1.37 Flow Rate: 1 mL/min Load: 2mg/mL x 5µL (10 µg) UV: 230 nm</p> <p>Peak 1: 9.81 min Peak 2: 12.26 min</p>	
<p>Analytical: 250 x 4.6 mm, 5µM Mobile Phase: 100/0.1w%, MeOH/NH<sub>4</sub>TFA + 5% H<sub>2</sub>O Selectivity: 1.29 Flow Rate: 1.0 mL/min Load: 8 mg/mL x 0.1mL (0.8 mg) UV: 254 nm</p>	
<p>Perp: 250 x 21.2mm, 5µM Mobile Phase: 100/0.1w%, MeOH/NH<sub>4</sub>TFA + 5% H<sub>2</sub>O Flow Rate: 20 mL/min Load: 7 mg/mL x 3mL (21 mg*) UV: 254 nm Throughput: 1.4 mg/g CSP/hr <i>*Note: higher loading causes peak one to split but does not affect recovery.</i></p>	

Advanced Separation Technologies Inc. presents this information as an aid to the successful use of these products. Exceptions are possible. All statements herein are expressions of opinion which we believe to be accurate and reliable, but are presented without guarantee or responsibility on our part. While Advanced Separation Technologies Inc. has used its best efforts to present useful instructions, no warranty expressed or implied is given with respect to this material or the products recommended herein.

©2003 Advanced Separation Technologies Inc.  
All rights reserved.  
Printed in USA.



*Advanced Separation Technologies Inc.*

World Headquarters: 37 Leslie Court, P. O. Box 297, Whippany, NJ 07981 USA  
Tel: (973) 428-9080 Fax: (973) 428-0152 E-mail [astecusa@aol.com](mailto:astecusa@aol.com)

UK and Ireland Sales Office: 1 Blake Street, Congleton, Cheshire CW12 4DA UK  
Tel: +44 (0) 1260 276276 Fax: +44 (0) 1260 290067 E-mail: [info@asteceuro.com](mailto:info@asteceuro.com)

Visit our website at [www.astecusa.com](http://www.astecusa.com)