



## Bio-Scale CHT10-I Hydroxyapatite Column

(12 mm x 88 mm)    Catalog Number 751-0025

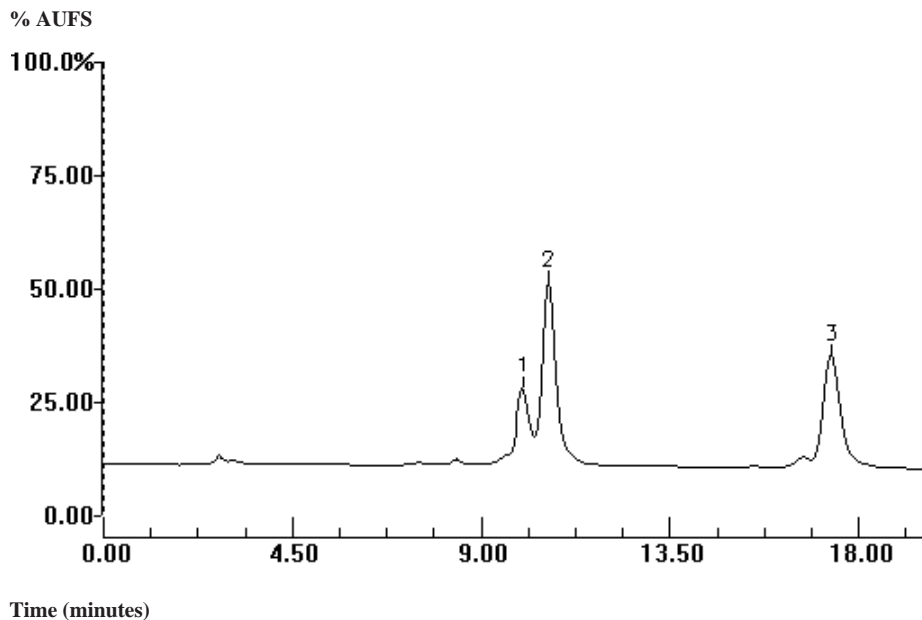
Serial Number: \_\_\_\_\_  
Batch Number: \_\_\_\_\_  
Resin Batch #: \_\_\_\_\_  
Date Tested: \_\_\_\_\_  
  
Checked By: \_\_\_\_\_

### Column Specifications

Efficiency (N):	$\geq 22,700$ / meter
Asymmetry (b/a):	0.85 - 2.50
Column Pressure (2.93 ml/min):	< 160 psi

### Typical Chromatogram\*

<b>Buffer A:</b>	10 mM Sodium Phosphate, pH 6.8	<b>Sample:</b>	Protein Standard for Cation Exchange Chromatography (Catalog Number 125-0562)
<b>Buffer B:</b>	500 mM Sodium Phosphate, pH 6.8	<b>Inj. Vol:</b>	100 $\mu$ l (0.6 mg Total Protein)
<b>Gradient:</b>	0-80% B in 10 minutes	<b>Peaks:</b>	1. Myoglobin 2. Ribonuclease A 3. Cytochrome C
<b>Flow Rate:</b>	2.93 ml/min		
<b>Detection:</b>	UV @ 280 nm, 0.20 AUFS		



\* The above chromatogram and test conditions are provided to help resolve any technical problems.