

Bio-Scale CHT20-I Hydroxyapatite Column

(15 mm x 118 mm) Catalog Number 751-0027

Serial Number: _____
Batch Number: _____
Resin Batch #: _____
Date Tested: _____

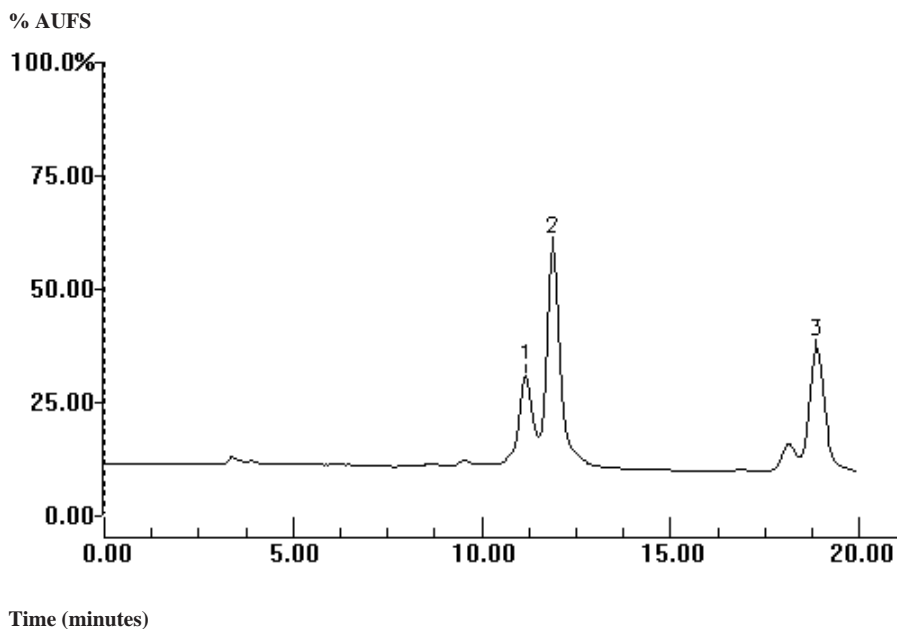
Checked By: _____

Column Specifications

Efficiency (N):	$\geq 17,000$ / meter
Asymmetry (b/a):	0.85 - 2.50
Column Pressure (4.59 ml/min):	< 160 psi

*Typical Chromatogram**

Buffer A:	10 mM Sodium Phosphate, pH 6.8	Sample:	Protein Standard for Cation Exchange Chromatography (Catalog Number 125-0562)
Buffer B:	500 mM Sodium Phosphate, pH 6.8	Inj. Vol:	100 μ l (0.6 mg Total Protein)
Gradient:	0-80% B in 20 minutes	Peaks:	1. Myoglobin 2. Ribonuclease A 3. Cytochrome C
Flow Rate:	4.59 ml/min		
Detection:	UV @ 280 nm, 0.10 AUFS		



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