

Bio-Scale CHT5-I Hydroxyapatite Column

(10 mm x 64 mm) Catalog Number 751-0023

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

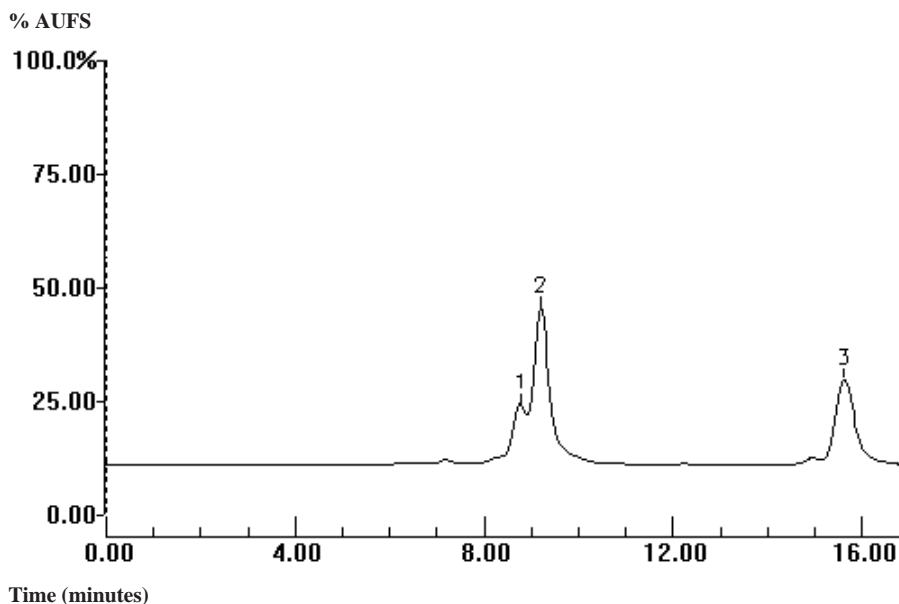
Checked By: _____

Column Specifications

Efficiency (N):	$\geq 20,300$ / meter
Asymmetry (b/a):	0.85 - 2.50
Column Pressure (2.04 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 10 minutes	Peaks:	1. Myoglobin 2. Ribonuclease A 3. Cytochrome C
Flow Rate:	2.04 ml/min		
Detection:	UV @ 280 nm, 0.30 AUFS		



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