

Bio-Scale DEAE5 Anion Exchange Column

(10 mm x 64 mm) Catalog Number 751-0033

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

Checked By: _____

Column Specifications

Efficiency (N):	$\geq 21,000$ / meter
Asymmetry (b/a):	0.85 - 1.50
Column Pressure (2.04 ml/min)	< 160 psi

*Typical Chromatogram**

Buffer A: 20 mM Tris, pH 7.6

Buffer B: 20 mM Tris + 1.0 M Sodium Chloride
pH 7.6

Gradient: 0% B in 6 minutes
0 - 50% B in 20 minutes

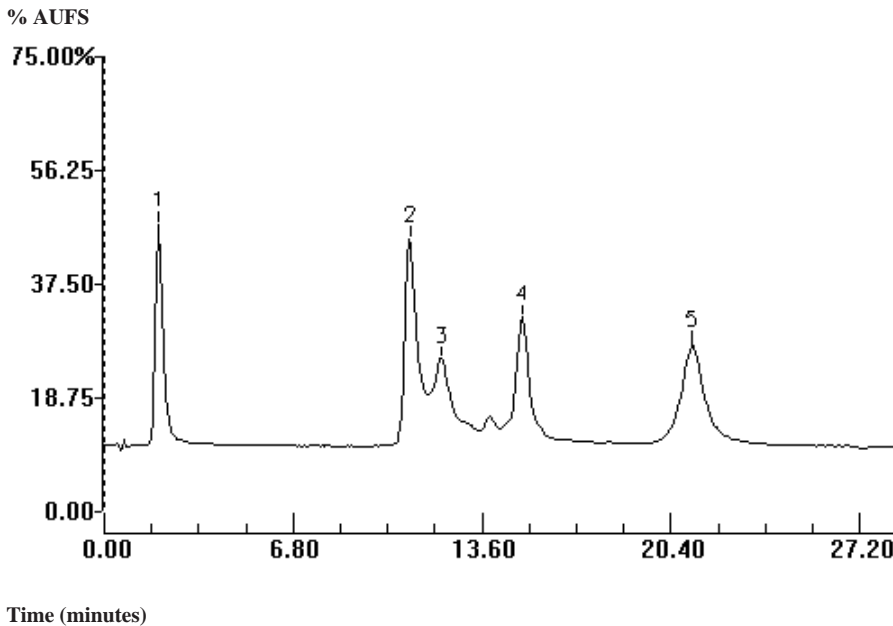
Flow Rate: 2.04 ml/min

Detection: UV @ 280 nm, 0.16 AUFS

Sample: Protein Standard for Anion
Exchange Chromatography
(Catalog Number 125-0561)

Inj. Vol: 250 μ l (2250 μ g Total Protein)

Peaks: 1. Myoglobin
2. Conalbumin - A
3. Conalbumin - B
4. Ovalbumin
5. Soybean Trypsin Inhibitor



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