

## Bio-Scale DEAE5 Anion Exchange Column

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

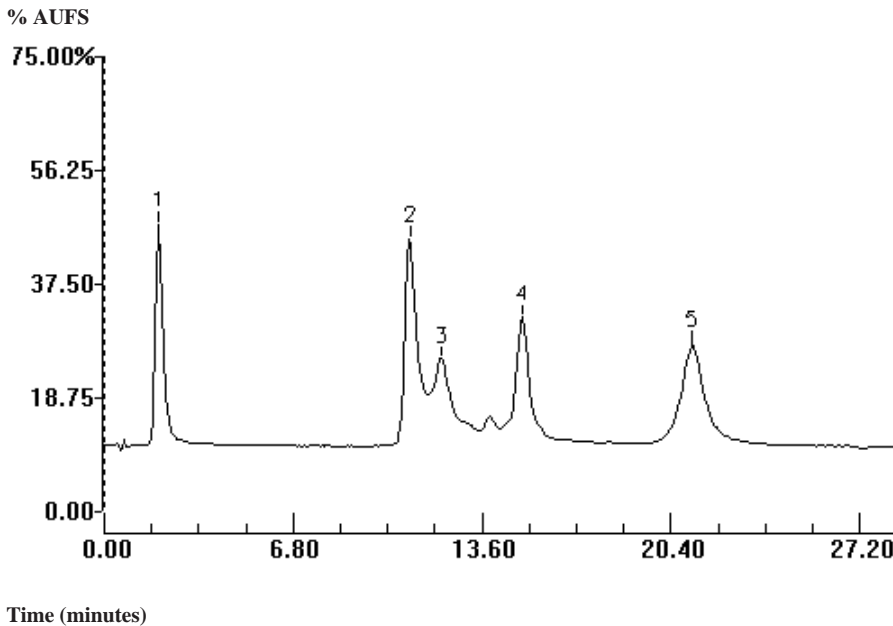
Serial Number: \_\_\_\_\_  
Batch Number: \_\_\_\_\_  
Resin Batch #: \_\_\_\_\_  
Date Tested: \_\_\_\_\_  
Checked By: \_\_\_\_\_

### 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\*

<b>Buffer A:</b>	20 mM Tris, pH 7.6	<b>Sample:</b>	Protein Standard for Anion Exchange Chromatography (Catalog Number 125-0561)
<b>Buffer B:</b>	20 mM Tris + 1.0 M Sodium Chloride pH 7.6	<b>Inj. Vol:</b>	250 $\mu$ l (2250 $\mu$ g Total Protein)
<b>Gradient:</b>	0% B in 6 minutes 0 - 50% B in 20 minutes	<b>Peaks:</b>	1. Myoglobin 2. Conalbumin - A 3. Conalbumin - B 4. Ovalbumin 5. Soybean Trypsin Inhibitor
<b>Flow Rate:</b>	2.04 ml/min		
<b>Detection:</b>	UV @ 280 nm, 0.16 AUFS		



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