

USING THE UNO S1 COLUMN IN THIS SEPARATION SCHEME ALLOWED:

- PURIFICATION OF CRUDE INSULIN UP TO THREE TIMES FASTER THAN A MONODISPERSED BEADED COLUMN
- FAST SEPARATION WITHOUT SACRIFICING RESOLUTION OR PURITY DUE TO THE UNO COLUMN'S UNIQUE CONTINUOUS BED

Purification of Insulin on UNO™ S1 Column

Insulin is a peptide hormone which stimulates anabolic reactions for carbohydrates, proteins, and fats resulting in lowered blood glucose. It is used therapeutically for treatment of diabetes mellitus and was the first recombinant protein on the market. Although insulin is well characterized and understood, it is often difficult to purify because of the harsh pH and denaturing conditions needed to solubilize and disaggregate the protein. A comparison of the UNO S1 (Bio-Rad catalog # 720-0021) and a popular monodispersed cation exchange column showed that it was advantageous to use the UNO S1 column for small scale insulin separation or analysis.

Experimental

The insulin used in this study was an acid precipitate in suspension. The insulin was solubilized in acidic buffer with urea. The insulin sample was applied separately to an UNO S1 and a monodispersed cation exchange column and eluted by linear gradients. Fractions were collected and neutralized with ammonium hydroxide to pH 7.0. Due to the viscosity of the buffers, both columns were run at half the maximum recommended flow rates. The separation performed on the UNO S1 was run at 3 ml/min (Figure 1). The separation performed on the monodispersed cation exchanger was performed at 1 ml/min (Figure 2). Fractions 43–50 and fractions 30–34 were pooled from the separation performed on the UNO S1 and the monodispersed column respectively and analyzed by reversed phase HPLC. Analysis performed on Hi-Pore® RP-304 reversed phase column (Bio-Rad catalog # 125-0551). The products from both columns were 98% pure. Figure 3 shows the HPLC chromatogram of the pooled fractions from the UNO S1 column after subtraction of the background UV from the buffer.

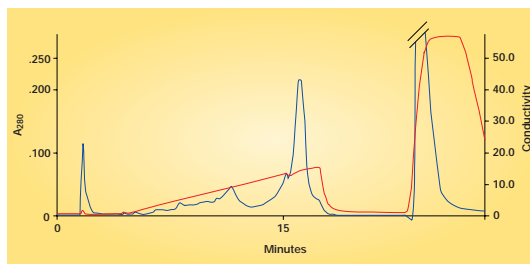


Fig. 1. Chromatogram of insulin purification on UNO S1 column. Buffer A: 0.25 M HOAc, 7.0 M urea. Buffer B: A + 1.0 M NaCl. Gradient: 0–30% B over 40 column volumes. Flow rate: 3 ml/min. Sample size: 0.25 ml.

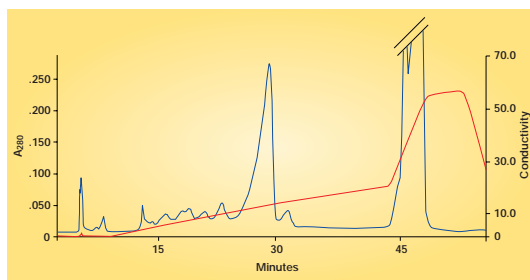


Fig. 2. Chromatogram of insulin purification on monodispersed cation exchange column. Conditions were as in Figure 1 except flow rate was 1 ml/min.

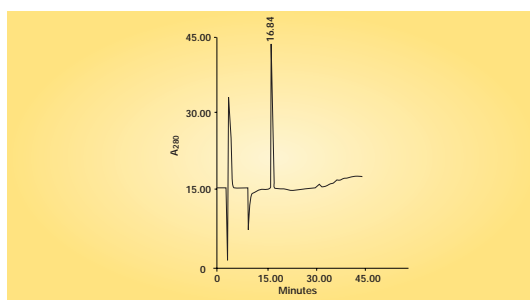


Fig. 3. Chromatogram of HPLC analysis on fractions 43–50 from UNO S1. Buffer A: 0.01% TFA. Buffer B: 95% Acetonitrile 5% buffer A. Flow rate: 1 ml/min.

BIO-RAD

**Bio-Rad
Laboratories**

Life Science
Group

Website www.bio-rad.com U.S. (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771
China 86-10-62051850 Denmark 45 39 17 99 47 Finland 358 (0)9 804 2200 France (01) 43 90 46 90 Germany 089 318 84-0
Hong Kong 852-2789-3300 India (91-11) 461-0103 Israel 03 951 4127 Italy 02-21609.1 Japan 03-5811-6270 Korea 82-2-3473-4460
The Netherlands 31 318-540666 New Zealand 64-9-4152280 Singapore 65-2729877 Spain (91) 661 70 85 Sweden 46 (0)8 627 50 00
Switzerland 01-809 55 55 United Kingdom 0800-181134