

THE UNO Q1 WAS CHOSEN FOR THIS SEPARATION BECAUSE OF ITS UNIQUE CONTINUOUS BED MATRIX WHICH ALLOWS:

- QUICK SCREENING FOR OPTIMUM ELUTION SCHEMES, EVEN WITH VISCOUS BUFFERS
- MINIMAL LOSS OF RESOLUTION AT FAST FLOW RATES, SINCE MASS TRANSFER EFFECTS ARE MINIMAL
- FAST ONE STEP PURIFICATION OF rIL-13 AFTER SAMPLE PREPARATION OF CELL PELLET

## Purification of rIL-13 Using an UNO™ Q1 Column

Interleukin 13 is a recently described cytokine produced by activated T cells. This protein is responsible for both positive and negative responses to regulatory molecules within the immune system. rIL-13 has been shown to modulate macrophage functions as well as play a direct role in regulation of proliferation and differentiation of hematopoietic stem cells. In addition, a recent study has indicated that rIL-13 inhibits human immunodeficiency virus type I production in blood derived macrophages. For this application, recombinant rIL-13 was overexpressed in *E. coli*. This report describes the isolation and purification of rIL-13 using an UNO Q1 (Bio-Rad catalog # 720-0001) column.

### Experimental

rIL-13 was obtained from the cell pellet in a two part process, sample preparation and chromatographic separation. See Figure 1 for a flow chart of sample preparation. Sample preparation involved suspending the cell pellet in 50 mM Tris, pH 8.0, 20 mM DTT with 350 µl of bacterial protease inhibitor (Sigma P8465). The suspension was then passed through a Microfluidics cell disruptor at 12,000 psi five times. The lysed cell suspension was then centrifuged, washed, and the rIL-13 was solubilized with 8 M urea. It was found that rIL-13 will only solubilize with a high concentration of chaotrope under reducing conditions. The final solubilized preparation was then filtered through a 0.45 µm filter and applied to an UNO Q1 column.

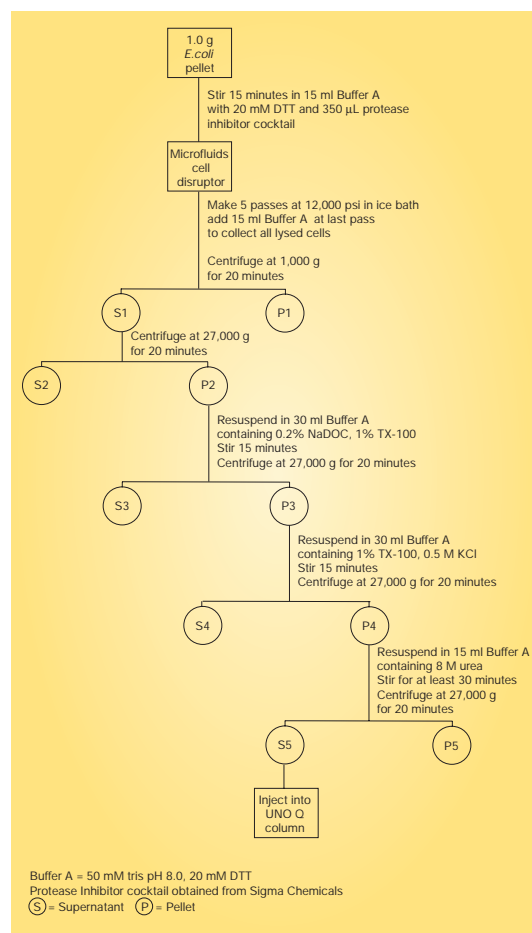


Fig. 1. Flow chart of sample preparation for rIL-13.

The rIL-13 was eluted by a linear gradient as described in Figure 2. Due to the viscosity of the buffers used in the experiment, the UNO Q1 column was run at 3 ml/min.

## Results

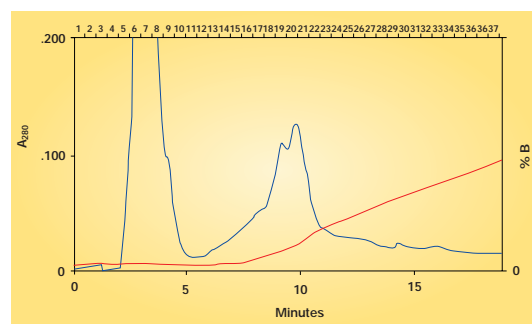
Figure 2 shows the elution profile from the UNO Q1 column. rIL-13 was found to elute from this column in the conductivity range of 1.75-3.55 ms/cm.

SDS-PAGE (Figure 3) analysis using a 15% Tris-HCl Ready Gel (Bio-Rad catalog #161-1157) indicated fractions 18–20 contained rIL-13 of at least 90% purity, while fraction 21 contained a small amount of impurities. A western blot of the gel was performed to identify the rIL-13 (data not shown). The apparent heterogeneity in the elution profile may arise from protein-protein interaction or folding intermediates, since by SDS-PAGE, the rIL-13 was nearly homogeneous.

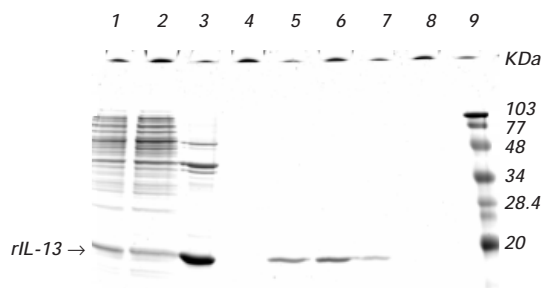
## References

- 1 Doherty, T. M., Kastelein R., Menon S., Andrade S., and Coffman, R. L., *J. Immunol.*, **151**, 7151, (1993).
- 2 Montaner, L. J., Doyle, A. G., Collin, M., Herbein G., Illei, P., James W., Minty, A., Caput D., Ferrara P., and Gordon, S., *J. Exp. Med.*, **178**, 743, (1993).
- 3 Jacobsen, S. E., Okkenhaug, C., Veiby, O. P., Caput D., Ferrara P., and Minty A., *J. Exp. Med.*, **180**, 75, (1994).
- 4 Lai, H. Y., Heslan, J. M., Poppema, S., Elliot, J. F., and Mosmann, T. R., *The Journal of Immunology*, **156**, 3166, (1996).

We gratefully acknowledge receipt of rIL-13 containing E. coli cell pellet constructed by Madeleine Huey working under Dr. John Roberts at UCSF liver transplant.



**Fig. 2.** Chromatogram of rIL-13 separated from fraction S5 on UNO Q1 column. Buffer A: 20 mM CHES, 8.0 M urea, pH 9.8. Buffer B: Buffer A: + 1.0 M NaCl. gradient was 0–20% B over 10 column volumes.



**Fig. 3.** SDS-PAGE of rIL-13 fractions from UNO Q1 column on 15% Tris-Glycine Ready Gel. Lane 1: cell pellet. Lane 2: lysed resuspension. Lane 3: S5 from Figure 1. Lane 4–7: fractions 18–21. Lane 8: fraction 41. Lane 9: MW standard (catalog #161-0305).

**BIO-RAD**

**Bio-Rad  
Laboratories**

Life Science  
Group

Website [www.bio-rad.com](http://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