

# UNO

## Use of an UNO Q1 Column Step in the Purification of Phosphoinositide 3-Kinase from Rat Liver

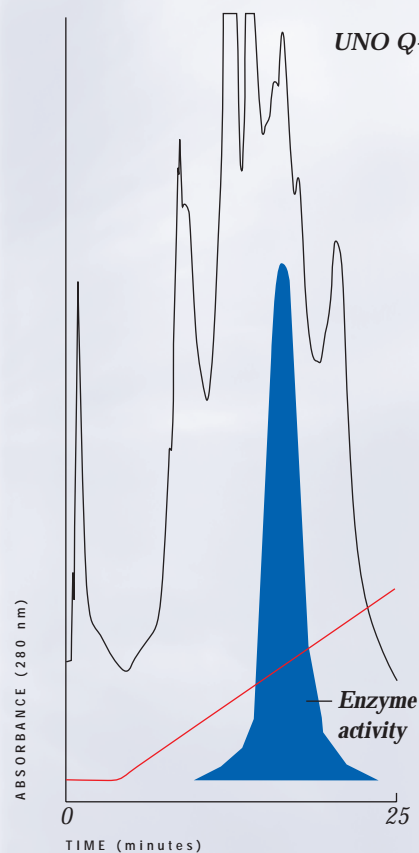
**The homogeneous, nonporous, Continuous Bed matrix allows:**

- Extremely fast mass transfer of proteins
- Minimal band-broadening during elution
- Unsurpassed resolution, even at high flow rates
- Separation times five-fold faster than conventional beaded columns

Polyphosphoinositides and their metabolites are important intracellular signals involved in the responses to a number of hormones and growth factors. The discovery of a phosphatidylinositol kinase that phosphorylates PI at the 3 position of the inositol ring uncovered a new pathway of PI metabolism and potential intracellular signals.

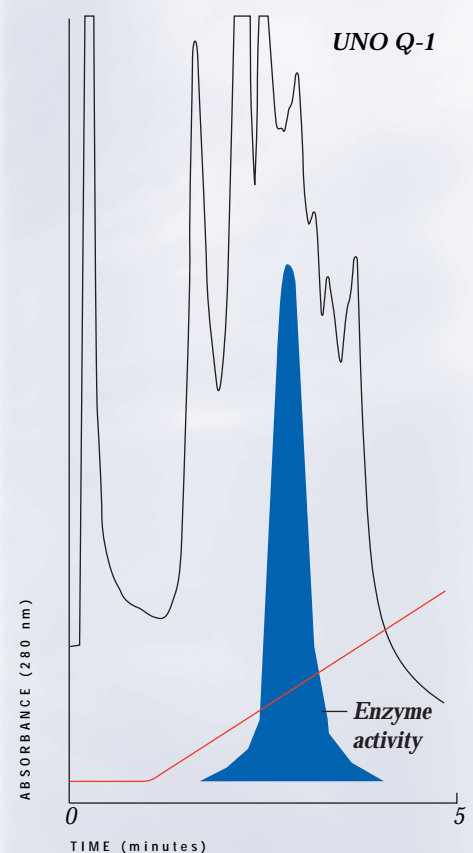
Phosphoinositide 3-kinase has been purified to homogeneity from rat liver and bovine thymus. The enzyme from rat liver is a heterodimer of 110 and 85 kDa proteins. Two forms of the 110 kDa protein exist, and have similar but not identical amino acid sequences. Recent studies have suggested that the 110 kDa protein is the catalytic protein and that the 85 kDa protein plays a regulatory role.

**Figure 1. Purification of Phosphoinositide-3-Kinase at 0.5 ml/min.**



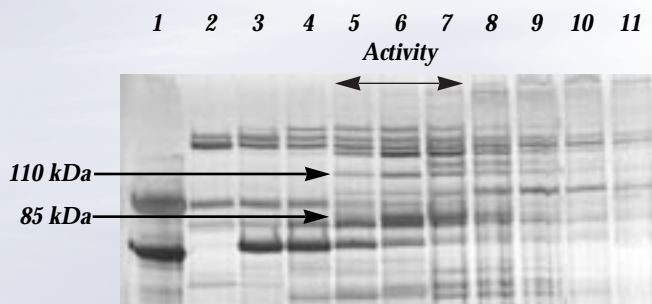
Buffer A: 40 mM TEA, pH 7.6  
Buffer B: A + 1 M KCl, pH 7.6  
Gradient: 0–35% B in 20 ml  
Flow rate: 0.5 ml/min  
System: FPLC®

**Figure 2. Purification of Phosphoinositide-3-Kinase at 2.5 ml/min.**



Buffer A: 40 mM TEA, pH 7.6  
Buffer B: A + 1 M KCl, pH 7.6  
Gradient: 0–35% B in 20 ml  
Flow rate: 2.5 ml/min  
System: FPLC®

**Figure 3. SDS-PAGE of PI-3K Active Fractions.**



Phosphoinositide 3-kinase has proven difficult to purify in an active form when overexpressed in various expression systems. Consequently, for continued biochemical studies of this phosphoinositide pathway, a routine purification protocol for phosphoinositide 3-kinase from rat liver has been established. The protocol involves multiple steps including size exclusion, hydroxyapatite, and ion-exchange chromatography.

The chromatograms (Figures 1 and 2) illustrate the use of an UNO Q1 column to partially purify a cytosolic phosphoinositide 3-kinase from a rat liver homogenate at a flow rate 5-fold higher than previously used with a conventional monodispersed beaded media.

Both the overall resolution and protein activity profile are maintained at the higher flow.

## Acknowledgement

Data courtesy of Dr. A. Couvillon, Beth Israel Deaconess Medical Center, Boston, MA.

FPLC is a trademark of Pharmacia

**BIO-RAD**

**Bio-Rad  
Laboratories**

Life Science  
Group

**Website** [www.bio-rad.com](http://www.bio-rad.com) **Bio-Rad Laboratories Main Office** 2000 Alfred Nobel Drive, Hercules, California 94547, Ph. (510) 741-1000, Fx. (510)741-5800  
**Also in:** **Australia** Ph. 02-9914-2800, Fx. 02-9914-2889 **Austria** Ph. (1)-877 89 01, Fx. (1) 876 56 29 **Belgium** Ph. 09-385 55 11, Fx. 09-385 65 54  
**Canada** Ph. (905) 712-2771, Fx. (905) 712-2990 **China** Ph. (86-10) 2046622, Fx. (86-10) 2051876 **Denmark** Ph. 39 17 9947, Fx. 39 27 1698  
**Finland** Ph. 90 804 2200, Fx. 90 804 1100 **France** Ph. (1) 43 90 46 90, Fx. (1) 46 71 24 67 **Germany** Ph. 089 31884-0, Fx. 089 31884-100  
**Hong Kong** Ph. 7893300, Fx. 7891257 **India** Ph. 91-11-461-0103, Fx. 91-11-461-0765 **Israel** Ph. 03 951 4127, Fx. 03 951 4129  
**Italy** Ph. 02-21609.1, Fx. 02-21609.399 **Japan** Ph. 03-5811-6270, Fx. 03-5811-6272 **The Netherlands** Ph. 0313 18-540666, Fx. 0313 18-542216  
**New Zealand** Ph. 09-443 3099, Fx. 09-443 3097 **Singapore** Ph. (65) 272-9877, Fx. (65) 273-4838 **Spain** Ph. (91) 661 70 85, Fx. (91) 661 96 98  
**Sweden** Ph. 46 (0) 8 627 50 00, Fx. 46 (0) 8 627 54 00 **Switzerland** Ph. 01-809 55 55, Fx. 01-809 55 00 **United Kingdom** Ph. 0800 181134, Fx. 01442 259118