

## Purification of elongation factors, EF-1 $\beta\gamma\delta$ and EF-1 $\beta\gamma$ , from *Xenopus Laevis* Oocytes Using UNO Q1 Column

**Incorporating UNO into this purification scheme resulted in:**

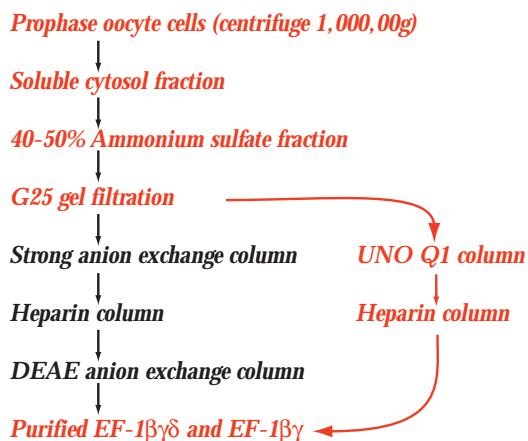
- **Reduction** of the number of purification steps
- **High degree of resolution** achieved at a flow-rate **3x faster** than a conventional beaded column

### Introduction

Elongation factor 1 (EF-1) mediates the elongation step of mRNA translation. The transfer of aminoacyl-tRNA to ribosomes mediated by hydrolysis of GTP is catalyzed by a GTP binding protein, EF-1. A guanine-nucleotide exchange complex, EF-1 $\beta\gamma\delta$ , replaces GDP by GTP on EF-1 $\alpha$ . A complex, purified from *Xenopus laevis* oocytes as a substrate for the meiotic and mitotic p34<sup>cdc2</sup> protein kinase, was shown to contain the guanine-nucleotide exchange activity. This complex is composed of three main proteins, p30, p36, and p47. This application describes the isolation of the EF-1 $\beta\gamma\delta$  and EF-1 $\beta\gamma$  complexes using UNO Q1 column.

### Purification

The traditional protocol to purify the EF-1 $\beta\gamma\delta$  and EF-1 $\beta\gamma$  complexes required six steps as outlined below.



The traditional separation protocol involved separation of the gel filtration desalted *Xenopus laevis* prophase oocyte sample on a strong anion exchange column (see Figure 1). The collected fraction contained both the EF-1 $\beta\gamma\delta$  (p47, p36 and p30) and EF-1 $\beta\gamma$  (p47, p30) complexes, a p40 contaminate and some high molecular weight contaminates (see Figure 2, Mono Q lane). The high molecular weight and p40 contaminates were

removed using a Heparin-Sepharose column (see Figure 2, Heparin lane). Fractionation of the heparin fraction on a DEAE column resulted in the separation of the EF-1 $\beta\gamma$  and EF-1 $\beta\gamma\delta$  complexes. (see Figure 2, DEAE lanes 1&3).

Purification of the EF-1 $\beta\gamma\delta$  and EF-1 $\beta\gamma$  complexes using the UNO Q1 column required only five steps. The gel filtration fraction was separated on the UNO Q1 column (see Figure 3) at 4.5 ml/min. This was accomplished at **3x** faster flow rate than that of the traditional strong anion exchange column separation, (1.5 ml/min). Note the high degree of resolution achieved on the UNO Q1 column.

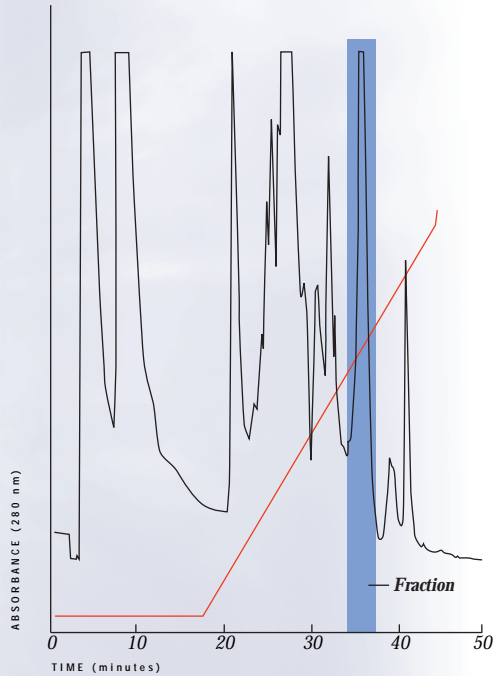
The collected fractions from the UNO Q1 separation were subjected to electrophoretic analysis. The EF-1 $\beta\gamma\delta$  complex was resolved into p47, p36 and p30 bands in fraction 16 and the EF-1 $\beta\gamma$  complex was resolved into p30 and p47 bands in fraction 14 (see Figure 4). The densitometry scan of fraction 16, showed that the p40 contaminate that was present in the Mono Q fraction was not present in the UNO fraction (data not shown). The high molecular weight contaminants in fraction 14 and 16 were removed using a heparin column (data not shown).

Incorporating the UNO Q1 column into the purification scheme of elongation factors from *Xenopus laevis* prophase oocytes resulted in improved purification and a reduction of the number of purification steps. This was achieved using an UNO Q1 high resolution separation at 4.5 ml/min.

### References:

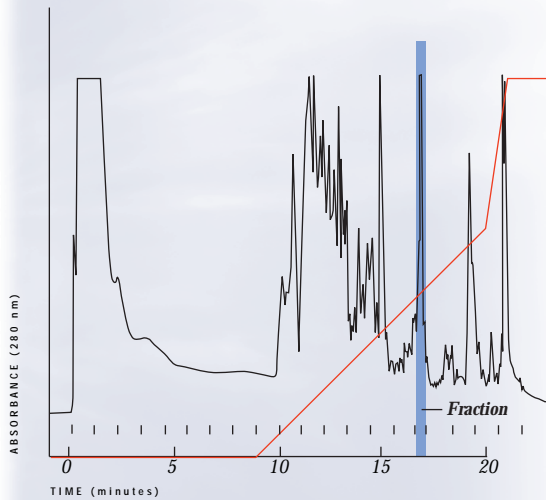
- Mulner-Lorillon O. et al., (1989), *FEBS Letters*, **251**, 219-224.
- Janssen, G. et al, (1991), *J. Biol. Chem.*, **266**, 23, 14885-14888.
- Mulner-Lorillon O. et al., (1994), *J. Biol. Chem.*, **269**, 31, 20201-20207.

**Figure 1. (NH<sub>4</sub>)<sub>2</sub>SO<sub>4</sub> Precipitate Fractionation of desalted *Xenopus laevis* oocytes on Strong Anion Exchanger.**



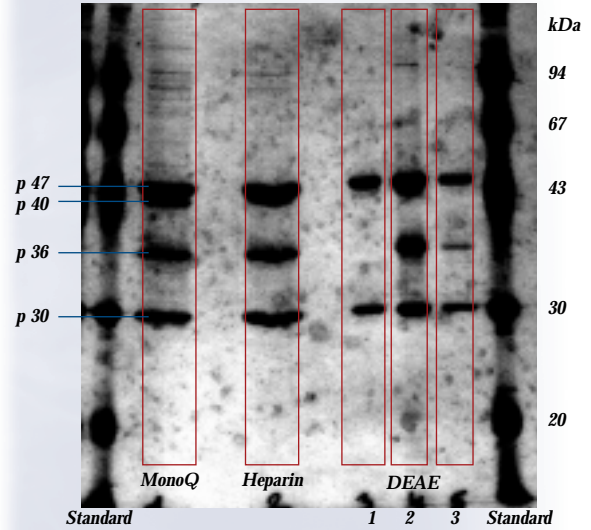
Protein load: 12 mg; Buffer A: 50 mM Tris, 1 mM DTT, 2 mM EDTA, pH 7.4; Buffer B: A + 1 M NaCl, Gradient: 0-70%B in 40 ml; Flow Rate: 1.5 ml/min System: FPLC®

**Figure 3. Fractionation of desalted *Xenopus laevis* oocytes extract on UNO Q1.**

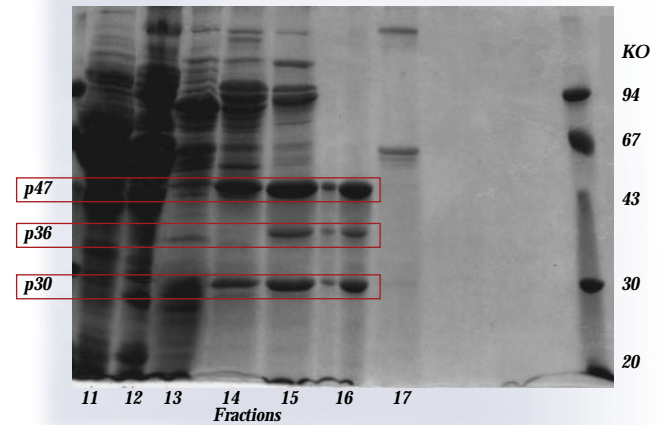


Protein load: 6 mg; Buffer A: 50 mM Tris, 1 mM DTT, 2 mM EDTA, pH 7.4; Buffer B: A + 1 M NaCl, Gradient: 0-60%B in 50 ml; Flow Rate: 4.5 ml/min; System: FPLC®

**Figure 2. SDS-PAGE of Mono Q, Heparin Sepharose, and DEAE column fractions containing EF-1βγ and EF-1βγδ.**



**Figure 4. SDS-PAGE of UNO Q1 Fractions.**



Data courtesy of R. Poulhe, Laboratoire de la reproduction, CNRS, Université Pierre et Marie Curie, Paris, France.



**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