

Preparative 2-D Electrophoresis System Purifies Recombinant Nuclear Proteins from Whole Bacterial Lysates

Gaëton Gagnon, A. Marie-Josée Sasseville, and Yves Raymond, Institute du Cancer de Montréal, Montréal, Québec Canada H2L 4M1

Introduction

This report describes the 2-step isolation of recombinant nuclear proteins from whole bacterial cell lysates. The Rotofor® preparative isoelectric focusing cell was used in the first step of purification. Subsequent purification on a molecular weight basis was accomplished using Bio-Rad's Model 491 prep cell. These 2 techniques, used in combination, completely isolated recombinant proteins from bacterial contaminants.

The nuclear lamins are a family of structural proteins in the range of 60 to 70 kD (Figure 2, lane L) that are found in eukaryotic nuclei. Type A lamins have nearly neutral isoelectric points and are composed of lamins A and C in mammalian cells. Lamins A and C share an extensive portion of their sequence and differ only by the length of their carboxy termini (McKeon et al. 1986, Fisher et al. 1986). Type B lamins have acidic isoelectric points. Nuclear lamins have been identified as targets for human autoantibodies, but the extent to which these reactions are specific to certain autoimmune diseases has not been completely clarified (Senecal and Raymond 1992).

Further studies of the specificity of autoantibodies require an abundant source of highly purified lamins A, B, and C of human origin. For this purpose, prokaryotic expression vectors were prepared harboring individual cDNAs for human lamins A and C (McKeon et al. 1986) and B (Pollard et al. 1990). Preliminary experiments showed that the bulk of bacterial proteins have isoelectric points below pH 7. During isoelectric focusing, most bacterial proteins focus in the acidic region of the pH gradient, well separated from recombinant human lamins A and C, which focus toward the alkaline end of the pH gradient. We therefore envisaged the use of the Rotofor cell as a primary fractionation step for preparative purification of these lamins from whole bacterial lysates.

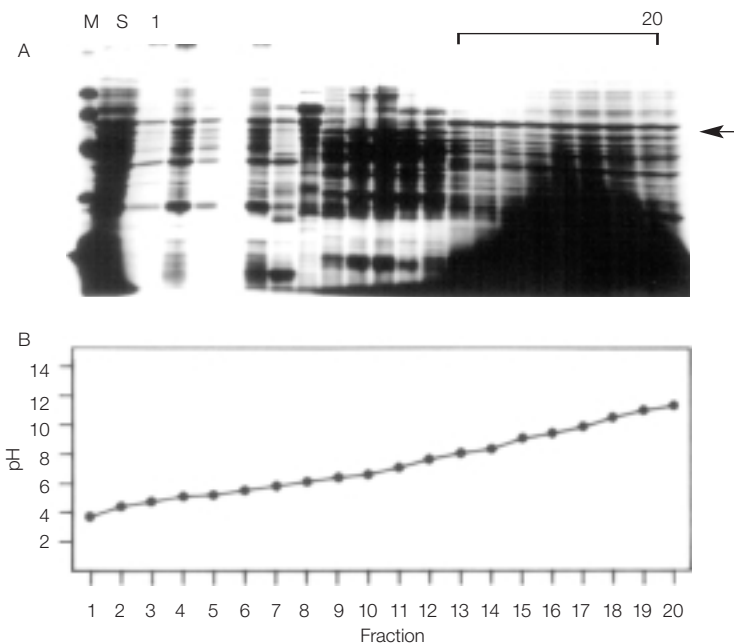


Fig. 1. Purification of recombinant human lamin C by preparative isoelectric focusing. A, 10% polyacrylamide gel of the starting material (whole bacterial lysate, lane S) and of Rotofor fractions 1–20 after staining with Coomassie Blue. The position of lamin C is indicated by the arrow. Fractions 13–20 were pooled for further purification on the Model 491 prep cell. B, pH gradient for Rotofor fractions 1–20.

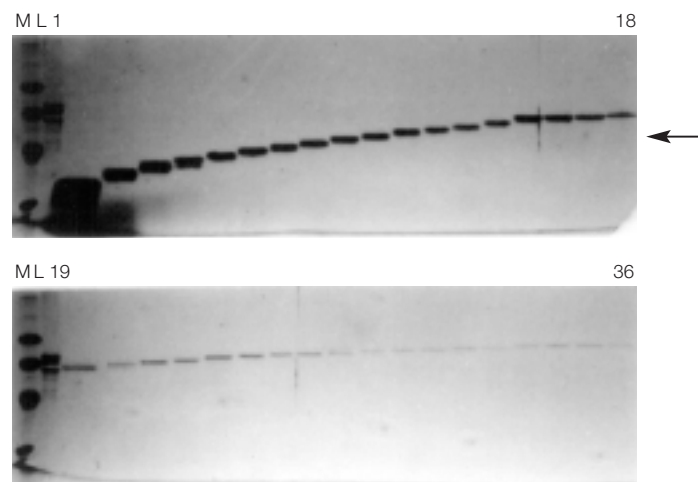


Fig. 2. Further purification of proteins. The contents of every third fraction eluted from the prep cell were analyzed on two 8% polyacrylamide gels. Lane M, molecular weight markers; lane L, rat liver nuclear lamins used as markers (lamins A, B, and C from top to bottom).

Methods

Protein Production in *E. coli*

DNA sequences coding for human lamin C were subcloned in the *NcoI*-*Bam*HI cloning site of a pET-11d translation vector bearing the bacteriophage T7 promoter (Studier et al. 1990). Plasmids were transformed into a BL21 (DE3)pLysS host (purchased from Novagen, Inc.) Bacteria were grown in Luria broth in the presence of carbenicillin at 100 µg/ml and chloramphenicol at 34 µg/ml until they reached an A₆₀₀ of 0.6. Synthesis of the recombinant lamin C was induced by the addition of IPTG at 0.4 mM and cultures were incubated for a further 3 hr at 37°C. Bacteria were harvested by centrifugation at 6,000 x g for 15 min and pellets were processed immediately or stored at -80°C.

Isoelectric Focusing on the Rotofor Cell

A bacterial pellet equivalent to 400 ml of the induced culture was dissolved in 7 ml of 10 M urea, extensively sonicated with a tip sonifier, and ultracentrifuged at 70,000 x g for 30 min. The supernatant, containing approximately 120 mg of protein, was adjusted to 8 M urea and 1% Bio-Lyte® ampholytes (pH range 3–10) in a final volume of 55 ml. Focusing was carried out at 12 W constant power for 4.5 hr at room temperature with cooling provided by running tap water. The initial conditions were 400 V and 30 mA. At equilibrium the conditions were 1,800 V and 6.5 mA. Twenty fractions of 2.5 ml each were collected, their pH measured, and their polypeptide composition analyzed by SDS-PAGE. See Figure 1.

Preparative Gel Electrophoresis on the Model 491 Prep Cell

Rotofor fractions containing lamin C identified by Coomassie Blue staining (Figure 1A) and immunoblotting (not shown) were pooled and dialyzed extensively against 10 mM Tris-HCl, pH 7.5, and then distilled water at 4°C. Proteins were concentrated by lyophilization. The dried protein, approximately 6 mg, was dissolved in 2 ml of SDS sample buffer and applied to the Model 491 prep cell for purification. The gel composition was 7%T/2.67%C, with a height of 10 cm and a gel tube size of 28 mm. The stacking gel composition was 4% with a height of 2.1 cm. Running conditions were 40 mA constant current (180–230 V) for 8–10 hr. After elution of Bromophenol Blue tracking dye, 2.5 ml fractions were collected. The polypeptide composition of every third fraction was analyzed by SDS-PAGE. Lamin C was identified by molecular weight (Figure 2) and antibody reactivity (not shown).

Results

The Rotofor cell was extremely efficient for initial purification of lamin C from bacterial proteins. The SDS-PAGE gel shown in Figure 1A demonstrates how the Rotofor cell separated the bulk of the contaminating proteins in the lysate from the recombinant lamin C-containing fractions. A similar separation profile was obtained for recombinant lamin A (not shown) which displays the same isoelectric point as lamin C.

Figure 1A demonstrates that Rotofor fractions containing lamin C also contained higher and lower molecular weight contaminants. Therefore, further purification by molecular weight was performed using the Model 491 prep cell. The efficiency of this apparatus at separating polypeptides is clearly illustrated in Figure 2. Lamin C was recovered in a small number of fractions, completely isolated from bacterial contaminants.

The isolation of lamin C by a method using denaturation by urea and ionic detergents is not a problem for the studies we are pursuing. First, nuclear lamins do not possess any enzymatic activity that could be irreversibly lost upon denaturation. Second, lamins, like other intermediate filament proteins, can be subjected to denaturation-renaturation cycles without loss of polymerization ability. And third, very efficient extraction methods exist whereby polypeptides can be completely recovered free of SDS, and we have used such preparations in immunofluorescence competition assays where the smallest trace of SDS would have adversely affected the antibodies under study (Collard et al. 1992).

In conclusion, a combination of isoelectric focusing on the Rotofor cell and molecular weight fractionation on the Model 491 prep cell allowed the complete purification of 1.5–2.0 mg of lamin C from a complex mixture of proteins equivalent to 400 ml of whole bacterial cultures in only 2 steps and with high yield.

References

- Collard JF et al., Redistribution of nuclear lamin A is an early event associated with differentiation of human promyelocytic leukemia HL-60 cells, *J Cell Sci* 101, 657–670 (1992)
- Fisher DZ et al., cDNA sequencing of nuclear lamins A and C reveals primary and secondary structural homology to intermediate filament proteins, *Proc Natl Acad Sci USA* 83, 6450–6454 (1986)
- McKeon FD et al., Homologies in both primary and secondary structure between nuclear envelope and intermediate filament proteins, *Nature* 319, 463–468 (1986)
- Pollard KM et al., In vitro posttranslational modification of lamin B cloned from a human T-cell line, *Mol Cell Biol* 10, 2164–2175 (1990)
- Senecal JL and Raymond Y, Autoantibodies to major and minor nuclear lamins are not restricted to autoimmune diseases, *Clin Immunol Immunopathol* 63, 115–125 (1992)
- Studier FW et al., Use of T7 RNA polymerase to direct expression of cloned genes, *Methods Enzymol* 185, 60–89 (1990)

Coomassie is a trademark of Imperial Chemical Industries PLC.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 507 6191
Canada (905) 712-2771 China (86-21) 63052255 Denmark 45 44 52-1000 Finland 358 (0)9 804 2200 France 01 47 95 69 65 Germany 089 318 84-177
Hong Kong 852-2789-3300 India (91-124) 6398112/113/114 Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460
Latin America 305-894-5950 Mexico 52 5 534 2552 to 54 The Netherlands 0318-540666 New Zealand 64-9-4152280 Norway 47-23-38-41-30
Russia 7 095 721 1404 Singapore 65-2729877 Spain 34 91 661 7085 Sweden 46 (0)8-55 51 27 00 Switzerland 061-717-9555 United Kingdom 0800-181134