

Fractionation of *Bovicola ovis* Homogenates Using the Mini Whole Gel Eluter—Effect of Buffer Composition on Elution Efficiency

Brian J. Shiell and Wojtek P. Michalski, CSIRO Division of Animal Health, Australian Animal Health Laboratory, Geelong, Victoria, Australia

Introduction

Amino acid sequence and composition analyses of proteins can be performed on samples derived from polyacrylamide gels. Protein electro-elution, the most commonly used method of recovering proteins from gels, can be performed on “cut-out” gel slices or using the Whole Gel Eluter recently introduced by Bio-Rad.¹⁻⁴ The Mini Whole Gel Eluter was used here to fractionate crude homogenate of the common sheep louse with the aim of characterizing immunogenic components.

Methods

Common sheep lice, *Bovicola ovis*, were harvested from infested animals, homogenized by ultrasonication, and the homogenate was clarified by centrifugation on a benchtop microfuge. Homogenate samples (1 mg protein) were separated by SDS-PAGE in 12% preparative well Tris-HCl Ready Gels (Bio-Rad). Protein fractions were immediately electro-eluted at 100 mA for 25 min using the Mini-Whole Gel Eluter. Three elution buffers: (A) 100 mM NH_4HCO_3 , 0.01% SDS, pH 8.5, (B) 10 mM phosphate buffer, 0.005% SDS, pH 7.8, and (C) 25 mM Tris, 19.2 mM glycine, 0.01% SDS, pH 8.3, were used to determine the elution efficacy. The fractions were analyzed in 12%, 15 well Tris-HCl Ready Gels (Bio-Rad) and proteins were visualized by

Coomassie® blue staining. Bio-Rad’s low molecular weight standards for SDS-PAGE are shown in Figure 1B, lane b.

Results and Discussion

Louse homogenate preparations are often rich in lipids and do not separate well on SDS-PAGE gels (Figure 1A, lane a). Individual proteins are poorly resolved and their extraction from individual “cut-out” bands is often inefficient. Several attempts were made to obtain amino acid sequence from various protein bands of louse homogenate using a variety of standard methods. This approach proved to be relatively ineffective and time-consuming.

The Mini Whole Gel Eluter was then used to fractionate crude louse homogenate. Samples of louse homogenate were separated on 12% preparative well gels and subjected to electro-elution. Two elution buffers, B and C, varying slightly in pH (7.8 v. 8.3) and in SDS content (0.005% v. 0.01%), were applied and in both experiments elution resulted in good recovery of a majority of proteins. It should be noted, however, that high molecular proteins did not elute well in buffer (B) containing low concentration of SDS (0.005%). The application of NH_4HCO_3 /SDS buffer (A) resulted in virtually no protein elution, possibly due to the low stability of ammonium carbonate buffer.

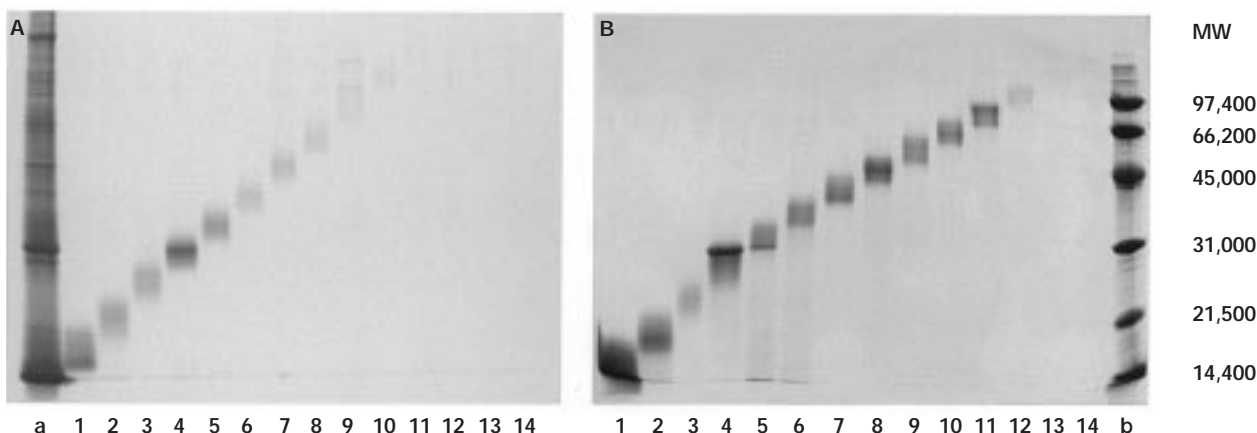


Fig. 1. SDS-PAGE profiles of protein fractions (components of sheep louse components) electro-eluted from mini SDS-PAGE gels using Mini Whole Gel Eluter. Gel A – electro-elution with buffer B: 10 mM phosphate buffer, 0.005% SDS, pH 7.8. Gel B – electro-elution with buffer C: 25 mM Tris, 19.2 mM glycine, 0.01% SDS, pH 8.3. Louse homogenate – gel A, lane a; Low molecular weight standards – gel B, lane b.

It was obvious that the buffer (C) containing Tris and a higher SDS concentration (0.01%) was a much better choice, as it resulted in the elution of virtually all proteins present in the louse homogenate ranging from <14,000 to >100,000 daltons. Increased SDS concentration also yielded approximately 2–3 fold greater amount of protein eluted in each of 14 fractions (Figure 1B).

Protein fractions electro-eluted on the Mini Whole Gel Eluter with buffer C, containing 0.01% SDS, pH 8.3, were transferred on Prosorb, desalted, and subjected to automated N-terminal amino acid sequencing. Of three major proteins analyzed, sequence information was obtained for two. One of the proteins contained a sequence that was similar to that of a reported allergen.

References

- 1 Andersen, P. and Heron, I., *J. Immunol. Meth.*, **161**, 29 (1993).
- 2 Dunbar, B. and Wilson, S. B., *Anal. Biochem.*, **216**, 227 (1994).
- 3 Andersen, P., Andersen, Å. B., Sørensen, A. L. and Nagai S., *J. Immunol.*, **154**, 3359 (1995).
- 4 Brandt, L., T. Oettinger, A., Holm, Andersen, Å. B. and P. Andersen, *J. Immunol.*, **157**, 3527 (1996).

Coomassie is a trademark of ICI.

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 86-10-62051850/51 **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** 39-02-216091 **Japan** 03-5811-6270 **Korea** 82-2-3473-4460
Latin America 305-894-5950 **Mexico** 514-2210 **The Netherlands** 0318-540666 **New Zealand** 64-9-4152280 **Russia** 7-095-4585822
Singapore 65-2729877 **Spain** 34-91-661-7085 **Sweden** 46 (0)8 627 50 00 **Switzerland** 01-809 55 55 **United Kingdom** 0800-181134