

January 01, 2004

**Keywords or phrases:**

His-Tagged Proteins, GFP,  $\beta$ -glucanase, Metal Affinity Chromatography, Membrane Adsorbers, Sartobind® Lab

# Purification of His-Tagged Proteins by Metal Affinity Chromatography with Sartobind® IDA Lab

Dr. Ricarda Busse<sup>1</sup>, Dr. Miyako Hirai<sup>1</sup>, Dr. Carsten Voss<sup>1</sup>, Dr. John Cashman<sup>2\*</sup>

1. Sartorius Stedim Biotech GmbH, August-Spindler-Straße 11, 37079 Göttingen, Germany

2. Sartorius UK Ltd., Blenheim Road, Epsom, KT19 9QQ, United Kingdom

\* Correspondence

Email: [john.cashman@sartorius.com](mailto:john.cashman@sartorius.com)

## Abstract

One of the practical methods to isolate recombinant proteins is to label them for purification by affinity chromatography. A common protein tag for this purpose is a peptide of six to eight histidine residues (His<sub>6</sub> or His<sub>8</sub>) cloned at the N- or C-terminus, which enables purification by metal chelate chromatography. This modality is based on the complexing of metal ions such as Ni<sup>2+</sup>, Cu<sup>2+</sup> or Co<sup>2+</sup>, immobilized on a chromatography matrix, by chelate formation through the imidazole ring of the amino acid histidine. The bound His<sub>6</sub>-tagged protein can subsequently be eluted from the matrix selectively with imidazole, which competitively binds to the metal complex. Sartobind® IDA utilizes a cellulosic membrane matrix, functionalized with iminodiacetic acid (IDA). This membrane therefore enables the immobilization of any suitable metal ion for metal affinity purification. In this study, Sartobind® IDA Lab units were evaluated for binding capacity, effectiveness, and handling in the purification of two hexa-histidine tagged proteins.

## Introduction

GFP is commonly used as a reporter protein. During the autocatalytic process, GFP forms a structure that emits green light at 508 nm if stimulated by light with wavelength of between 396 and 475 nm (absorbance maxima). Therefore, it is relatively easy to follow GFP during processing by means of fluorescence light. However, it becomes unstable or loses its fluorescent under certain conditions. GFP may be intracellular when produced in *E. coli* and therefore cells must be disrupted prior to separation.

The gene of  $\beta$ -glucanase is a fusion of glucanase genes from *Bacillus macerans* and *Bacillus amyloliquefaciens* [Borris *et al.*, 1990]. This fusion was cloned into a pET20b+ vector, while the signal sequence of  $\beta$ -glucanase was preserved. Transcription in *E. coli* was carried out under the IPTG-inducible T7 promoter. A secretion cassette is located on the plasmid. It consists of the gene for the bacterial release protein (kil) under the control of the weak stationary phase promoter *fic* [Miksch *et al.*, 1997]. The moderate expression of kil allows the secretion of the glucanase from the periplasm into the culture medium. The isolation and purification of  $\beta$ -glucanase can therefore be carried out after removal of *E. coli* from the fermentation medium.

Sartobind® IDA Lab offers the possibility to purify virtually any his-tagged target protein by affinity chromatography. Supplied uncharged, the user is free to choose the metal ion of their choice, to achieve optimum performance. Furthermore, the integral Luer connectors enable convenient use with syringes, while UNF 10-32 adapters included with the devices provide flexibility for operation by peristaltic pump or liquid chromatography systems. In this study, we demonstrate the performance of Sartobind® IDA Lab units (Figure 1) for metal affinity purification, using  $\text{Cu}^{2+}$  to isolate both GFP and  $\beta$ -glucanase carrying His<sub>6</sub> tags at their N-termini.



**Figure 1:** Sartobind® IDA Lab Membrane Adsorber Unit

## Materials and Methods

### Production of His<sub>6</sub>-GFP

BL21 *E. coli* were transfected with pHis<sub>6</sub>-GFP. Cells were cultivated in 300 mL medium (10 g/L peptone, 10 g/L yeast extract, 20 g/L glycerol, 10 g/L NaCl, 100 mg/L ampicillin) in a 1 L baffled flask, inoculated with 1 mL of starter culture. The culture was incubated with shaking at 37 °C overnight, with gene expression induced by IPTG after 2 hours. The cells were isolated by centrifugation, the resulting pellet (23.3 g) suspended in approximately 10 mL of lysis buffer (50 mM  $\text{NaH}_2\text{PO}_4$ , 300 mM NaCl, 10 mM imidazole, pH 8.0) and 25 mg lysozyme dissolved in 1 mL buffer was added (final concentration = 1 mg/mL). Cells were disrupted in an ice bath using pulsed ultrasound and the resulting lysate centrifuged. After removal of 5 mL for analysis, the supernatant was passed through a 0.2  $\mu\text{m}$  filter and the filtrate was taken forward for purification.

### Production of His<sub>6</sub>- $\beta$ -glucanase (His<sub>6</sub>-bgl)

BL21 (DE3) *E. coli* cells transfected with pET-bgl-His<sub>6</sub>-sec were cultivated under the same conditions as for His<sub>6</sub>-GFP production. The cells were harvested by centrifugation at 11,000 g for 30 min. A portion of the supernatant was passed through a 0.2  $\mu\text{m}$  filter and the filtrate was used for purification.

### Preparation of Sartobind® IDA Lab Unit

Each Sartobind® IDA Lab unit containing 9.3 cm<sup>2</sup> adsorption area (3x layers of 3.1 cm<sup>2</sup> membrane discs) was flushed with 10 mL of deionized water, which was slowly pumped from bottom to top until the unit was vented completely. Then the remaining water was applied to the top of the unit and allowed to flow through by gravity. The following solutions were passed through the unit in the same way:

1. 0.5 M  $\text{CuSO}_4$   
(charging buffer)
2. 10 mM imidazole in 25 mM potassium phosphate buffer, pH 8 (equilibration buffer 1)
3. 25 mM potassium phosphate buffer, pH 8  
(equilibration buffer 2)

### Protein Purification

The filtered lysate of His<sub>6</sub>-GFP (11 mL) and cell-free supernatant of His<sub>6</sub>-bgl (80 mL) were applied to the Sartobind® IDA Lab units and the flow through fractions collected. The units were washed twice with 10 mL of washing buffer (50 mM  $\text{NaH}_2\text{PO}_4$ , 300 mM NaCl, 20 mM imidazole, pH 8.0). Both fractions were collected and pooled. The proteins were eluted by applying 2x 3 mL of elution buffer (50 mM  $\text{NaH}_2\text{PO}_4$ , 300 mM NaCl, 250 mM imidazole, pH 8.0). Each eluate fraction was retained separately and all samples were stored at -20 °C until analysis.

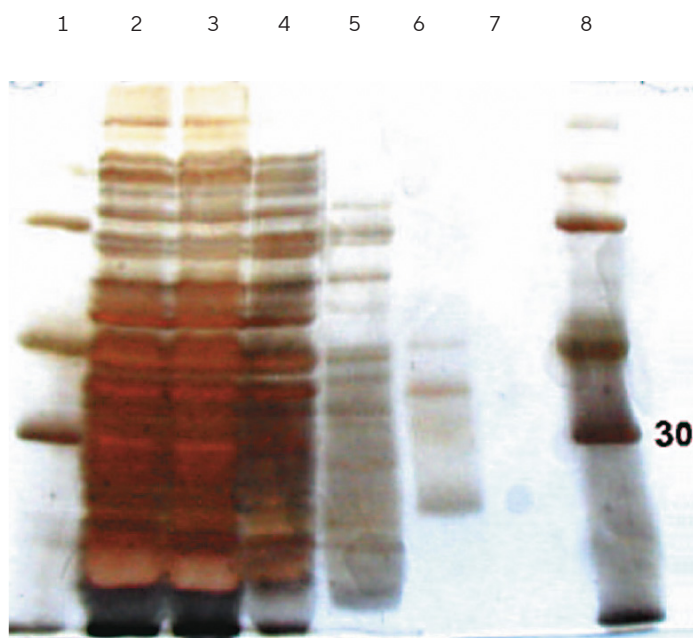
## Protein Analysis

Lysate | cell-free samples, and flow through, wash and eluate fractions were analyzed by polyacrylamide gel electrophoresis (PAGE) to assess purity. Protein activity was also assessed by fluorescence estimation for GFP or enzymatic activity of  $\beta$ -glucanase.

## Results and Discussion

### His<sub>6</sub>-GFP

Cell lysate (Figure 2, lanes 2 and 3) and Sartobind® IDA Lab flow through fraction (lane 4) show a broad range of intracellular proteins. Only the bands of abundant proteins are visible in the wash fractions (lane 5). Eluate 1 (lane 6) shows a clear band below 30 kDa, corresponding to His<sub>6</sub>-GFP with an expected molecular weight of 27.7 kDa. The two other bands are intracellular proteins which were also bound by the membrane adsorber – a common occurrence in purifications from whole cell lysates. Eluate 2 (lane 7) showed no detectable proteins, indicating that all bound proteins were recovered in the first elution step.



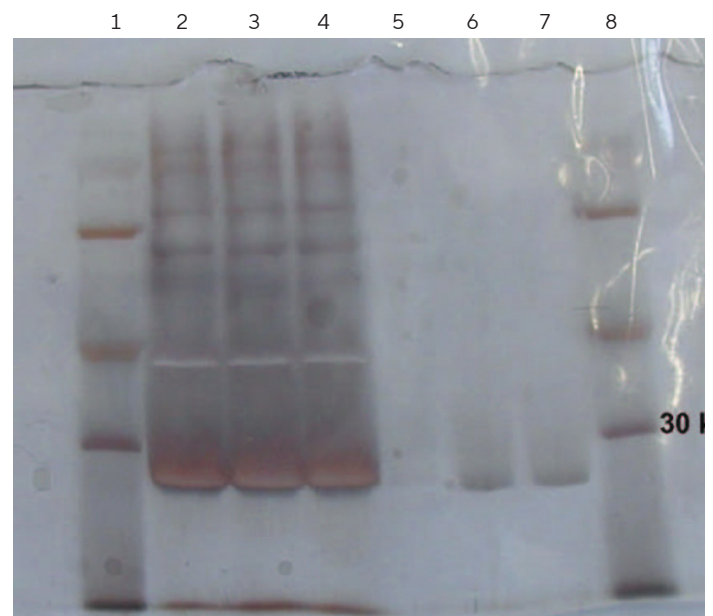
**Figure 2:** Polyacrylamide Gel Analysis of His<sub>6</sub>-GFP Purification Samples

Lane 1: Molecular weight standards  
Lane 2: Lysate  
Lane 3: Filtered lysate  
Lane 4: Flow through  
Lane 5: Wash fractions (pooled)  
Lane 6: Eluate 1  
Lane 7: Eluate 2  
Lane 8: Molecular weight standards

### His<sub>6</sub>- $\beta$ -glucanase

The samples of cell culture supernatant and Sartobind® IDA Lab flow through were concentrated prior to PAGE. These samples show a distinct, broad band migrating to <30 kDa (Figure 3, lanes 2-4), which corresponds to a high quantity of  $\beta$ -glucanase. The high quantity of this target protein in the flow through suggests that the amount of  $\beta$ -glucanase exceeded the binding capacity of the Sartobind® IDA Lab unit. The pooled wash fractions (lane 5) show no detectable protein, indicative of stable binding of the target molecule to the membrane adsorber. Each eluate (lanes 6 and 7) displays a single band of the target protein. This indicates that the membrane adsorber unit was completely saturated with the target protein, in keeping with the previously mentioned detection of a high quantity of  $\beta$ -glucanase in the flow through. No additional proteins were detected in the eluates.

The protein concentration in eluates 1 and 2 was 232 and 127  $\mu$ g/mL, respectively, corresponding to a total protein content of 1,077  $\mu$ g in the 2x 3 mL fractions. This places the binding capacity of Sartobind® IDA Lab for  $\beta$ -glucanase at approximately 1.1 mg, although we assume a third eluate would have yielded further protein and therefore this represents the minimum capacity of the unit (118  $\mu$ g/cm<sup>2</sup> or 4.3 mg/mL, where 1 mL membrane = 36.4 cm<sup>2</sup>).



**Figure 3:** Polyacrylamide Gel Analysis of His<sub>6</sub>-bgl Purification Samples

Lane 1: Molecular weight standards  
Lane 2: Lysate  
Lane 3: Filtered lysate  
Lane 4: Flow through  
Lane 5: Wash fractions (pooled)  
Lane 6: Eluate 1  
Lane 7: Eluate 2  
Lane 8: Molecular weight standards

## Conclusion

Our results show that, when charged with Cu<sup>2+</sup>, Sartobind® IDA Lab units are suitable for the purification of both intra- and extra-cellular proteins. The units exhibited high binding capacities for both His<sub>6</sub>-GFP and His<sub>6</sub>-β-glucanase, and both proteins were eluted at high purity.

## References

Borris, R., Buettner, K. and Maentsaelae, P. (1990). Structure of the beta-1,3-1,4-glucanase gene of *Bacillus macerans*: homologies to other beta-glucanases. *Molecular & general genetics* 222 (2-3), 278-283.

Miksch, G., Fiedler, E., Dobrowolski, P., Fries, K. (1997). The *kil* gene of the ColE1 plasmid of *Escherichia coli* controlled by a growth-phase-dependent promoter mediates the secretion of a heterologous periplasmic protein during the stationary phase. *Arch. Microbiol.* 167 (2+3), 143-150.

This Application Note is made available courtesy of Friehs K., Bielefeld University: Isolierung rekombinanter Proteine mit Membranmodulen. Bericht.

## Note


Literature published up to c.2022 may reference the use of Sartobind® MA, which is a name previously used for the Sartobind® Lab membrane adsorbers. When these products were renamed, there was no change made to fit, form or function. Therefore, results collected and methods established using Sartobind® MA remain valid also for Sartobind® Lab.

### Germany

Sartorius Lab Instruments GmbH & Co. KG  
Otto-Brenner-Straße 20  
37079 Göttingen  
Phone +49 551 308 0

### USA

Sartorius Corporation  
565 Johnson Avenue  
Bohemia, NY 11716  
Phone +1 631 254 4249  
Toll-free +1 800 635 2906

 For further information, visit  
[www.sartorius.com](http://www.sartorius.com)