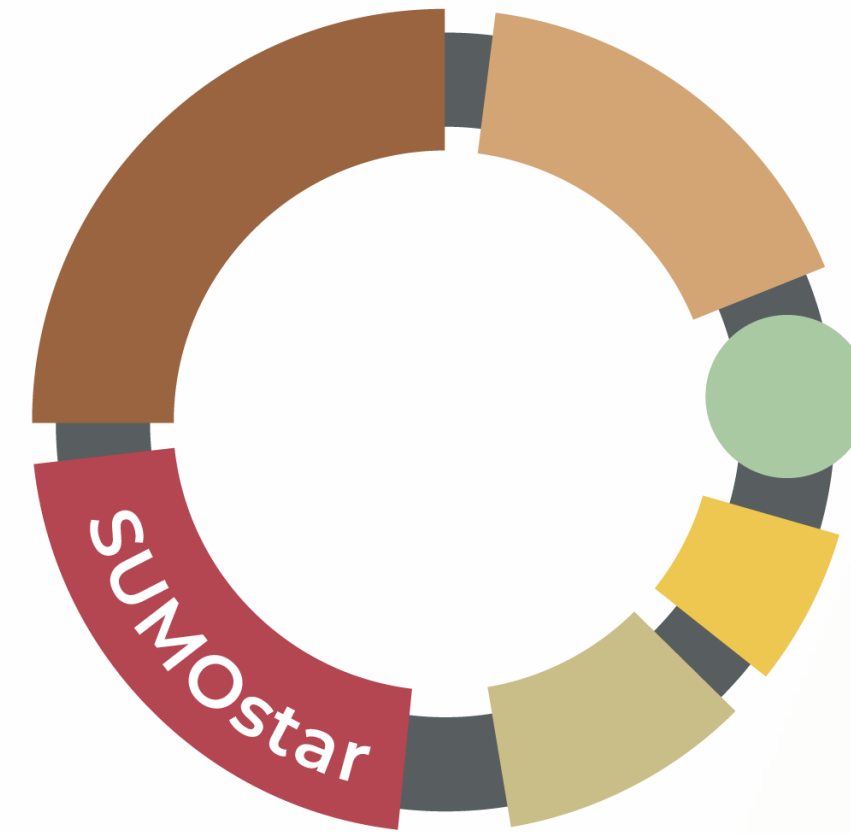


# Manufacturing Single Chain Antibodies, Nanobodies with SUMO Expression Platform



LifeSensors Inc.  
271 Great Valley Parkway  
Malvern PA 19355

610-644-8845  
bd@lifesensors.com  
www.lifesensors.com

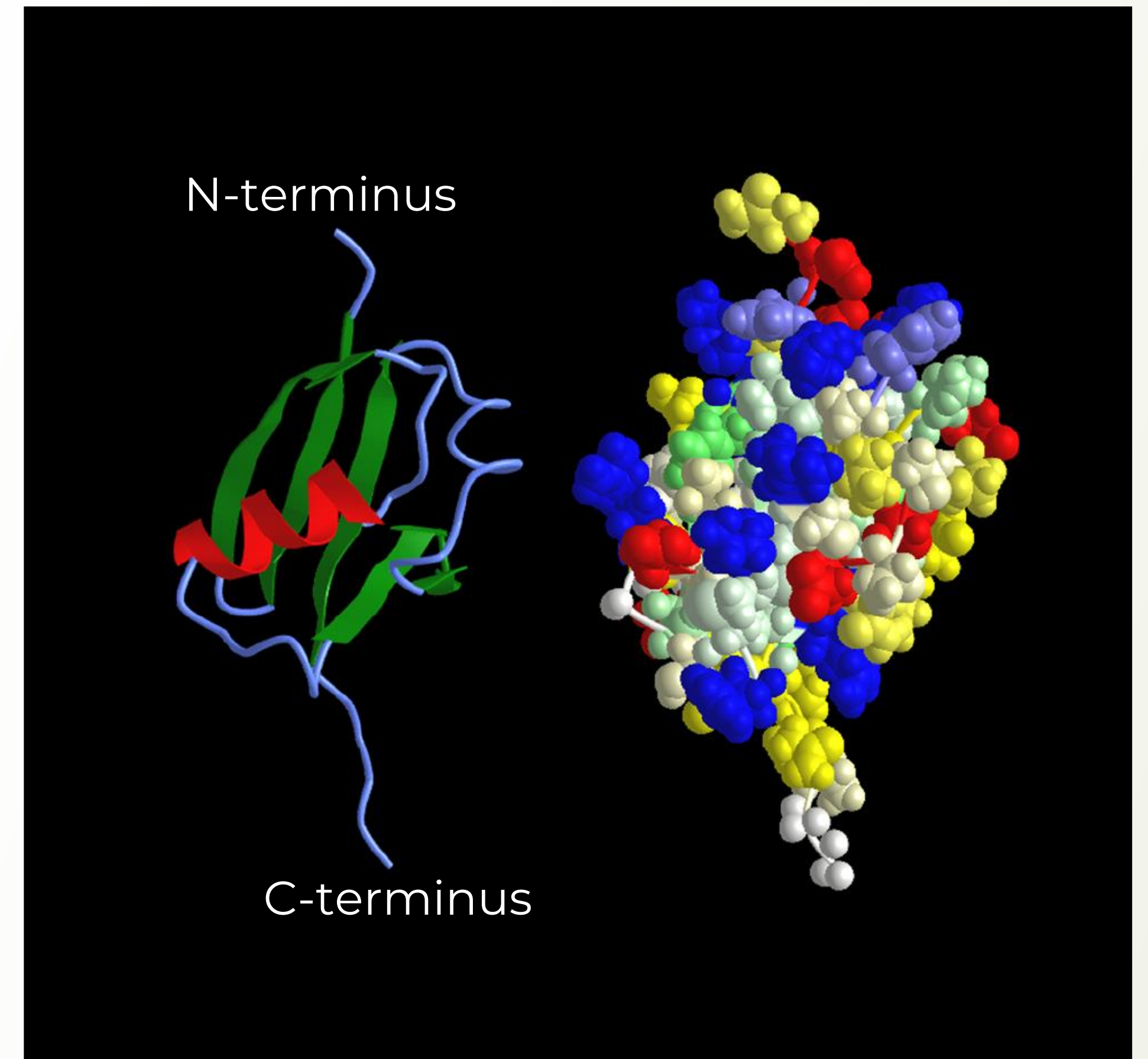
# SUMO Advantage

- Nanobodies or engineered single chain antibodies don't express well in E.coli
- Clinical programs are abandoned due to poor expression
- More than 90 proteins expressed in E.coli have been approved by FDA †
- SUMO system dramatically enhances expression of soluble nanobodies and decreases the cost of goods

† Full-length recombinant antibodies from E.coli --- clinical evaluation. MABS, 14, 1, e2111748 (20 pages)

# How Does SUMO Help Enhance Nanobody Expression?

- Nature designed SUMO to act as a chaperone for proteins
- SUMO, or Small Ubiquitin-like MOdifier, is a member of the ubiquitin family
- Flexible N-terminal region followed by a compact ball-shaped ubiquitin-like fold
- SUMO improves solubility due to hydrophilic outer surface and hydrophobic core



# SUMO Platform for Nanobodies

## PROBLEMS

Unstable protein,  
low expression

Insoluble Nanobodies

Miss-folded nanobodies

High cost of goods



## SOLUTIONS

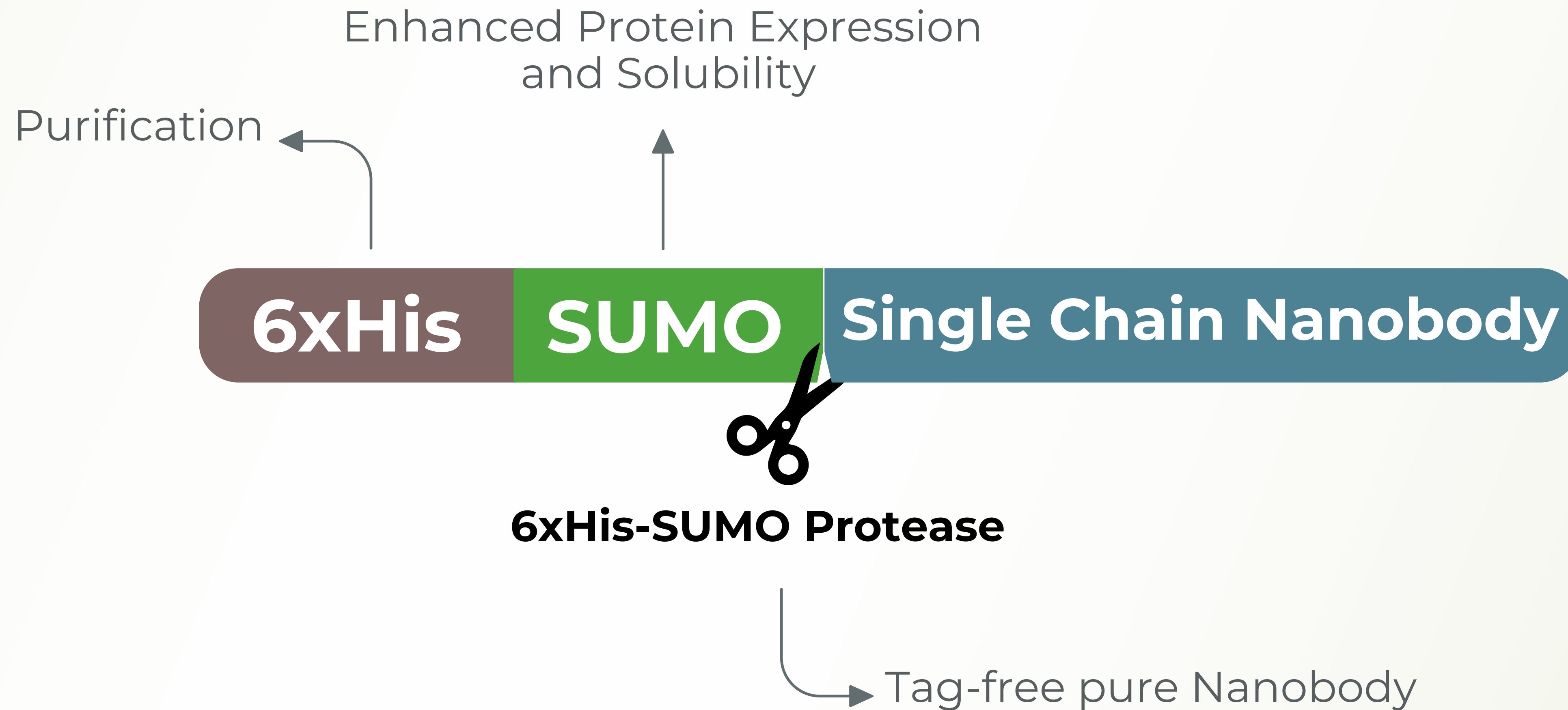
SUMO chaperoning,  
enhanced expression

SUMO derived solubility

SUMO-driven folding

High yield, saves cost

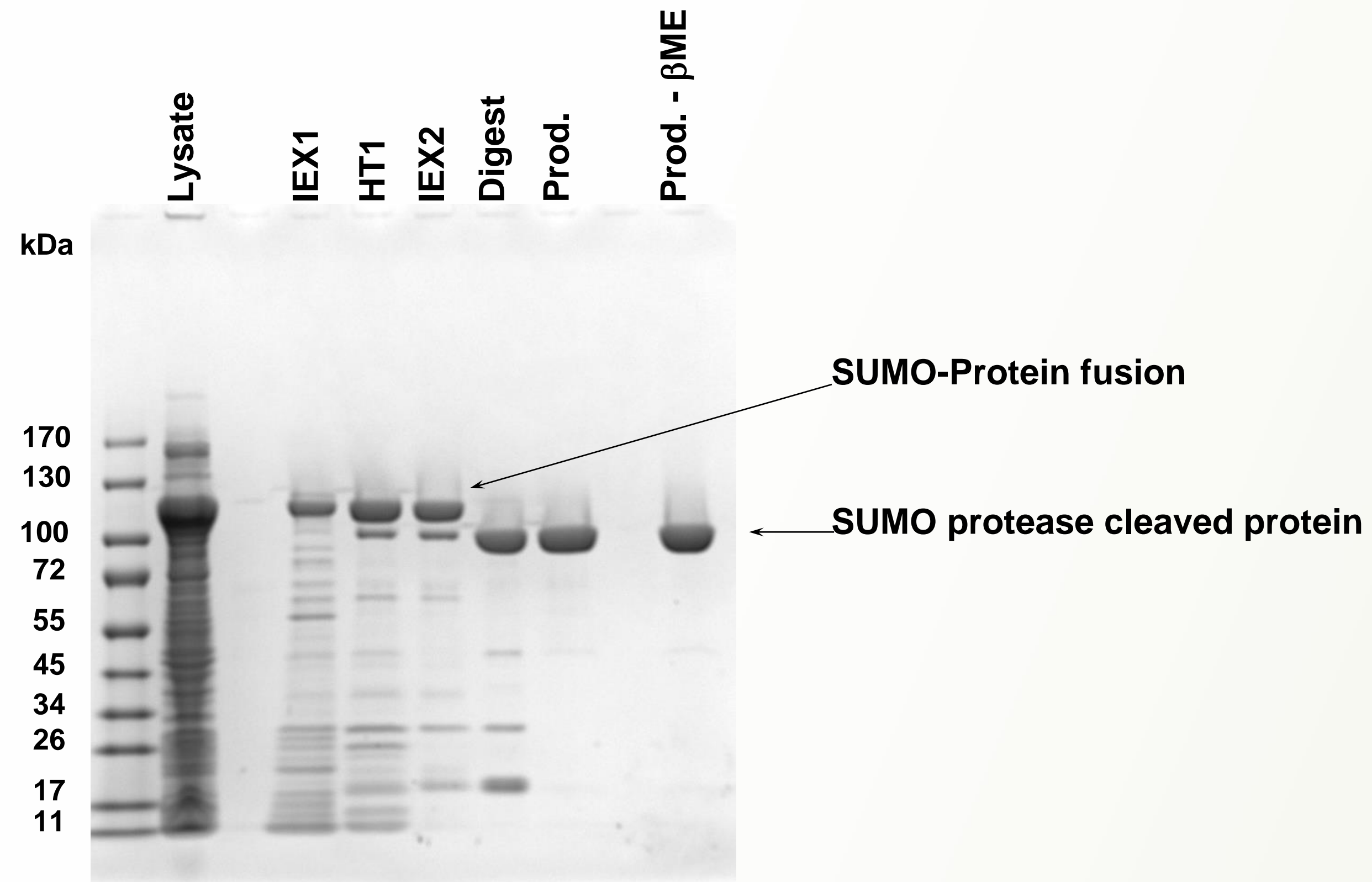
# SUMO Platform For Native Nanobodies





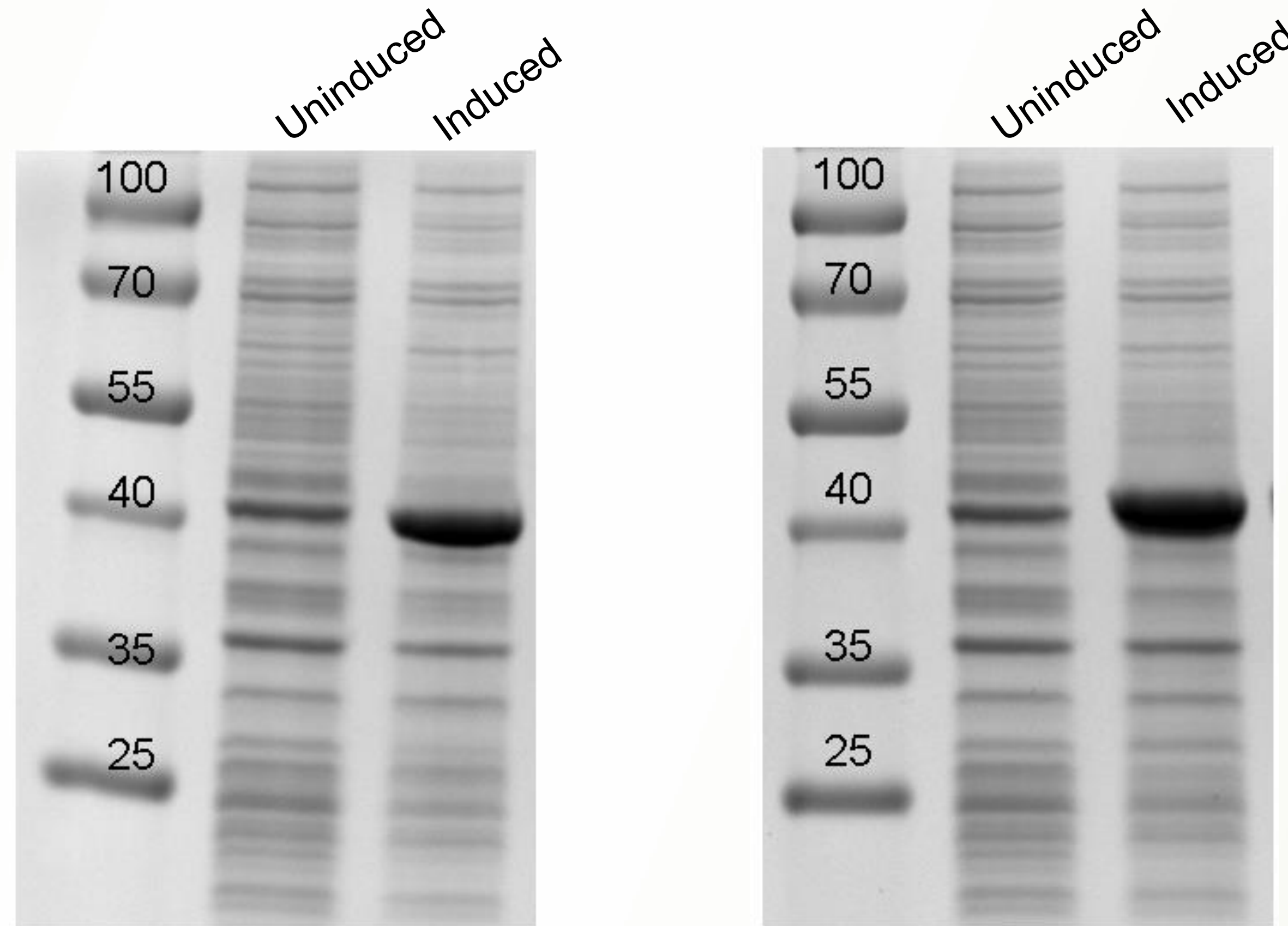
# Expression of Candidate Protein in *E. coli*

- Improved expression of Protein from 8% TCP to >20% TCP
- Purification reduced from 5 column steps to 4
- Overall yield improved from ~16% to >40%





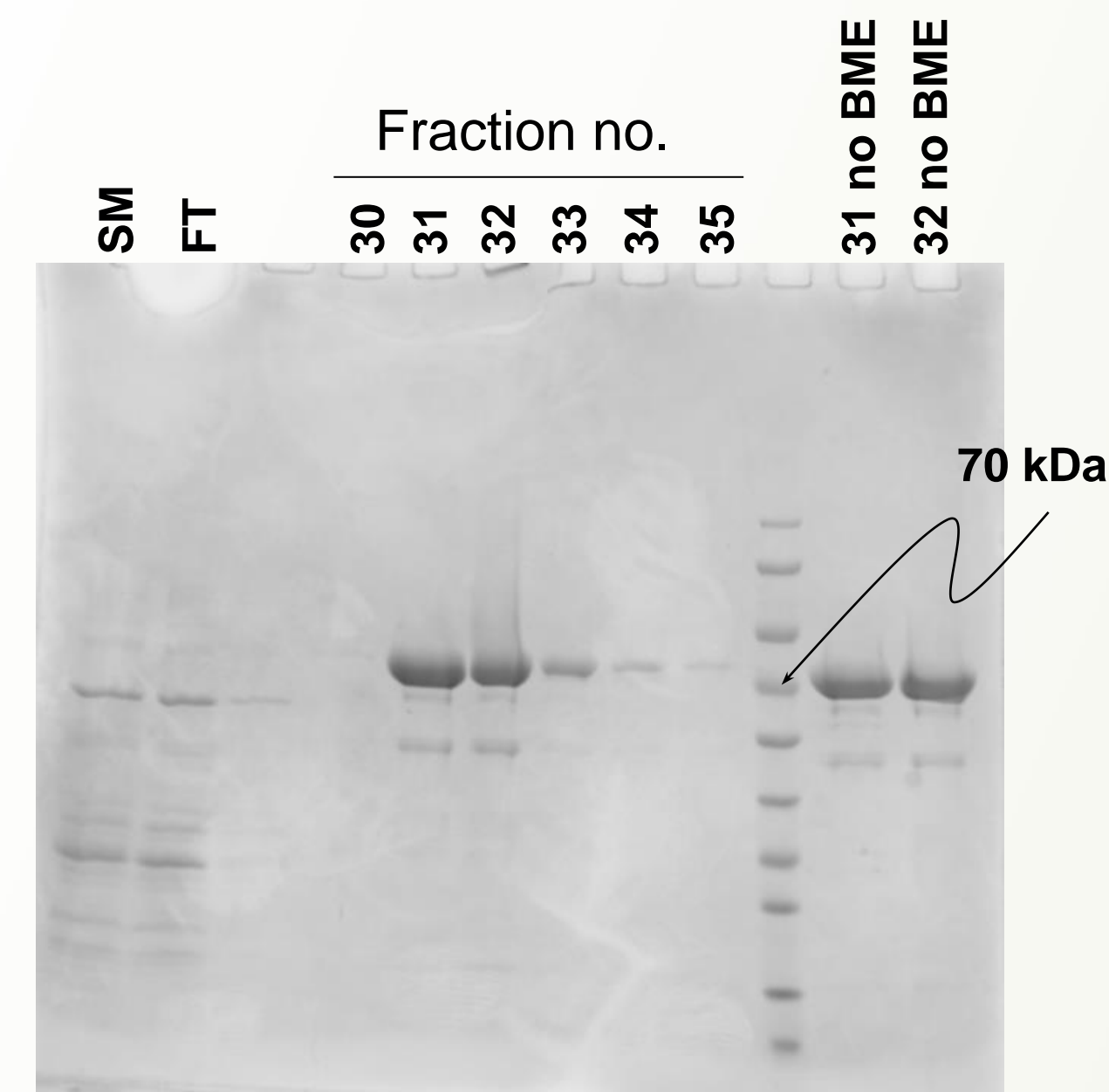
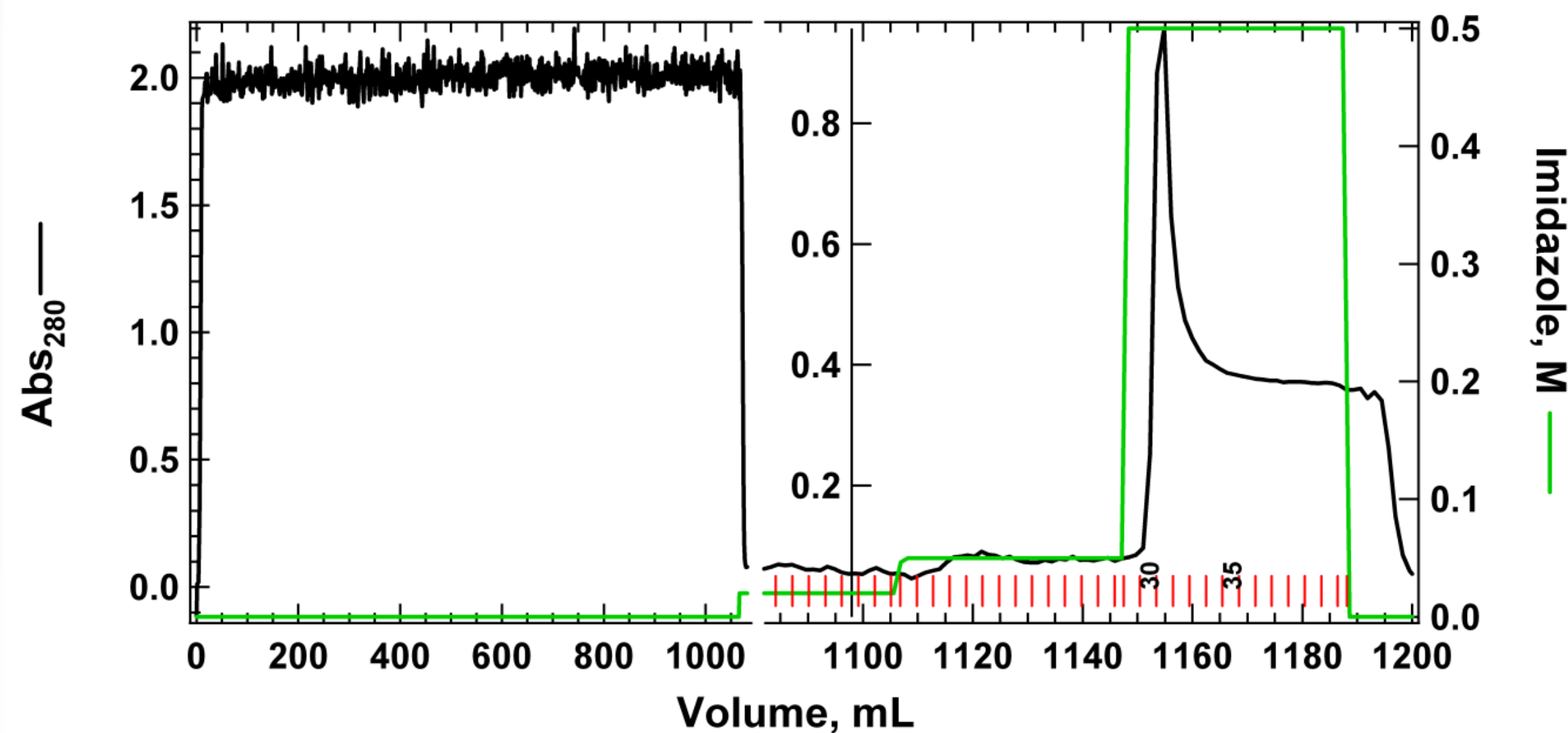
# Expression of single chain antibodies as SUMO fusions in *E. coli*



# scFv<sub>2</sub> IMAC purification



## 0.8 M Arg hypertonic periplasmic extract



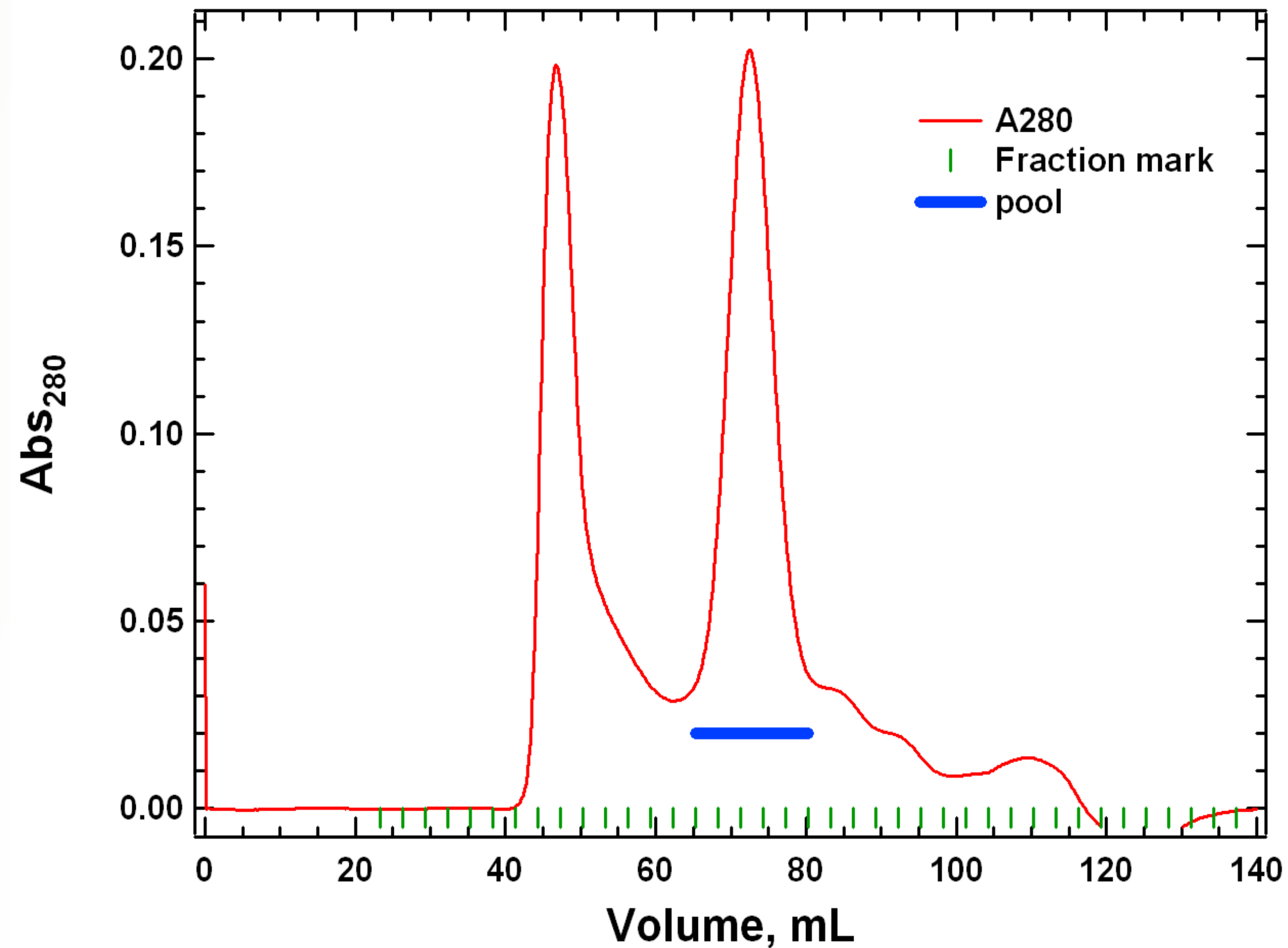
Coomassie blue stained gel.

8  $\mu$ L of 3 mL fraction analyzed





## SEC analysis of SUMO-scFv<sub>2</sub> (S200 10/30)

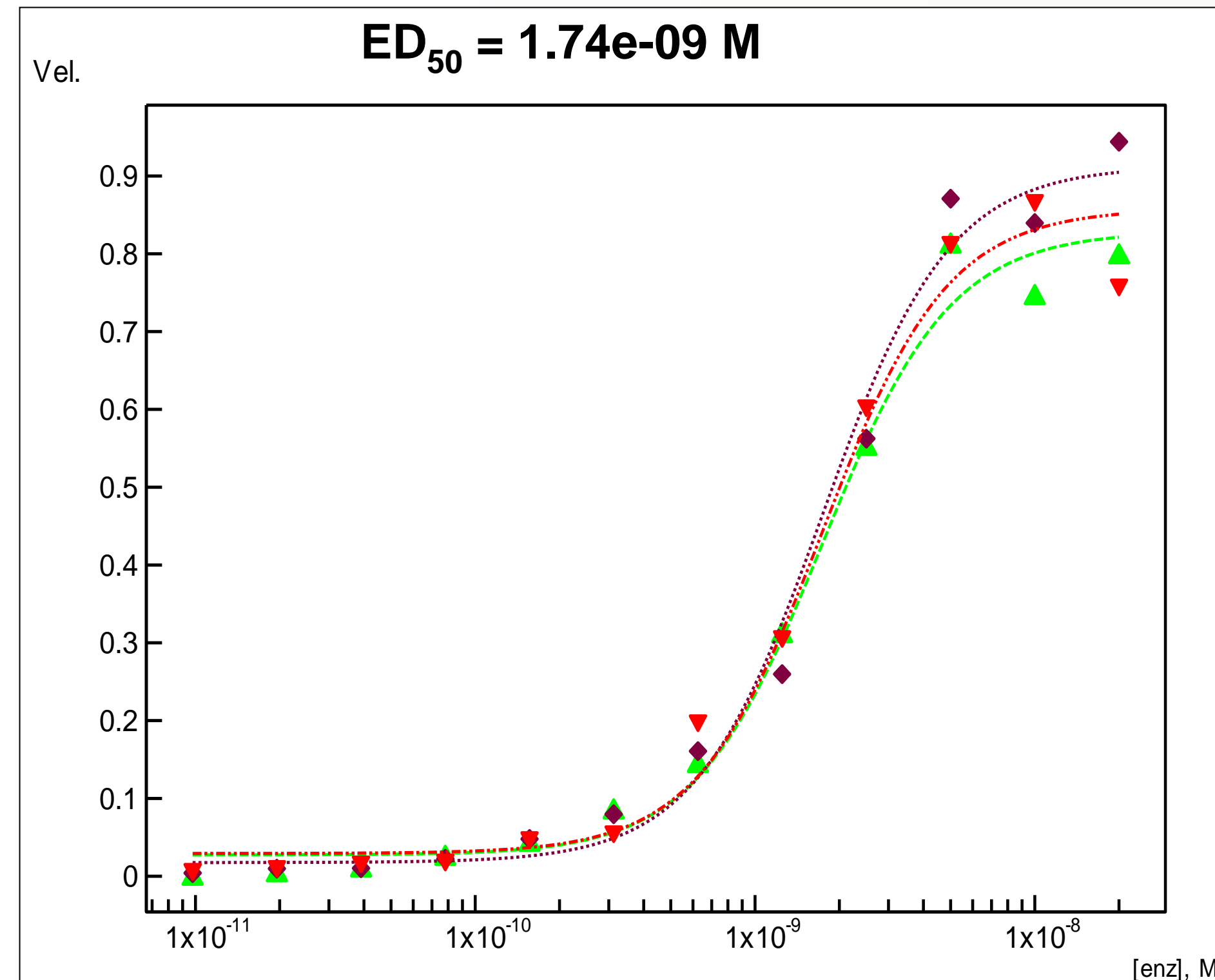
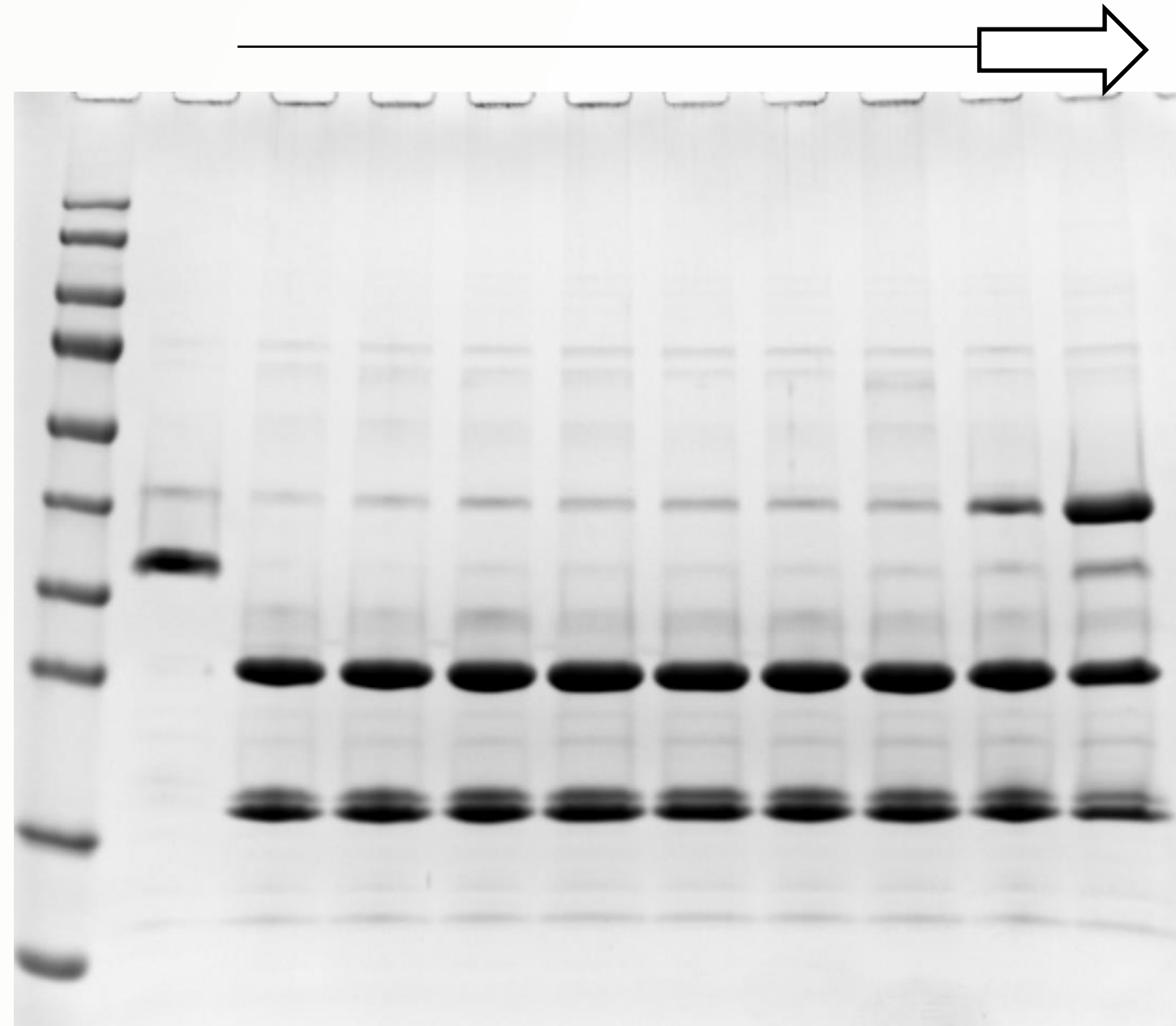


*SUMO-scFv<sub>2</sub> elution profile consistent with dimer formation*



# Remarkably Efficient Removal of SUMO by SUMO-protease

Decreasing conc. of SUMO protease



10 ug SUMO-GFP was incubated with 0.1-1.00 unit of [SUMO protease](#) for 37 for one hour. SDS PAGE was stained with commissive blue. Data shows that LifeSensors SUMO protease is highly active. See next slide for comparison ion between different proteases



# Comparison Between SUMO Protease and Commonly used Tag Removal Enzymes

Protease	Cleavage site	Location	Residual AAs	pH range/ optimum	Chaotrope Sensitivity	Salt Sensitivity	Enz: target	\$/mg
TEV, AcTEV, ProTEV	ExxYxQ/(G/S)	N	G/S	5.5-8.5/7	2 M urea	≤ 0.1 M	1-3% (wt/wt)	36 <sup>1</sup>
		C	ExxYxQ					
Thrombin (fIIa)	LVPRG	N	G	6-9	≤ 0.1 M urea	≤ 0.15 M	1-10% (wt/wt)	3.5 <sup>2**</sup>
		C	LVPR					
Factor Xa (fXa)	IEGR/x	N	none	6-9	≤ 0.1 M urea	≤ 0.15 M	1-10% (wt/wt)	26.9 <sup>3*</sup>
		C	IEGR					
Enterokinase (EK)	DDDDK/x	N	none	7-8	N/A	N/A	0.1% (wt/wt)	8.2 <sup>4**</sup>
		C	DDDDK					
<b>SUMO proteases</b>	<b>No sequence preference</b>	<b>N only</b>	<b>Encode desired seq</b>	<b>6-9</b>	<b>0.1 M GnCl &lt; 2 M Urea</b>	<b>&lt; 0.5 M</b>	<b>0.1% (wt/wt)</b>	<b>10<sup>5</sup></b>

- <https://www.fishersci.com/shop/products/promega-protev-plus-protease-2/p-4389612>
- <https://www.sigmaaldrich.com/US/en/product/sigma/ge27084601>
- [https://www.neb.com/products/p8010-factor-xa-protease#Product%20Information\\_Properties%20&%20Usage](https://www.neb.com/products/p8010-factor-xa-protease#Product%20Information_Properties%20&%20Usage)
- <https://www.neb.com/products/p8070-enterokinase-light-chain#Product%20Information>
- <https://lifesensors.com/product/sp4010-sumo-protease-1/>

\*Assumes a **6 hour** reaction as opposed to a 1 hour reaction.

\*\*Assumes a **16 hour** reaction as opposed to a 1 hour reaction

Yields are based upon shake flask growth and are unoptimized.

# Examples of Nanobodies Expressed as SUMO

- **Fibronectin type III domain based monobodies:** Monobodies were expressed and purified using SUMO fusion system in E.coli (Gorman K et al, Curr Protoc Chem Biol. 2018; Hussain M et al, Biochemistry. 2018)
- **Fab fragments:** Cytosolic expression of functional fab fragments in Escherichia coli using a novel combination of dual SUMO expression cassette and EnBase® cultivation mode (Rezaie F et al, J Appl Microbiol. 2017)
- **Single-chain variable fragment (scFv):** High-level expression of an anti-VEGF165 scFv in E. coli achieved by SUMO fusion (Sillen M et al, Int J Mol Sci. 2020)
- **Nanobodies (Nbs):** Nbs against a GPCR, PAI and SARS-CoV-2 RBD were expressed and purified in E.coli (Koehl A et al, Nature. 2019; Sillen M et al, Int J Mol Sci. 2020; Shi Z et al, Structure. 2022; Sillen M et al, J Thromb Haemost. 2020; Sillen M et al, Int J Mol Sci. 2021)
- **VHH antibodies:** Single-Domain VHH antibodies against FMDV proteins and Chikungunya E2 protein were expressed and purified with SUMO fusion system (Li H et al, Vaccines (Basel). 2021; Li Q et al, J Nanobiotechnology. 2022; Wang D et al, BMC Vet Res. 2015)
- **Heavy-chain-only antibodies (HCAbs):** HCAbs against influenza HA protein were produced in E.coli fused with SUMO.

# SUMO Advantage

- **Dramatic Enhancement of Nanobodies yield as SUMO-fusion**
- **Application of the most efficient SUMO protease to generate native proteins**
- **More than 90 proteins expressed in E.coli have been approved by FDA †**
- **Application of SUMO system decreases the cost of goods**

† Full-length recombinant antibodies from E.coli --- clinical evaluation. MABS, 14, 1, e2111748 (20 pages)

# Thank You

We are your partner in Manufacturing Difficult to Express Proteins

## Contact Us!

<b>Research &amp; Product Inquiries</b>	<b>R&amp;D</b>	<b><a href="mailto:info@lifesensors.com">info@lifesensors.com</a></b>	<b>610-644-8845 (ext 339)</b>
<b>Custom Service &amp; Assays</b>	<b>BD</b>	<b><a href="mailto:bd@lifesensors.com">bd@lifesensors.com</a></b>	<b>610-644-8845 (ext 310)</b>