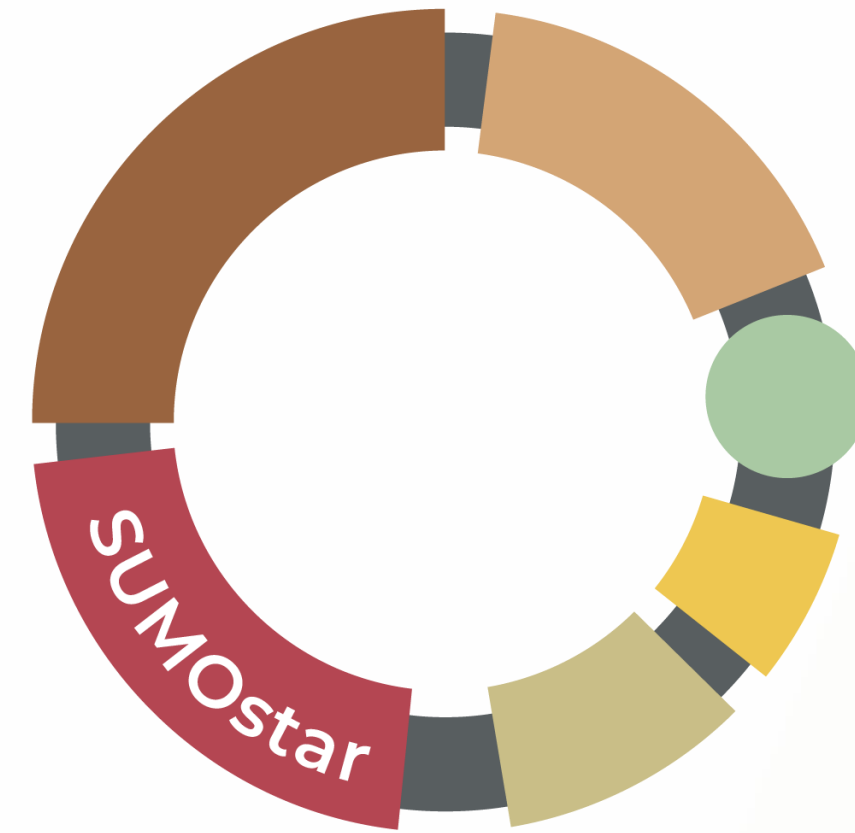


# Therapeutic Proteins Manufacturing with SUMO Expression Platform



LifeSensors Inc.  
271 Great Valley Parkway  
Malvern PA 19355

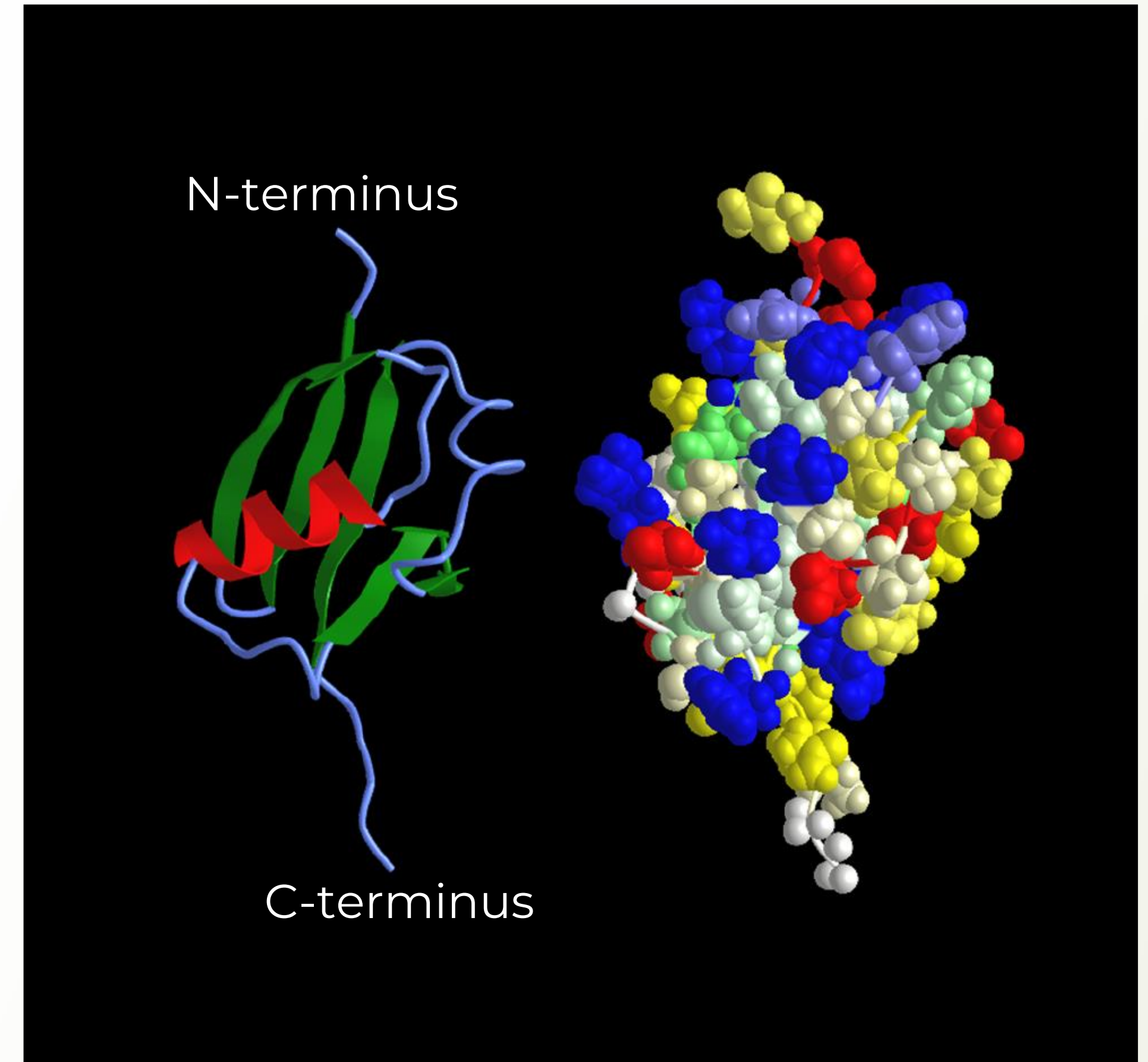
610-644-8845  
bd@lifesensors.com  
www.lifesensors.com

# SUMO Advantage

- Efficient translation of proteins is key to the success of therapeutics, gene and cell therapy and mRNA therapy
- Poor expression of engineered genes leads to failure in the clinic
- SUMO system dramatically enhances expression, and provides unique features not possible with traditional systems
- More than 150 references of therapeutic proteins expressed as SUMO-fusion in E.coli and mammalian cells

# What is SUMO?

- SUMO, or Small Ubiquitin-like modifier, is a member of the ubiquitin family
- SUMO is *not involved* in targeting proteins for degradation
- Flexible N-terminal region followed by a compact ball-shaped ubiquitin-like fold
- Nature designed SUMO to act as a chaperone for proteins
- SUMO improves solubility due to hydrophilic shell and hydrophobic core



# Benefits of the SUMO Platform

## PROBLEMS

Unstable protein,  
low expression

Specific N-terminus required

Insoluble protein

High cost of goods



## SOLUTIONS

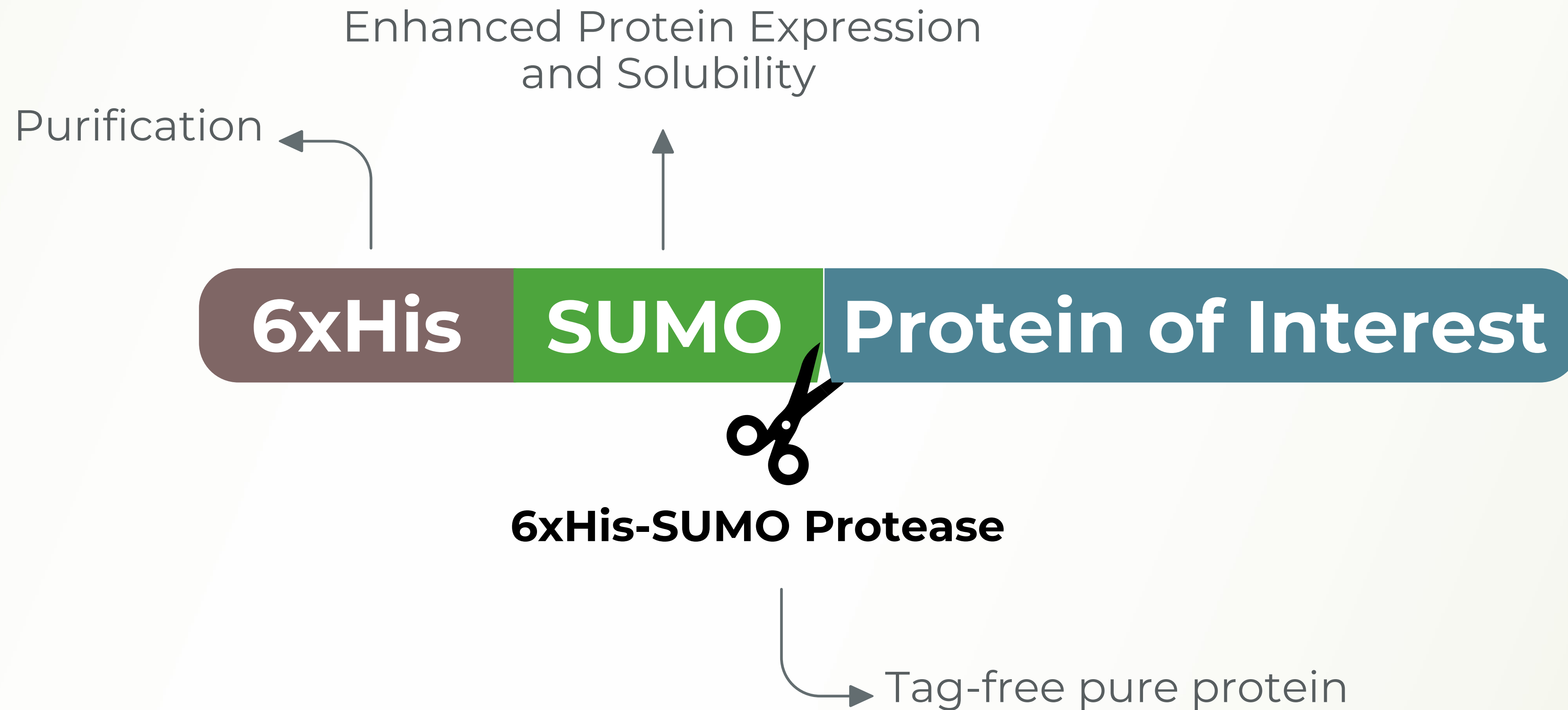
SUMO chaperoning,  
enhanced expression

SUMO-specific processing

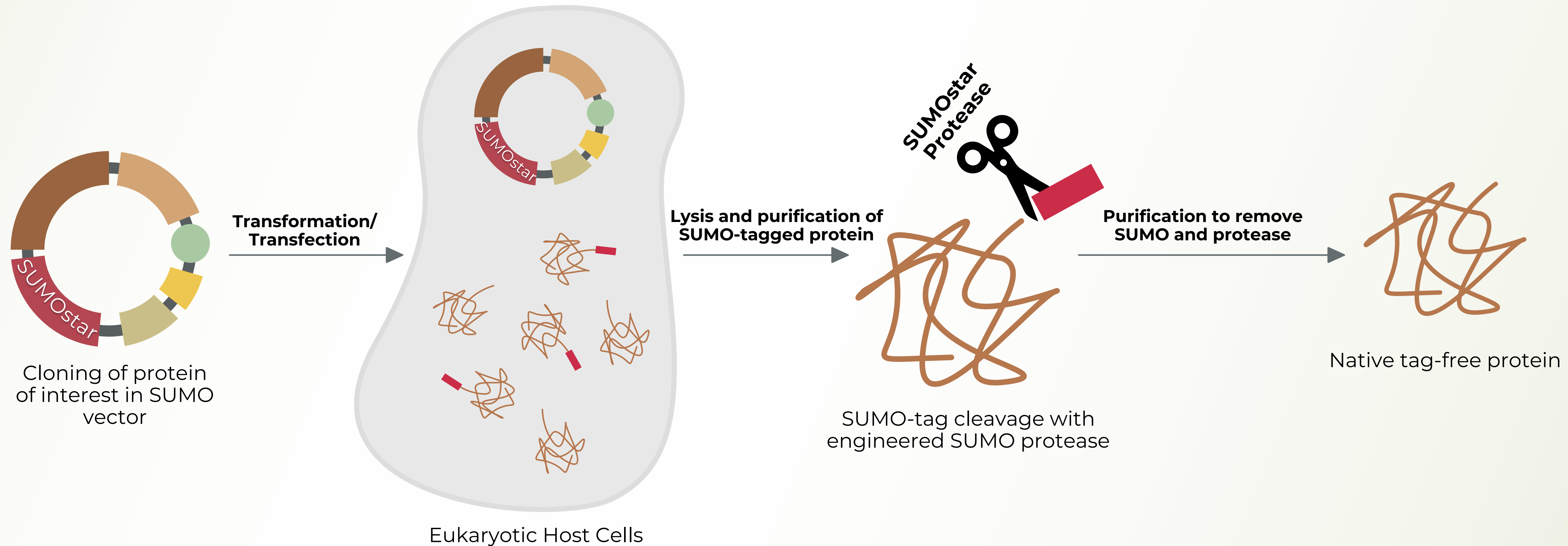
SUMO-driven folding

High yield, saves cost

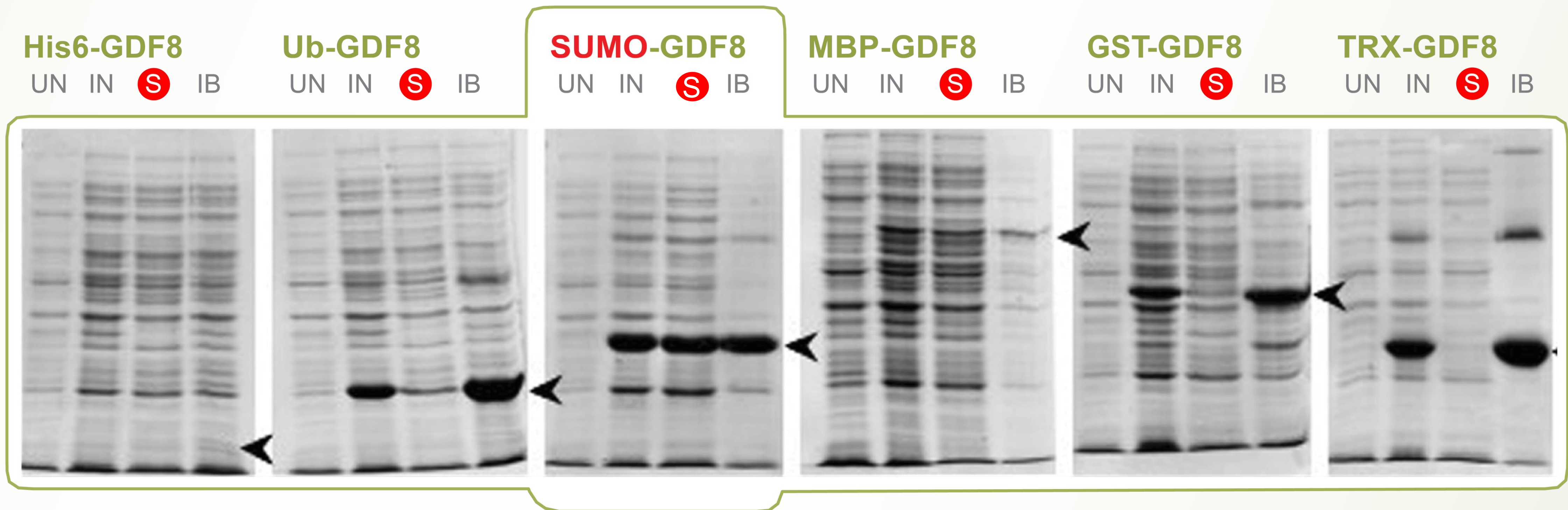
# SUMO Fusion Tag Platform For E.coli



# Protein Expression and Purification Process Using SUMO-tag



# SUMOpro<sup>®</sup> fusion enhances expression and solubility in *E. coli*



Conditions: UN – Uninduced, IN – Induced, S – Soluble Fraction, IB – Inclusion Bodies

Marblestone et. al, Protein Sci. 2006 Jan;15(1):182-9

# Therapeutic Cytokines/chemokines/GFs in the clinic

Therapeutic Protein/Polypeptide	Production Method	Company	Indication
<b>G-CSF (Filgrastim/Neupogen, Sargramostim (Leukine))</b>	Recombinant (E.coli, Yeast)	<a href="#">Amgen</a> , <a href="#">Sanofi</a>	low neutrophil count
<b>Erythropoietin (Epoetin alfa/Epogen, Procrit)</b>	Recombinant (E.coli, yeast)	<a href="#">Pfizer</a> , <a href="#">Amgen</a>	Chronic Renal Failure
<b>Interferon alfa-2b (Intron A)</b>	Recombinant (CHO), HEK293	<a href="#">Biogen (Avonex)</a> , <a href="#">EMD Serono/Merck (Rebif)</a> , <a href="#">Sigma</a>	Multiple Sclerosis
<b>RANTES (regulated on activation, normal T cell expressed and secreted)</b>	Recombinant (E.coli)	<a href="#">StemCell Tech</a>	<b>In development</b>
<b>IP-10 (interferon-inducible protein 10)</b>	Recombinant (E.coli)	<a href="#">Sigma</a>	<b>In development</b> (potent inhibitor of angiogenesis and displays thymus-dependent anti-tumor effects)
<b>SDF-1 (stromal cell-derived factor 1)</b>	Recombinant (E.coli)	<a href="#">Acro Biosystems</a>	<b>In development</b>
<b>Eotaxin</b>	Recombinant (HEK293)	<a href="#">Acro Biosystems</a>	<b>In development</b>
<b>MCP-1 (monocyte chemoattractant protein-1)</b>	Recombinant (E.coli)	<a href="#">R&amp;D Systems</a>	<b>In development</b>
<b>Platelet-derived growth factor (PDGF)</b>	Recombinant (E.coli, yeast)	<a href="#">BioLegend</a> , <a href="#">Regeneron (becaplermin)</a>	diabetic neuropathic ulcers
<b>Insulin-like growth factor 1 (IGF-1)</b>	Recombinant (E.coli, yeast)	<a href="#">Ipsen Biopharmaceuticals (Mecasermin/Increlex)</a>	growth failure and short stature in children with severe primary IGF-1 deficiency
<b>Epidermal growth factor (EGF)</b>	Recombinant (E.coli, HEK293)	<a href="#">BPS</a> , <a href="#">Abcam</a>	Skin care, In development for clinical use



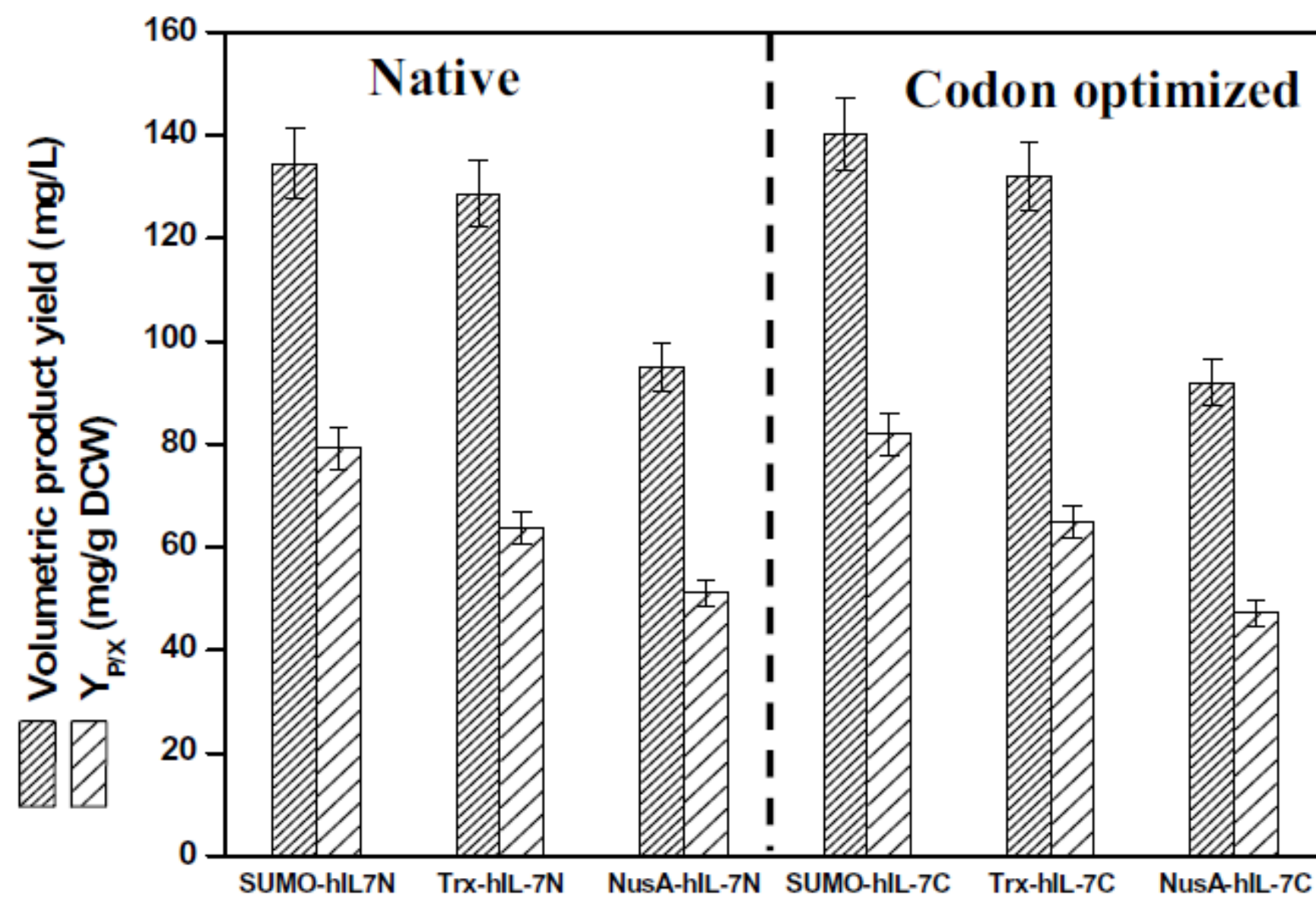
# Examples of SUMO-Enhanced Production of Therapeutic Proteins

	Peptide Name	Host	Yield (mg/L)
<b>Cytokines</b>	Human interleukin-7 (hIL-7) <sup>1</sup>	E.coli	135-140 mg/L
	TNF-alpha <sup>2</sup>	E.coli	97 mg/L
	IL-36 <sup>3</sup>	E.coli	10 mg/L
	IL-6 <sup>4</sup>	Nicotiana benthamiana leaves	18.5 mg/kg fresh weight
<b>Growth Factors</b>	Human Epidermal Growth Factor (hEGF) <sup>5,6</sup>	E.coli	281 mg/L
	Heparin-binding epidermal growth factor (HB-EGF) <sup>7,8</sup>	E.coli	104 mg/L
	FGF20 <sup>9</sup>	E.coli	-
	FGF21 <sup>10</sup>	E.coli	3.9 mg/L

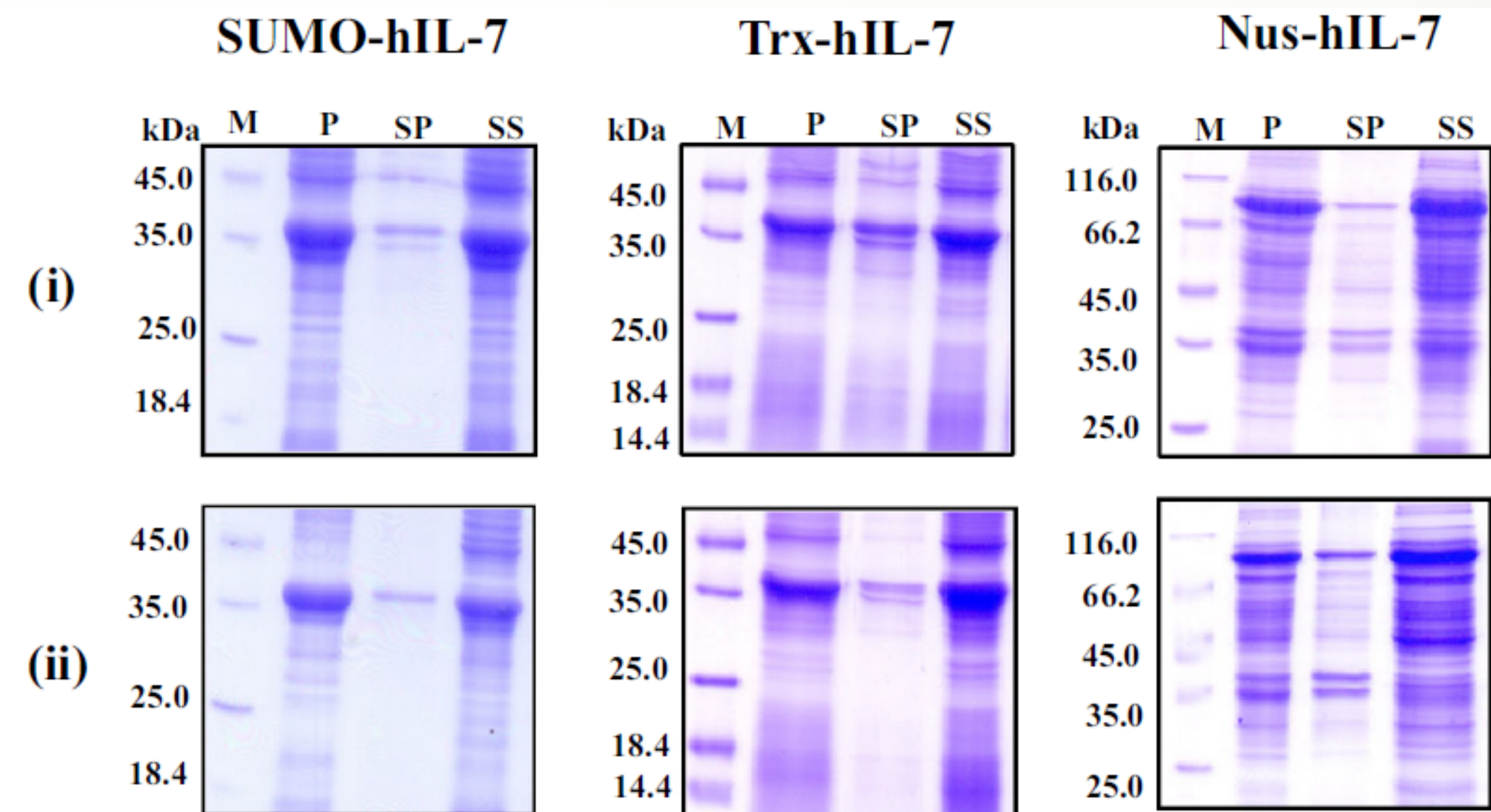
1. Devi N et al, Appl Microbiol Biotechnol, 2016
2. Hoffman A et al, Protein Expr Purif., 2010
3. Clancy DM et al, FEBS Open Bio. 2016
4. Islam MR et al, Plant Biotechnol J. 2019
5. Su Z et al, Protein Pept Lett. 2006;
6. Ma Y et al, Appl Microbiol Biotechnol, 2016
7. Lu, W. et al, Mol Biotechnol, 2010;
8. Ferreira AS et al, Sci Rep. 2022
9. Niu J et al Drug Deliv. 2018
10. Xu P et al, PLoS One. 2016

# Superior SUMO fusion enhanced the production of hIL-7 in E.coli

## Comparison Between SUMO, Trx and Nus Fusion



**Fig. 4** Shake flask expression studies of native and codon-optimized hIL-7 gene with different fusion tags (Trx, SUMO and NusA). **a, b** Growth profile ( $OD_{600}$ ) of native and codon-optimized hIL-7 gene,

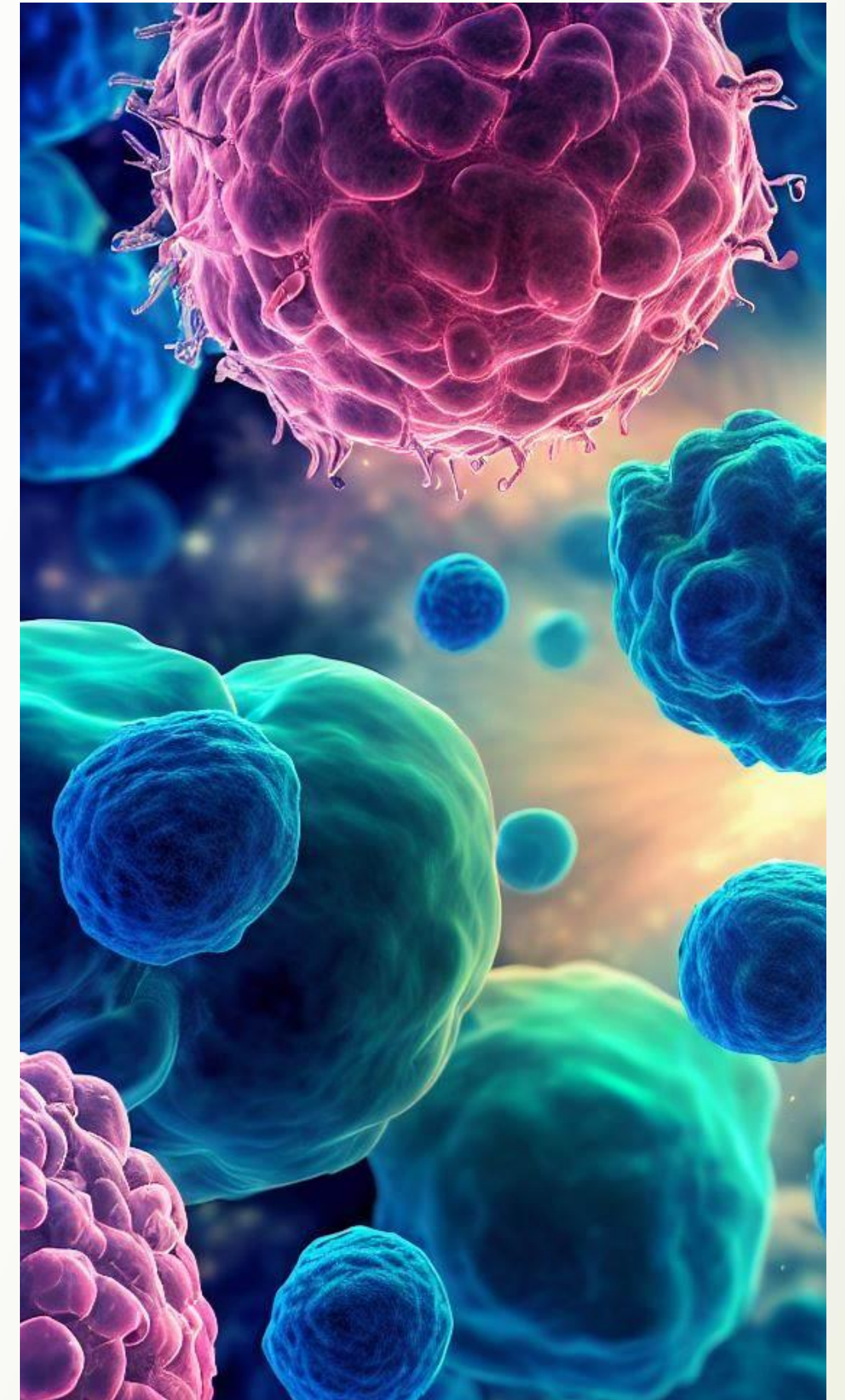


respectively. **c** Volumetric and specific product yield. **d** SDS-PAGE analysis (i) native and (ii) codon optimized

Devi, N., Adivitiya & Khasa, Y.P. *Appl Microbiol Biotechnol* **100**, 9979–9994 (2016).

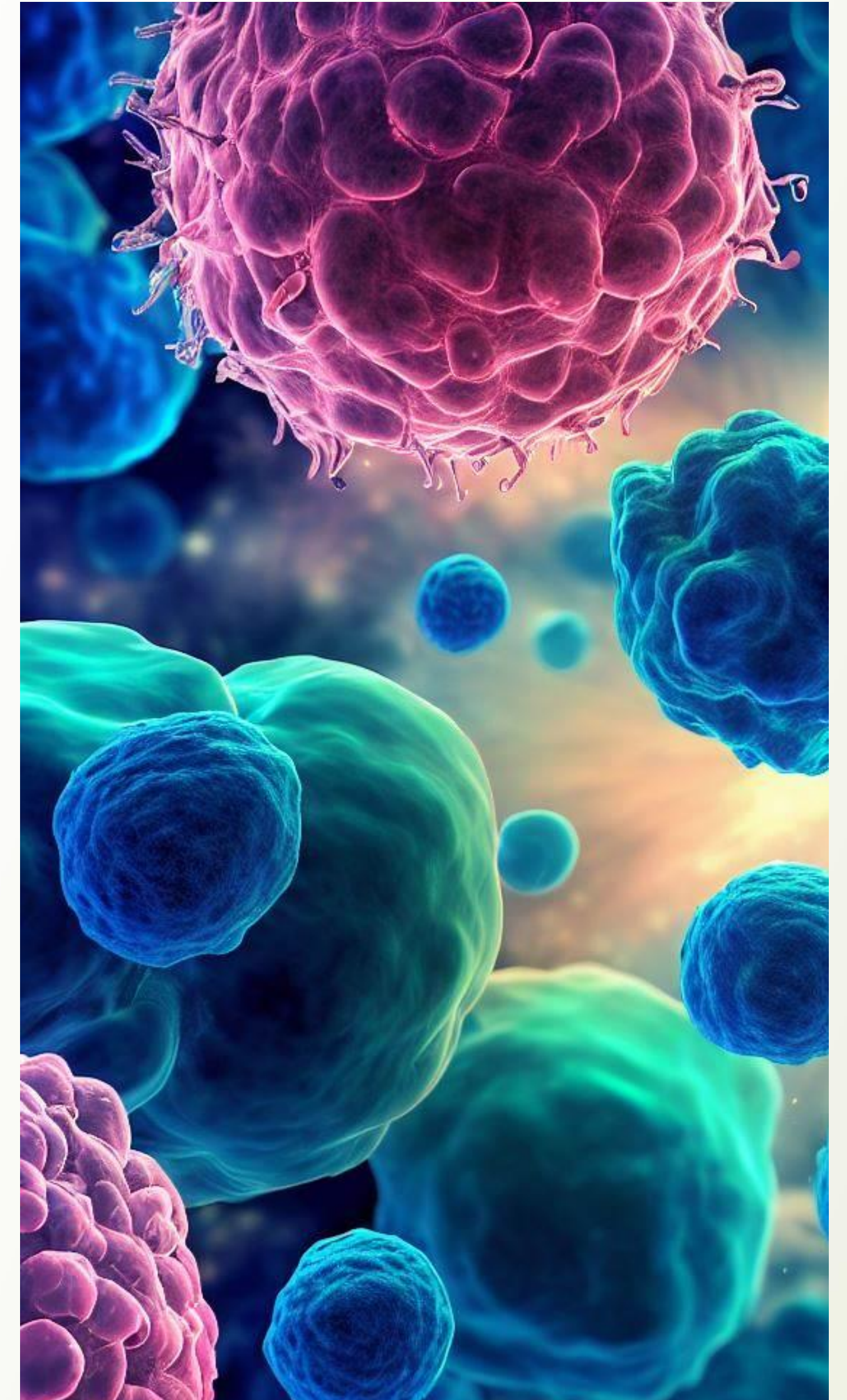
<https://doi.org/10.1007/s00253-016-7683-5>

# Challenges in expression of Cytokines, Chemokines and Growth Factors, Solutions offered by SUMO System



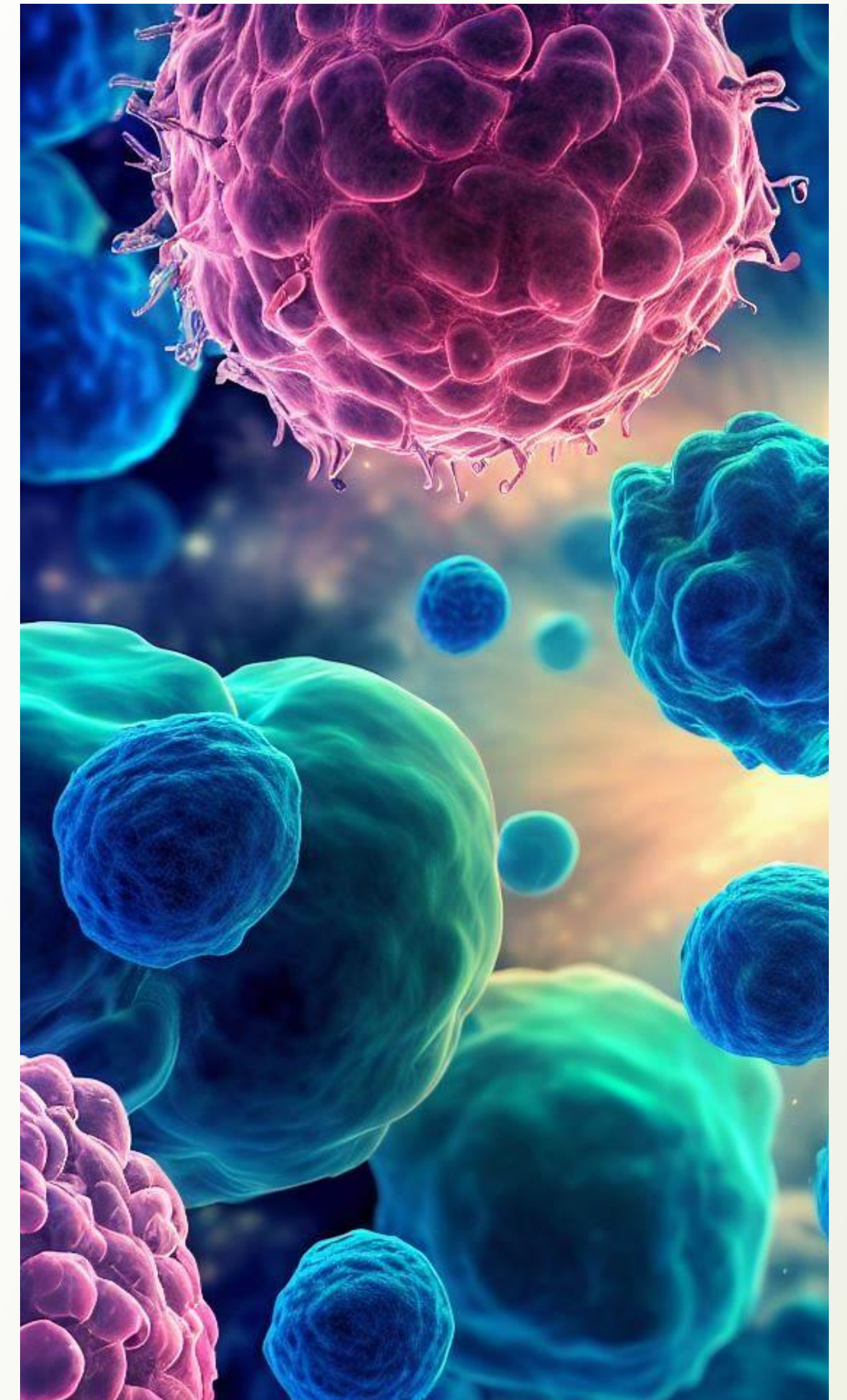
# Cytokines

- Families of cytokines expressed as SUMO-fusion, just bullets, no more than five bullets per slide.
- Need 1-2 real data slides, that will show impact of SUMO, no western blots or research centric data, just stained gel or manufacturing yield per liter
- Important references, 2-3, small font, at the end of the page, such as reviews can be listed



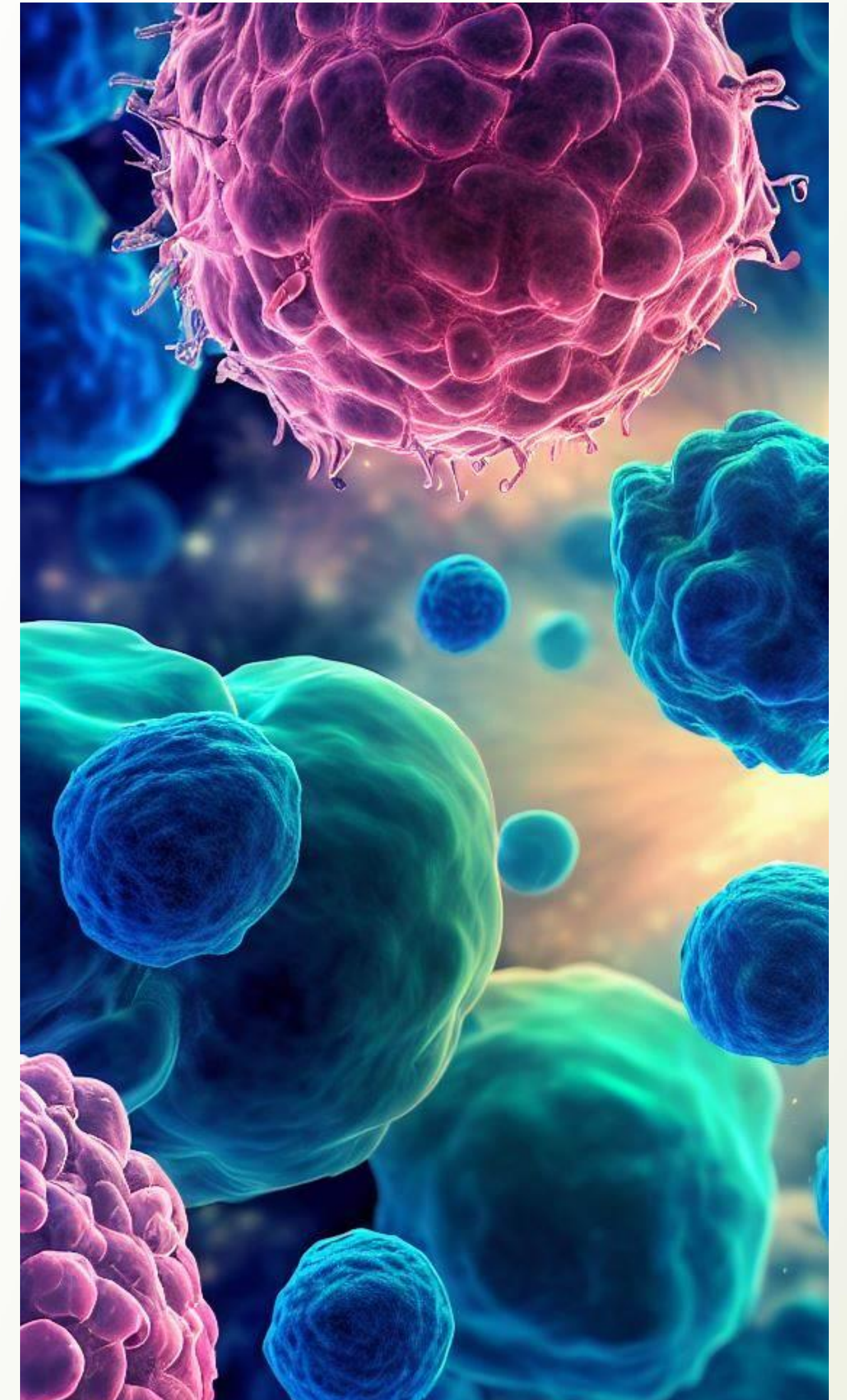
# Chemokines

- One line of intro re Chemokines
- Families of Chemokines expressed as SUMO-fusion, just bullets
- Need 1-2 real data slides, that will show impact of SUMO, no western blots or research centric data, just manufacturing
- Important references, 2-3, small font, at the end of the page, such as reviews can be listed



# Growth factors

- One line of intro re growth factor clinical value
- Families of growth factors expressed as SUMO-fusion, just bullets
- Need 1-2 real data slides, that will show impact of SUMO, no western blots or research centric data, just manufacturing
- Important references, 2-3, small font, at the end of the page, such as reviews can be listed



# Advantages of SUMO Mediated Enhanced Expression in Mammalian Cells

- Chemokines, cytokines and growth factors and other proteins are expressed as precursor proteins.
- SUMO facilitates expression of mature protein or peptide with desired N-terminus that is extremely important for biological activity
- Manufacturing proteins with stabilizing N-termini restores biological activity and improves yield
- Post-translational modifications are preserved with enhanced yield and biological activity
- Native SUMO is cleaved in mammalian cells, LifeSensors has developed an engineered SUMO that preserves SUMO chaperoning properties to enhanced expression while keeping the SUMO tag intact
- The engineered SUMO for mammalian cell is called SUMOstar. SUMOstar fusion can only be cleaved with SUMOstar protease

# SUMOstar, Preserving SUMO Enhanced Production in Eukaryotic Cells.

Generating Post-translationally Modified Proteins

Proteins that require post-translational modification, need to be produced in eukaryotic cells

## PROBLEM

Eukaryotic cells have endogenous SUMO proteases

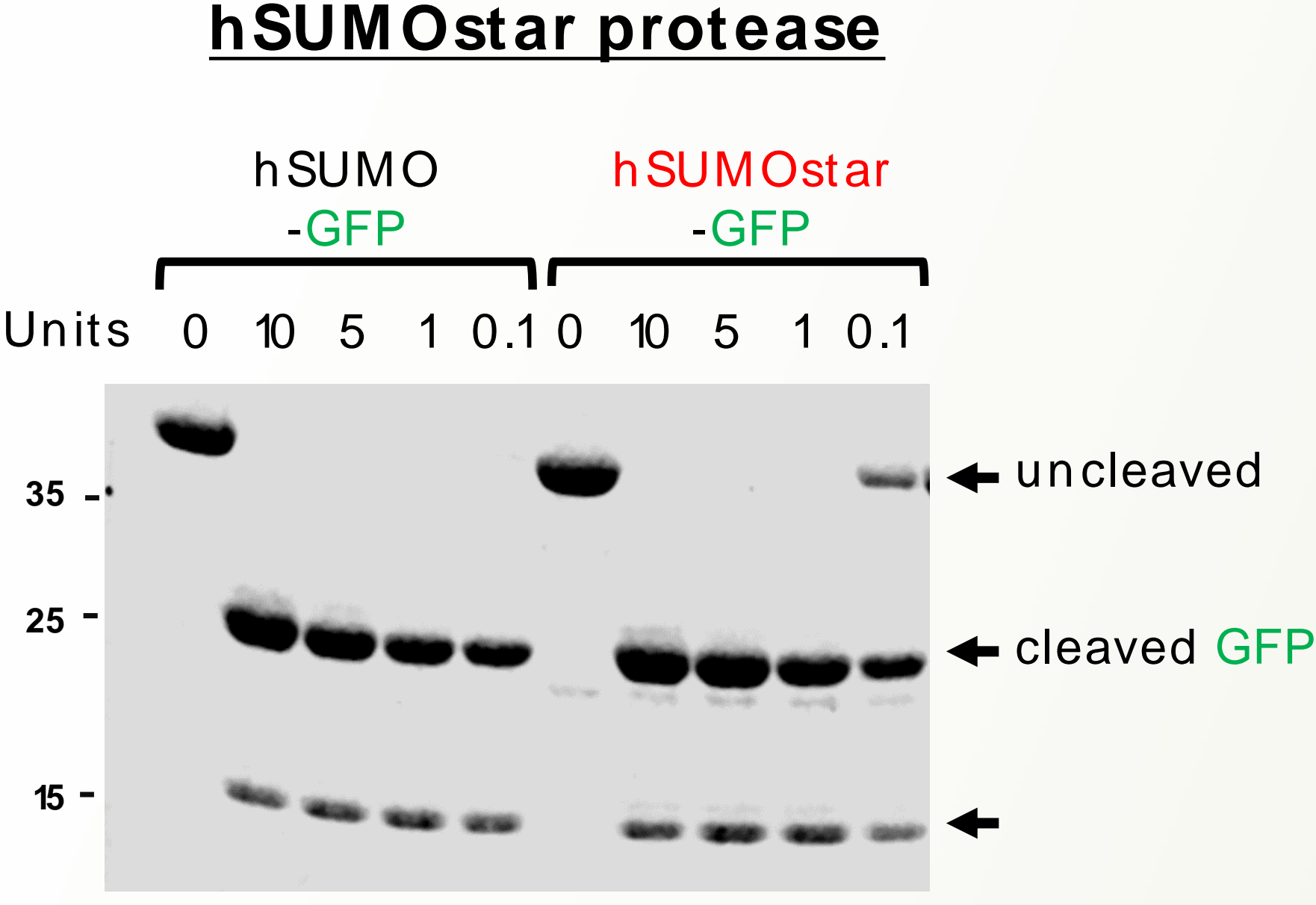
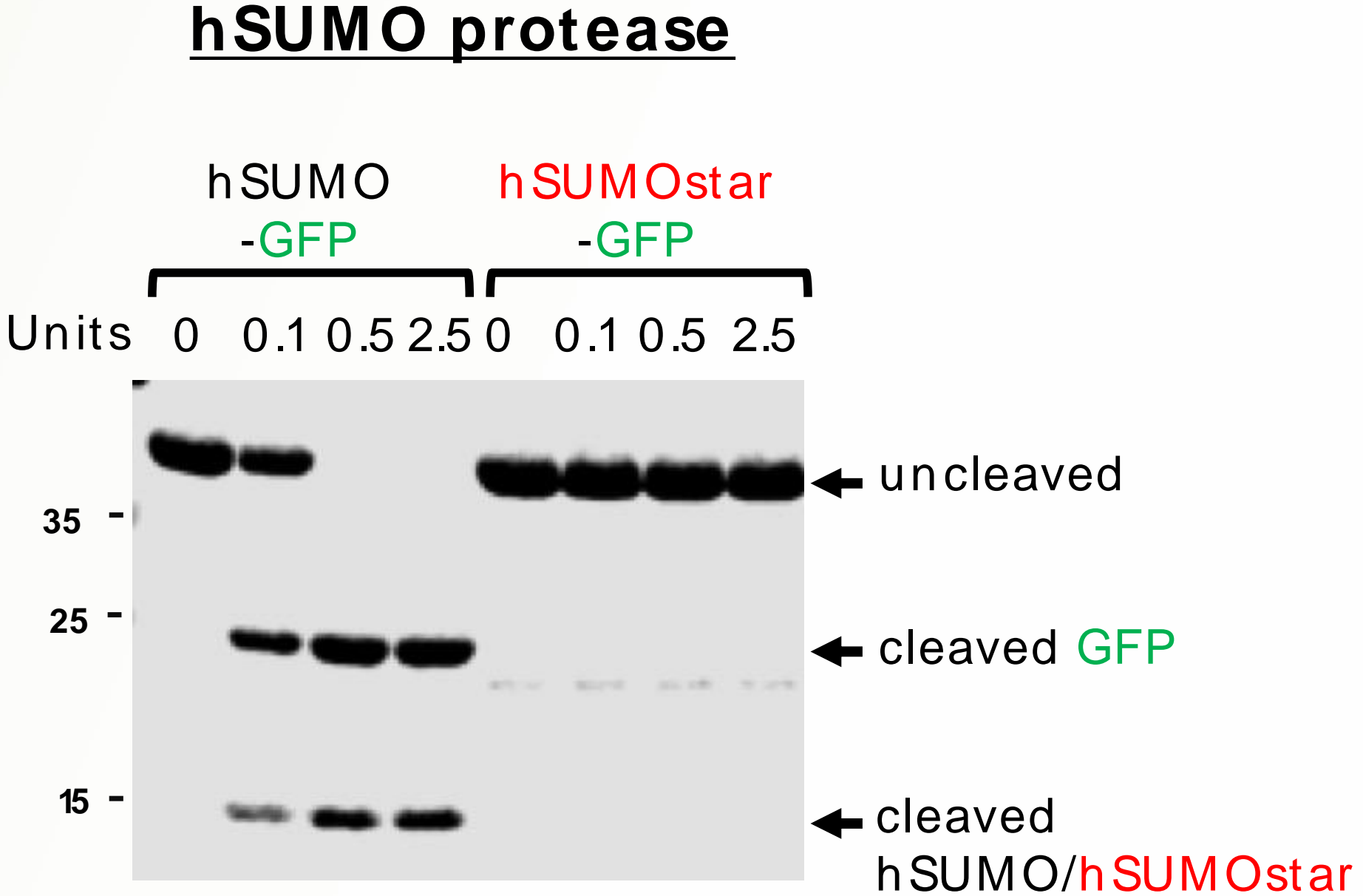
## SOLUTION

Engineered SUMOstar tag & SUMOstar protease

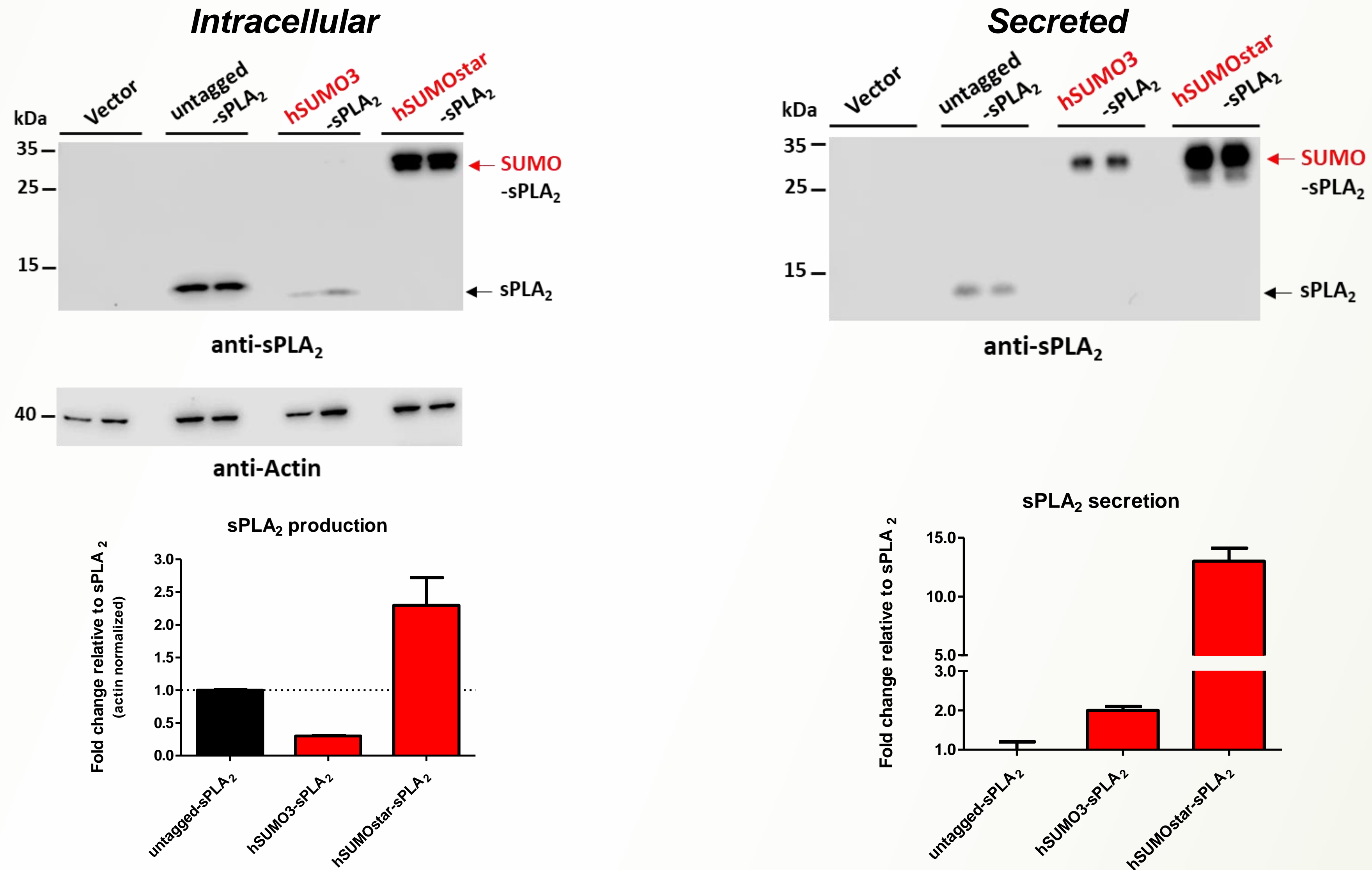




# hSUMOStar Protease Efficiently Cleaves the hSUMOStar Tag



# Human SUMOstar Dramatically Enhances sPLA<sub>2</sub> Production & Secretion in HEK293 cells



# Advantages of SUMO Mediated Enhanced Expression in Mammalian Cells, CAR-T and Gene Therapies

- Chimeric Antigen Receptors (CAR) are engineered genes that do not express well.
- Fusion with SUMO dramatically enhances their expression level and therapeutic function
- CAR system can be incorporated with human SUMO or hSUMOstar

# SUMOstar for Eukaryotic Expression

- Human SUMO and human SUMOstar dramatically enhances expression, and provides unique features not possible with traditional systems
- More than 150 publications describe therapeutic proteins expressed as SUMO or SUMOstar in E.coli and mammalian cells

# Thank You

We are your partner in Manufacturing Difficult to Express Proteins

## Contact Us!

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