



Eukaryotic Recombinant Protein Expression

LifeSensors' SUMO technologies are proven to dramatically enhance expression and solubility of heterologous proteins. Our SUMO proteases efficiently remove the SUMO fusion tags, leaving no residues behind giving you the unaltered protein you designed. The His₆-tags make purification fast and simple. Please contact us to see how our SUMO technology can benefit you.

E.coli



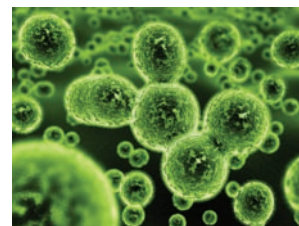
Mammalian



Insect



Yeast



Contact Us
To Learn More:
info@lifesensors.com

To order, please call
610-644-8845 or visit
www.lifesensors.com/sumostar



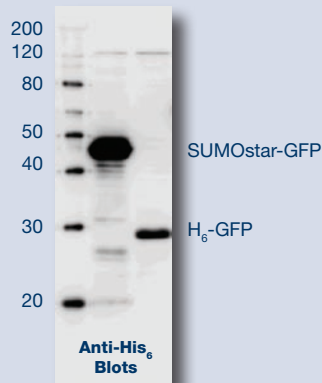
New Tag Technology for High-Level Expression of Recombinant Proteins in Eukaryotes

- Based on proven prokaryotic recombinant protein expression system
- Enhanced proper protein folding
- Improves solubility
- Cleavage yields native protein
- Rapid protein purification
- *E. coli*, Yeast, Insect and Mammalian cell expression systems

Express your protein in eukaryotic cells for proper post-translational modification! Fusion of heterologous proteins to the **S**mall **U**biquitin-like **M**odifier, SUMO, is a proven technology for enhancing the expression and solubility of proteins in *Escherichia coli*. Unlike native SUMO, the LifeSensors SUMOstar tag is not cleaved by endogenous desumoylases in eukaryotes improving folding, solubility, and yield of fused proteins. Our engineered SUMOstar desumoylase precisely and efficiently removes the SUMOstar tag following purification, generating the native N-terminus of the protein-of-interest.

Eukaryotic Expression

GFP expression in Insect Cells



PLA2 expression in Human HEK293

