



Buy One Try One!

Buy any one product, and receive any other product of equal or lesser value for free. August 16th - September 30th.

Color Key

SUMO PRO VECTORS

The original SUMOpro vectors offer increased expression and solubility for Prokaryotic systems. Benefits also include the use of highly specific SUMO Protease.

SUMO STAR VECTORS

A modified form of the original SUMOpro vectors that enables the use of SUMO for both Eukaryotic and Prokaryotic systems.

SUMO KITS

SUMO kits are a convenient package complete with SUMO Protease (1,000 units), Vector (Circular), Antibody and Control Protein.

SUMO Vectors and Systems

Catalog #	Product Name	Fusion Tag	Promoter	Selection Marker	System (Host)
E.coli					
1001K (Kan)	pE-SUMO Vector (E. coli)	His6-SUMO	T7	Kanamycin	E. coli
1001A (Amp)	pE-SUMO Vector (E. coli)	His6-SUMO	T7	Ampicillin	E. coli
1011 (Kan)	pE-SUMO3 Vector (E. coli)	His6-SUMO3	T7	Kanamycin	E. coli
1016 (Amp)	pE-SUMO3 Vector (E. coli)	His6-SUMO3	T7	Ampicillin	E. coli
LIC1000A	SUMOpro Protein and Peptide Expression Kit (Amp) LIC	His6-SUMO	T7	Ampicillin	E. coli
LIC1000K	SUMOpro Protein and Peptide Expression Kit (Kan) LIC	His6-SUMO	T7	Kanamycin	E. coli
1000A	SUMOpro Protein and Peptide Expression Kit (Amp)	His6-SUMO	T7	Ampicillin	E. coli
1000K	SUMOpro Protein and Peptide Expression Kit (Kan)	His6-SUMO	T7	Kanamycin	E. coli
1010A	SUMOpro-3 Protein and Peptide Expression Kit (Amp)	His6-HuSUMO3	T7	Ampicillin	E. coli
1010K	SUMOpro-3 Protein and Peptide Expression Kit (Kan)	His6-HuSUMO3	T7	Kanamycin	E. coli
1100A	SUMOstar Protein and Peptide Expression Kit (Amp)	His6-SUMOstar	T7	Ampicillin	E. coli
1100K	SUMOstar Protein and Peptide Expression Kit (Kan)	His6-SUMOstar	T7	Kanamycin	E. coli
LIC1001A	pE-SUMO LIC Vector (E. coli) Amp	His6-SUMO	T7	Ampicillin	E. coli
LIC1001K	pE-SUMO LIC Vector (E. coli) Kan	His6-SUMO	T7	Kanamycin	E. coli
1101 (Kan)	pE-SUMOstar Vector (E. coli)	His6-SUMOstar	T7	Kanamycin	E. coli
1106 (Amp)	pE-SUMOstar Vector (E. coli)	His6-SUMOstar	T7	Ampicillin	E. coli

Insect



LIC3100	Insect Intracellular SUMOstar Expression Kit LIC	His6-SUMOstar	Polyhedrin	Ampicillin	Insect
LIC3105	Insect Secretory SUMOstar Expression Kit LIC	gp67-His6-SUMOstar	Polyhedrin	Ampicillin	Insect
3100	Insect Intracellular SUMOstar Expression Kit	His6-SUMOstar	Polyhedrin	Ampicillin	Insect
3105	Insect Secretory SUMOstar Expression Kit	gp67-His6-SUMOstar	Polyhedrin	Ampicillin	Insect
LIC3106	pI-secSUMOstar LIC Vector (Insect-Secretory)	His6-SUMOstar	Polyhedrin	Ampicillin	Insect
LIC3101	pI-SUMOstar LIC Vector (Insect-Intracellular)	His6-SUMOstar	Polyhedrin	Ampicillin	Insect
3106	pI-secSUMOstar Vector (Insect-Secretory)	His6-SUMOstar	Polyhedrin	Ampicillin	Insect
3101	pI-SUMOstar Vector (Insect-Intracellular)	His6-SUMOstar	Polyhedrin	Ampicillin	Insect

SUMO Vectors and Systems

Catalog #	Product Name	Fusion Tag	Promoter	Selection Marker	System (Host)
Mammalian					
7110	Mammalian Intracellular SUMOstar Expression Kit	His6-SUMOstar	CMV	Ampicillin	Mammalian
7120	Mammalian Secretory SUMOstar Expression Kit	His6-secSUMOstar	CMV	Ampicillin	Mammalian
7121	pM-secSUMOstar (Mammalian-Secretory)	kappa secretory signal-His6-SUMOstar	CMV	Zeocin	Mammalian
7111	pM-SUMOstar (Mammalian-Intracellular)	His6-SUMOstar	CMV	Ampicillin (bacteria); Neomycin (mammalian)	Mammalian
Yeast					
2150	Pichia Secretory SUMOpro-3 Expression Kit	α -mating factor-His6-HuSUMO3	Alcohol Oxidase	Zeocin	Yeast (<i>P. pastoris</i>)
2151	pP-SUMO3 vector (<i>P.pastoris</i>)	α -mating factor-His6-huSUMO3	Alcohol Oxidase	Zeocin	Yeast (<i>P. pastoris</i>)
2101	pY-SUMOstar Vector (<i>S. cerevisiae</i>)	His6-SUMOstar	CUP	TRP	Yeast (<i>S. cerevisiae</i>)
2100	Saccharomyces Intracellular SUMOstar Expression Kit	His6-SUMOstar	CUP	TRP	Yeast (<i>S. cerevisiae</i>)

Protease, Protein, & Antibodies

PROTEASE

Catalog #	Item Name	Source	MW	Family	Buffer	Physical State	Purity	Alternate Name(s)	Fusion Tag	Concentration	Amount
4010	SUMO Protease 1	E.Coli	26.8 kDa	Cysteine Protease	Tris Buffer	Liquid	>95%	Ulp1 core	His6	10 U/ μ l	1000 units
4020	SUMO Protease 2	E.Coli	28.0 kDa	Cysteine Protease	Hepes Buffer	Liquid	>95%	SEN2 core	His6	10 U/ μ l	1000 units
4110	SUMOstar Protease 1	E.Coli	26.8 kDa	Cysteine Protease	Tris Buffer	Liquid	>95%		His6	10 U/ μ l	1000 units

SUMO PROTEINS

Catalog #	Item Name	Description
P6008	His6-huSUMO3	Recombinant His6-tagged Human SUMO3
P6004	His6-SUMO	Recombinant His6-tagged Yeast SUMO
P6012	His6-SUMOstar	Recombinant His6-tagged SUMOstar
P6006	huSUMO3 (untagged)	Recombinant untagged Human SUMO3
P6002	SUMO (Untagged)	Recombinant untagged Yeast SUMO
P6010	SUMOstar (untagged)	Recombinant untagged SUMOstar

ANTIBODIES

Catalog #	Item Name	Physical State	Specificity	Source Organism	Amount
AB7030	Anti-huSUMO3 antibody	Liquid	HuSUMO3	Chicken, affinity purified IgY	50 μ l
AB7002	Anti-SUMO/SUMOstar antibody	Liquid	SUMO (Smt3); SUMOstar	Chicken, affinity purified IgY	50 μ l

