



**pE-SUMO Vector (gp67, pPolh, Amp, 5761bp)**

6x His Tag Sdal

Met Gly His His His His His His His Gly Ser Leu Gln Asp  
 AATAATTTTGTTTAACTTTAAGAAGGAGATATACC **ATG** GGT CAT CAC CAT CAT CAT CAC GGG TCC CTG CAG GAC  
 TTATTTAAAACAAATTGAAATTCCTCTCTATATGG **TAC** CCA GTA GTG GTA GTA GTA GTA GTG CCC AGC GAC GTC CTG

Ser Glu Val Asn Gln Glu Ala Lys Pro Glu Val Lys Pro Glu Val Lys Pro Glu Thr His Ile  
 TCA GAA GTC AAT CAA GAA GCT AAG CCA GAG GTC AAG CCA GAA GTC AAG CCT GAG ACT CAC ATC  
 AGT CTT CAG TTA GTT CTT CGA TTC GGT CTC CAG TTC GGT CTT CAG TTC GGA CTC TGA GTG TAG  
**SUMOpro Fusion Tag**

Asn Leu Lys Val Ser Asp Gly Ser Ser Glu Ile Phe Phe Lys Ile Lys Lys Thr Thr Pro Leu  
 AAT TTA AAG GTG TCC GAT GGA TCT TCA GAG ATC TTC TTC AAG ATC AAA AAG ACC ACT CCT TTA  
 TTA AAT TTC CAC AGG CTA CCT AGA AGT CTC TAG AAG AAG TTC TAG TTT TTC TGG TGA GGA AAT

Arg Arg Leu Met Glu Ala Phe Ala Lys Arg Gln Gly Lys Glu Met Asp Ser Leu Arg Phe Leu  
 AGA AGG CTG ATG GAA GCG TTC GCT AAA AGA CAG GGT AAG GAA ATG GAC TCC TTA AGA TTC TTG  
 TCT TCC GAC TAC CTT CGC AAG CGA TTT TCT GTC CCA TTC CTT TAC CTG AGG AAT TGC AAG AAC

Tyr Asp Gly Ile Arg Ile Gln Ala Asp Gln Ala Pro Glu Asp Leu Asp Met Glu Asp Asn Asp  
 TAC GAC GGT ATT AGA ATT CAA GCT GAT CAG GCC CCT GAA GAT TTG GAC ATG GAG GAT AAC GAT  
 ATG CTG CCA TAA CTT TAA GTT CGA CTA GTC CGG GGA CTT CTA AAC CTG TAC CTC CTA TTG CTA

Ile Ile Glu Ala His Arg Glu Gln Ile Gly Gly BsaI  
 ATT ATT GAG GCT CAC CGC GAA CAG ATT **GG** AGGTTGAGACC ACTAGTGGTACC GGTCTCA **CTAGA**  
 TAA TAA CTC CGA GTG GCG CTT GTC TAA **CCTCCA** ACTCTGG TGATCACCATGG CCAGAGTGATC **T**  
**SUMO Protease 1 Cleavage Site** **BsaI** **MCS**

**Multiple Cloning Site (MCS)**

pE-SUMO vector has to be linearized with BsaI or Eco31I restriction enzyme

Gly Gly NotI

**GG** XbaI<sup>1</sup> BamHI SacI Sall HindIII EagI XhoI  
**CCTCCA** **CTAGA** GGATCC GAATTC GAGCTCC GTCGAC AAGCTT GCGGCCG ACTCGAG CACCA  
 T CCTAGG CTTAAG CTCGAGG CAGCTG TTCGAA CGCCGGCG TGAGCTC GTGGT

