

Polylinker Map (pP- α SUMO3 Vector, AOX1, Zeo)
Vector Size: 3840bp



Met Arg Phe Pro Ser Ile Phe Thr Ala Val Leu Phe Ala Ala Ser Ser Ala Leu
 ACTAATTATTCGAAACG **ATG** AGA TTT CCT TCA ATT TTT ACT GCT GTT TTA TTC GCA GCA TCC TCC GCA TTA
 TGATTAATAAGCTTTGC TAC TCT AAA GGA AGT TAA AAA TGA CGA CAA AAT AAG CGT CGT AGG AGG CGT AAT

Alpha Secretion signal

Ala Ala Pro Val Asn Thr Thr Thr Glu Asp Glu Thr Ala Gln Ile Pro Ala Glu Ala Val Ile Gly
 GCT GCT CCA GTC AAC ACT ACA ACA GAA GAT GAA ACG GCA CAA ATT CCG GCT GAA GCT GTC ATC GGT
 CGA CGA GGT CAG TTG TGA TGT TGT CTT CTA CTT TGC CGT GTT TAA GGC CGA CTT CGA CAG TAG CCA

Tyr Ser Asp Leu Glu Gly Asp Phe Asp Val Ala Val Leu Pro Phe Ser Asn Ser Thr Asn Asn Gly
 TAC TCA GAT TTA GAA GGG GAT TTC GAT GTT GCT GTT TTG CCA TTT TCC AAC AGC ACA AAT AAC GGG
 ATG AGT CTA AAT CTT CCC CTA AAG CTA CAA CGA CAA AAC GGT AAA AGG TTG TCG TCT TTA TTG CCC

Leu Leu Phe Ile Asn Thr Thr Ile Ala Ser Ile Ala Ala Lys Glu Glu Gly Val Ser Leu Glu Lys
 TTA TTG TTT ATA AAT ACT ACT ATT GCC AGC ATT GCT GCT AAA GAA GAA GGG GTA TCT CTC GAG AAA
 AAT AAC AAA TAT TTA TGA TCA TAA CGG TCG TAA CGA CGA TTT CTT CTT CCC CAT AGA GAG CTC TTT

6x His Tag

SdaI

Arg Gly His His His His His His Gly Ser Leu Gln Glu Glu Lys Pro Lys Glu Gly Val Lys Thr
 AGA GGT CAT CAC CAT CAT CAT CAC GGG TCC CTG CAG GAG GAG AAG CCC AAG GAG GGT GTG AAG ACA
 TCT CCA GTA GTG GTA GTA GTA GTG CCC AGG GAC GTC CTC CTC TTC GGG TTC CTC CCA CAC TTC TGT

Human SUMO3 Fusion Protein

Glu Asn Asp His Ile Asn Leu Lys Val Ala Gly Gln Asp Gly Ser Val Val Gln Phe Lys Ile Lys
 GAG AAT GAC CAC ATC AAC CTG AAG GTG GCC GGG CAG GAC GGC TCC GTG GTG CAG TTC AAG ATC AAG
 CTC TTA CTG GTG TAG TTG GAC TTC CAC CGG CCC GTC CTG CCG AGG CAC CAC GTC AAG TTC TAG TTC

Arg His Thr Pro Leu Ser Lys Leu Met Lys Ala Tyr Cys Glu Arg Gln Gly Leu Ser Met Arg Gln
 AGG CAC ACG CCG CTG AGC AAG CTG ATG AAG GCC TAC TGC GAG AGG CAG GGC TTG TCA ATG AGG CAG
 TCC GTG TGC GGC GAC TCG TTC GAC TAC TTC CGG ATG ACG CTC TCC GTC CCG AAC AGT TAC TCC GTC

Ile Arg Phe Arg Phe Asp Gly Gln Pro Ile Asn Glu Thr Asp Thr Pro Ala Gln Leu Glu Met Glu
 ATC AGA TTC AGG TTC GAC GGG CAG CCA ATC AAT GAA ACT GAC ACT CCA GCA CAG CTG GAG ATG GAG
 TAG TCT AAG TCC AAG CTG CCC GTC GGT TAG TTA CTT TGA CTG TGA GGT CGT GTC GAC CTC TAC CTC

Asp Glu Asp Thr Ile Asp Val Phe Gln Gln Gln Thr Gly Gly **BsmBI**
 GAC GAG GAC ACC ATC GAC GTG TTC CAG CAG CAG ACG **GG** AGGTAGAGAC GATCTGC CGTCTCA **CTAGA**
 CTG CTC CTG TGG TAG CTG CAC AAG GTC GTC GTC TGC **CCTCCA** TCTCTG CTAGACG GCAGAGTGATC **T**

SUMO Protease 2 cleavage site

BsmBI

Multiple Cloning Site (MCS)

pP- α SUMO3 vector has to be linearized with BsmBI restriction enzyme

