

GP78

Cat. # UB310

Background: GP78 is a ligase that functions in endoplasmic reticulum-associated degradation (ERAD). This enzyme construct is composed of the GP78 cytosolic domain (residues 309-643) with an N-terminal GST tag.

Application: Useful in conjunction with E1 and E2 for the study of enzyme kinetics, screening inhibitors and selectivity profiling.

Product Information

Molecular Weight:	72 kDa
Purity:	>80% by SDS-PAGE
Quantity:	20 µg
Physical State:	Liquid, 45 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 10% glycerol, 3 mM DTT
Species:	Human
Host:	E. coli
Tag:	N-terminal GST tag
Storage/Stability:	>6 months at -80°C

References

1. Fang S. *et al.*, *Proc. Natl. Acad. Sci. U.S.A.* **98** (25),14422-14427 (2001)
2. Zhang T. *et al.*, *PLoS One* **10** (3), e0118448 (2015)
3. Joshi V. *et al.* *Front. Cell Neurosci.* **11**, 259 (2017)

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2014 LifeSensors, Inc. All Rights Reserved