

**Background**

GP78 is an E3 ubiquitin ligase involved in endoplasmic reticulum-associated degradation (ERAD). This construct consists of the cytosolic domain of GP78 (residues 309–643) fused to an N-terminal GST tag.

**Alternate Names**

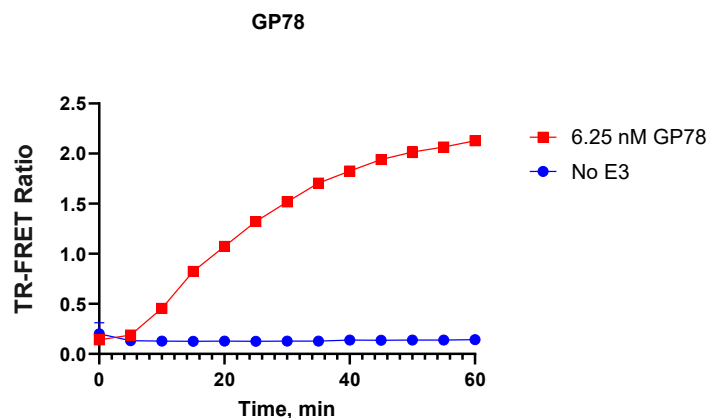
Autocrine Motility Factor Receptor (AMFR), E3 Ubiquitin-Protein Ligase AMFR, RING Finger Protein 45 (RNF45), RING-Type E3 Ubiquitin Transferase AMFR, AMF Receptor, SPG89

**Application(s)**

Ubiquitin ligation reactions

**Product Specifications**

<b>Affinity tag</b>	N-terminal GST tag
<b>Purity</b>	> 80% by SDS-PAGE
<b>Molecular Weight</b>	72 kDa
<b>Quantity</b>	25 µg
<b>Species</b>	Human
<b>Expression System</b>	<i>E. coli</i>
<b>Physical State</b>	Liquid
<b>Buffer</b>	45 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 10% glycerol, 3 mM DTT
<b>Activity</b>	A typical enzyme concentration of 2-100 nM is used for in vitro conjugation, depending on assay conditions
<b>Stability &amp; Storage</b>	> 6 months at -80°C. Avoid repeated freeze/thaw cycles

**Product QC**

**Activity Assay of GP78.** GP78 at a concentration of 6.25 nM was tested in a TR-FRET assay for 60 minutes, yielding a signal-to-background ratio of 15-fold.

**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)