

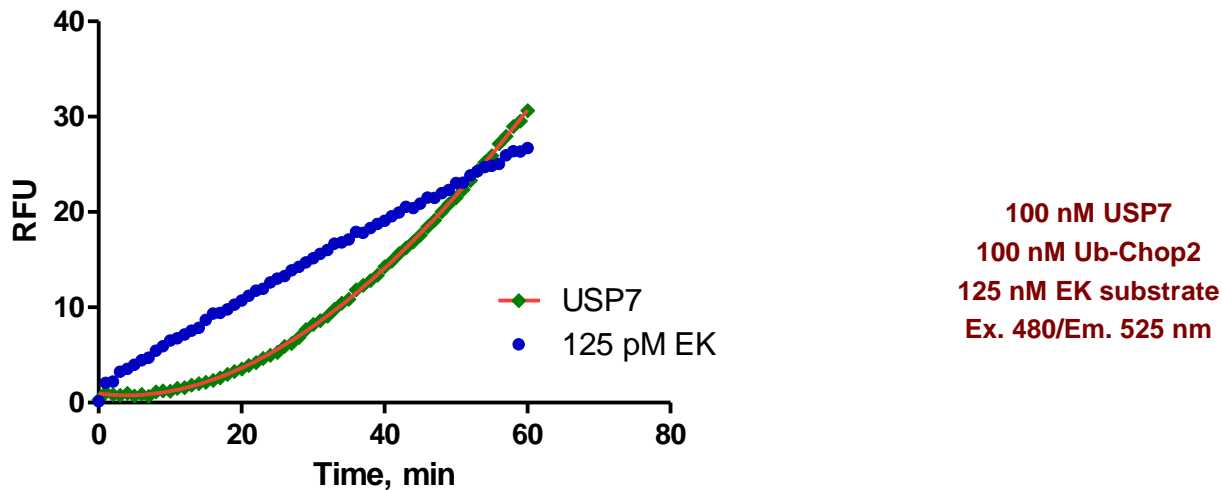
Enterokinase catalytic domain

Cat. # PR11EK

Background: Enterokinase is a serine protease with specificity for cleavage after the lysine residue in the sequence Asp-Asp-Asp-Asp-Lys↓. LifeSensors enterokinase CD is produced recombinantly in *E. coli*, purified, and refolded for use as a counterscreen with the UBL-CHOP2 assay suite (Nicholson, Leach et al. 2008).

Application: Cleavage after a FLAG™ sequence; Counterscreen in CHOP2 assay

Product Information	EC 3.4.21.9
Accession No:	P98073
Residues:	785 to 1019
Purity:	N/A
Molecular Weight:	26,246.8 Da
Physical State:	liquid
Storage:	-80° C. Avoid repeated freeze/thaw cycles



References

Nicholson, B., C. A. Leach, et al. (2008). "Characterization of ubiquitin and ubiquitin-like protein isopeptidase activities." *J Prot. Sci.* **17**: 1035-1043.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2015 LifeSensors, Inc. All Rights Reserved