

Human recombinant ubiquitin K33

Cat. # SI206

Background: Human recombinant ubiquitin containing a single lysine at position 33 is produced in *E. coli* and purified to >95% homogeneity. This protein contains no extraneous tags.

Application: For use in conjunction with the E3LITE Customizable Ligase Assay and E2 Profiling Kit to investigate conjugation/ligation enzyme activity and/or drug discovery. Can also be used for conventional ubiquitin pathway research.

Product Information

Purity:	≥ 95% by SDS-PAGE
Molecular Weight:	8,732.9 Da
Physical State:	Liquid (PBS)
Quantity:	1mg at 4mg/mL
Solubility:	>30mg/ml
Storage:	-80° C. Avoid repeated freeze/thaw cycles

References

1. Wilkinson, K. D. (2000). "Ubiquitination and deubiquitination: targeting of proteins for degradation by the proteasome." *Semin Cell Dev Biol* 11(3): 141-8.
2. Ciechanover, A. (2003) "The ubiquitin proteolytic system and pathogenesis of human diseases: a novel platform for mechanism-based drug targeting." *Biochem Soc Trans* 31(2): 474-81.
3. Marblestone, J.G., Kumar, K.G., Eddins, M.J., Leach C.A., Sterner, D.E., Mattern, M.R., Nicholson, B. (2010) Novel Approach for Characterizing Ubiquitin E3 Ligase Function. *J Biomol Screen* (In Press).

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2009 LifeSensors, Inc. All Rights Reserved