

Parkin C457S Mutant (His₁₀-SUMO)

Cat. # UB323

Background: Encoded by the PARK2 gene, the E3 ligase Parkin is part of the multi-protein E3 complex that encodes substrate proteins for degradation in the Ubiquitin-Proteasome Pathway¹. The precise function of this protein is unknown; however, mutations in this gene are known to cause a familial form of Parkinson's disease known as autosomal recessive juvenile Parkinson's disease (AR-JP). Parkin is described to be necessary for mitophagy (autophagy of mitochondria). The C457S mutant is a proposed activating mutant².

Application: For use in investigating and research of the Parkin and PINK1 pathway and/or drug discovery.

Product Information

Purity:	≥ 95% by SDS-PAGE
Molecular Weight:	64 kDa
Physical State:	Liquid, 50 mM Tris, pH 7.5, 0.15 M NaCl, 10% Glycerol
Species:	Human
Source:	<i>E. coli</i>
Quantity:	25 µg
Tag:	His ₁₀ -SUMO
Storage:	-80° C. Avoid repeated freeze/thaw cycles

References

1. Wang, X.-L. *et al.* Parkin, an E3 Ubiquitin Ligase, Plays an Essential Role in Mitochondrial Quality Control in Parkinson's Disease. *Cell Mol Neurobiol* **41**, 1395–1411 (2021).
2. Ordureau, A. *et al.* Quantitative Proteomics Reveal a Feedforward Mechanism for Mitochondrial PARKIN Translocation and Ubiquitin Chain Synthesis. *Molecular Cell* **56**, 360–375 (2014).

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