

Parkin Δ UBL (His₁₀-SUMO)

Cat. # UB319

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 Pathway. 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). Phosphorylation of Parkin's UBL at Ser65 by PINK1 kinase activates Parkin. The UBL domain has been removed from this from this construct.

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

Product Information

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

References

1. Sauv , V. *et al.* A Ubl/ubiquitin switch in the activation of Parkin. *EMBO J* **34**, 2492–2505 (2015).
 2. Quinn, P. M. J., Moreira, P. I., Ambr sio, A. F. & Alves, C. H. PINK1/PARKIN signalling in neurodegeneration and neuroinflammation. *Acta Neuropathologica Communications* **8**, 189 (2020).
-

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