

Untagged Parkin E3 Ligase

Cat. # UB318

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).

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

Product Information

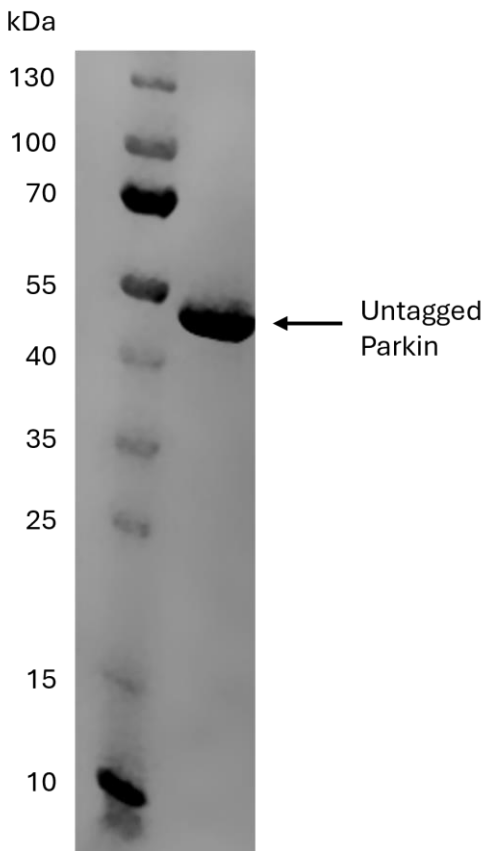
Species:	Human
Source:	<i>E. coli</i>
Purity:	≥ 95% by SDS-PAGE
Molecular Weight:	51.6 kDa
Tag:	None (Untagged)
Quantity:	25 µg
Physical State:	Liquid, 50 mM Tris, pH 7.5, 0.15 M NaCl, 10% Glycerol
Quantity:	25 µg
Storage:	-80° C. Avoid repeated freeze/thaw cycles

References

1. Ge, P., Dawson, V. L. & Dawson, T. M. PINK1 and Parkin mitochondrial quality control: a source of regional vulnerability in Parkinson's disease. *Molecular Neurodegeneration* **15**, 20 (2020).
 2. Quinn, P. M. J., Moreira, P. I., Ambrósio, 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

Data



SDS-Page Analysis of purified Untagged Parkin. Four μg of the protein was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue

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