

Parkin (human) polyclonal antibody

Parkin is a ubiquitin-protein ligase (E3), and mutations in the Parkin gene appear to be closely associated with autosomal recessive juvenile Parkinson's disease. Parkin has been localised to Lewy bodies in patients with sporadic Parkinson's disease, and in peri-nuclear aggresome-like inclusions in cells over-expressing wild-type Parkin where proteasomal activity has also been inhibited⁴. Both types of inclusion also contain ubiquitin, α -synuclein, proteasomal subunits and heat shock proteins. A recent paper suggests that loss of proteasomal activity leads to Parkin accumulation with a concomitant reduction in E3 ligase activity resulting in Lewy body formation⁵. Parkin is selectively recruited from the cytosol to damaged mitochondria to trigger their autophagy (mitophagy).

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-PW9365-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application IHC, IP, WB

Formulation Liquid. In PBS.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 399-415 (CE339QARWEAASKETIK412) of human parkin. A single amino acid substitution (A405E) exists in both the rat and mouse homologues.

Species Reactivity Human

UniProt ID O60260

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#)

.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)