

Parkin [pS101], Phosphospecific Antibody

Catalog no. 480026

(See product label for lot information)

Product Description

Affinity purified rabbit polyclonal antibody

Clone/PAD: N/A
Isotype: N/A
Qty: 100µl

Formulation

10 mM HEPES (pH 7.5), 150 mM NaCl,
100 µg per ml BSA and 50% glycerol.

Purification Method:

Prepared from rabbit serum by affinity purification via sequential chromatography on phospho- and dephosphopeptide affinity columns.

Validation

See www.invitrogen.com/antibodies for protocols

Validated for use in WB.

WB: 1:1000

Reactivity

This product had been directly tested for reactivity with Human.

Immunogen

Phosphopeptide corresponding to amino acid residues surrounding the phospho-Ser101 of human parkin.

Storage

2-8°C for up to 1mo, -20°C for long term storage. Avoid repeated freezing and thawing.

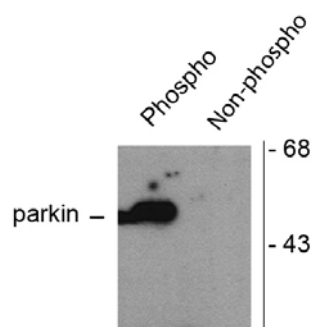
Expiration Date

Expires one year from date of receipt when stored as instructed.



Background

Parkin is an E3 ligase in the ubiquitin-proteasome system. Hereditary Parkinson's disease is most commonly caused by mutations in the parkin gene and is characterized by the progressive loss of dopaminergic neurons and the presence of Lewy bodies in the substantia nigra (Jenner et al., 1992). Recent evidence suggests that phosphorylation of parkin at Ser101 may have an important regulatory role on its E3 ubiquitin ligase activity (Yamamoto et al., 2005).



Western blot of HEK293 cells transfected with Parkin WT (Phospho) and Parkin S101 mutant (non-phospho) showing the phospho-specific immunolabeling of the ~ 52 k parkin protein. The immunolabeling is absent in the parkin S101 mutant.

References

- Jenner P, Dexter DT, Sian J, Schapira AH, Marsden CD (1992) Oxidative stress as a cause of nigral cell death in Parkinson's disease and incidental Lewy body disease. *Ann Neurol.* 32 Suppl: S82-7.
- Yamamoto A, Friedlein A, Imai Y, Takahashi R, Kahle PJ, Haass C (2005) Parkin phosphorylation and modulation of its E3 ubiquitin ligase activity. *J Biol Chem.* 280(5):3390-9

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com