

# PARK7 recombinant monoclonal antibody (007)

PARK7 (DJ-1) is a protein with antioxidant properties that plays a protective role against oxidative stress and cell death. It is highly relevant in neurodegenerative disease research, particularly in Parkinson's disease, where mutations in PARK7 are linked to early-onset forms of the condition.

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

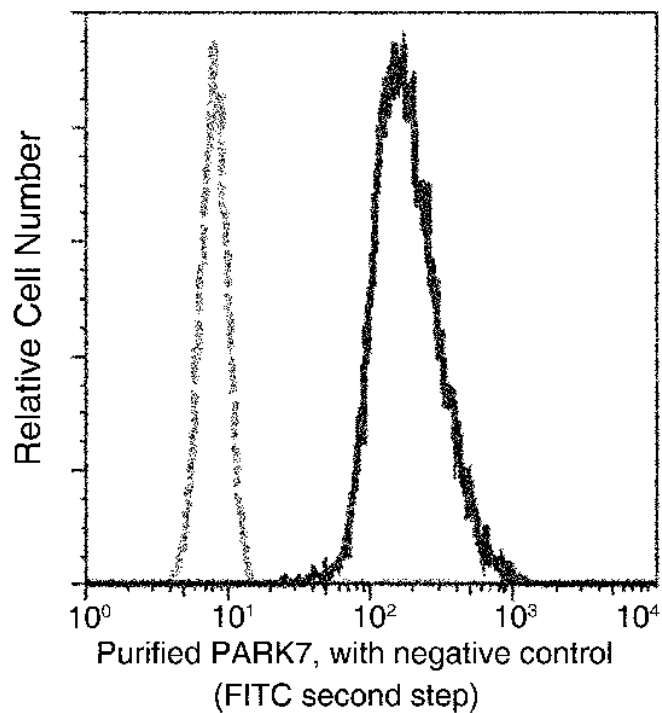
## Ordering Information

[Order Online »](#)

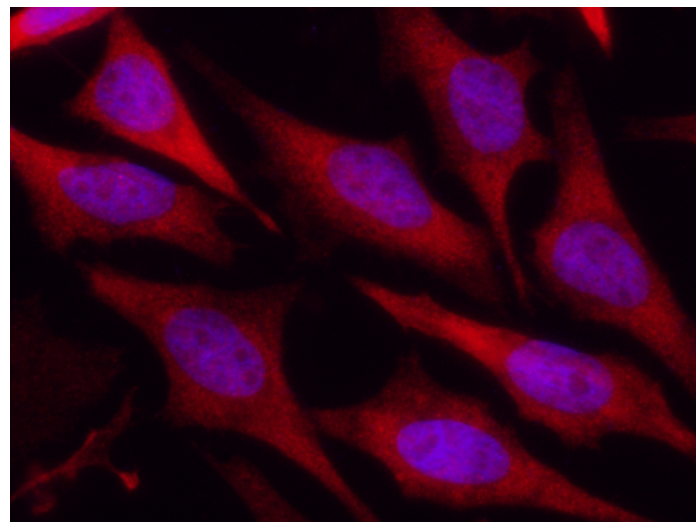
ENZ-ABS1038-0100	100µl
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Human PARK7 expression on HeLa cells. The cells were treated according to manufacturer's manual (BD Pharmingen™ Cat. No. 554714), stained with purified anti-Human PARK7, then a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.



Immunofluorescence staining of Human PARK7 in HeLa cells. Cells were fixed with 4% PFA, permeabilized with 0.3% Triton X-100 in PBS, blocked with 10% serum, and incubated with rabbit anti-Human PARK7 monoclonal antibody (1:60) at 37°C 1 hour. Then cells were stained with the Alexa Fluor® 594-conjugated Goat Anti-rabbit IgG secondary antibody (red) and counterstained with DAPI (blue). Positive staining was localized to cytoplasm.

## Handling & Storage

**Handling** Avoid freeze-thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** DJ-1

**Application** Flow Cytometry, ICC, IF, IP, WB

**Clone** 007

**Formulation** Liquid. 0.2 µm filtered solution in PBS. Preservative free. Carrier-free.

**Host** Rabbit

**Immunogen** Recombinant human PARK7 / DJ-1.

**Isotype** IgG

**Purity Detail** Protein A purified.

**Recommendation Dilutions/Conditions** Flow Cytometry (1:25-1:100)Immunocytochemistry / Immunofluorescence (1:20-1:100)Western Blot (1:1'000-1:5'000)Immunocytochemistry / Immunofluorescence (1:300-1:10'000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Human

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



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

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

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

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

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

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

Last modified: October 7, 2024