

DJ-1 monoclonal antibody (3E8)

Researchers cloned DJ-1 protein, consisting of 189 amino acids, from a HeLa cell cDNA library as a oncogene which weakly transformed NIH3T3 cells. Cotransfection with ras or c-myc enhances transformation by DJ-11. DJ-1 expression is induced by growth signals, and it is translocated from the cytoplasm to the nucleus during the S phase without the presence of any nuclear localization signals. In addition to its potential involvement in ras-mediated signal transduction pathways, DJ-1 is known to regulate transcription of androgen-responsive genes, and is sensitive to oxidative stress. DJ-1 is ubiquitously expressed in various human tissues including the brain, with high expression in astrocytes of the frontal cortex and substantia nigra. Homozygous mutations in the DJ-1 gene have been shown to cause early-onset autosomal recessive Parkinson's disease.

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

Citations: 6

[View Online »](#)

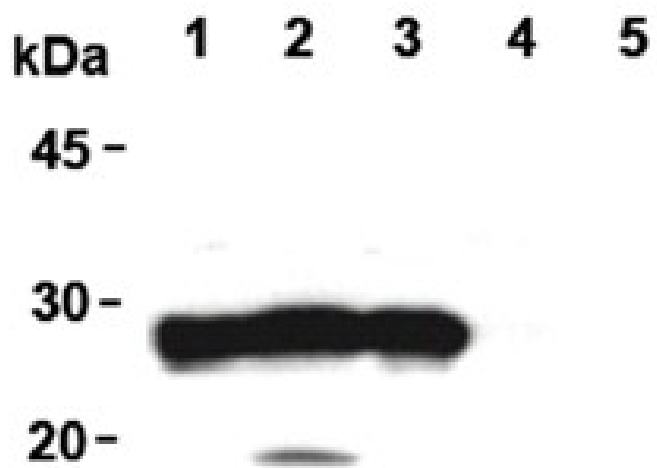
Ordering Information

[Order Online »](#)

ADI-KAM-SA100-E	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Jurkat (1), Raji (2), HeLa (3), NIH/3T3 (4) and PC12 (5) using DJ-1 mAb (3E8).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PARK7, Parkinson disease protein 7
Application	WB
Application Notes	Detects bands of ~28kDa by Western blot.
Clone	3E8
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol.
GenBank ID	BAA09603
Host	Mouse
Immunogen	Recombinant human DJ-1.
Isotype	IgG1
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1-10 µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	Q99497
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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)