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

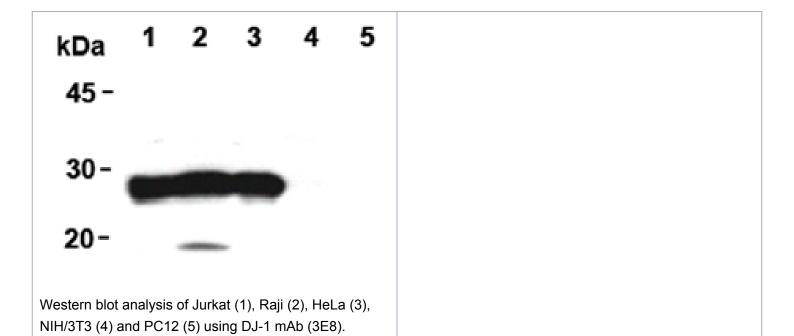
Order Online »

ADI-KAM-SA100-E

100µg

Manuals, SDS & CofA

View Online »



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 NotesDetects 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.

lsotype lgG1

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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com