GAD65/67 monoclonal antibody (9A6)

GAD, glutamic acid decarboxylase, is an integral membrane protein that catalyzes the formation of gamma-aminobutyric acid (GABA), which is an inhibitory neurotransmitter. The biological functions of GAD and GABA extend beyond regulation of neurotransmission to include effects on the immune system and modulation of cell proliferation, protein synthesis, and metabolism.

This antibody is covered by our Worry-Free Guarantee.

Citations: 12

View Online »

Ordering Information

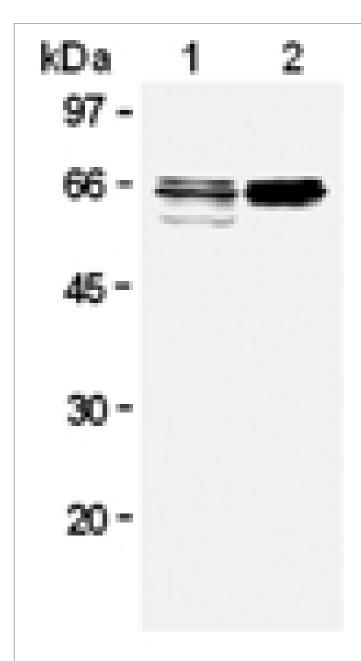
Order Online »

ADI-MSA-225-E

100µg

Manuals, SDS & CofA

View Online »



Western blot analysis of mouse (1) and rat (2) brain extracts, probed with GAD65/67, mAb (9A6).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC (PS), WB

Application Notes Detects bands of ~65kDa and ~67kDa, (corresponding to

GAD65 and GAD67, respectively) by Western blot.

Clone 9A6

Formulation Liquid. In PBS containing 50% glycerol.

Host Mouse

Immunogen Recombinant human GAD65.

lgG1

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections, 1-

10μg/ml)Western Blot (1μg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined

individually for each application.

Species Reactivity Human, Mouse, Rat

UniProt ID Q05329

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

ENZO LIFE SCIENCES, INC.

May 29,942024

usa@enzolifesciences.com

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com