Agrin monoclonal antibody (Agr-131)

Agrin is an essential extracellular matrix component which promotes clustering of nicotinic acetylcholine receptors (nAChRs) and other proteins at the neuromuscular junction during development. Agrin, MuSK, and Rapsyn are all essential components for AChR aggregation, through an unknown mechanism. The C-terminal region of agrin is released into the medium, interacts with receptors on the muscle surface and induces AChR aggregation.

This antibody is covered by our Worry-Free Guarantee.

Citations: 4

View Online »

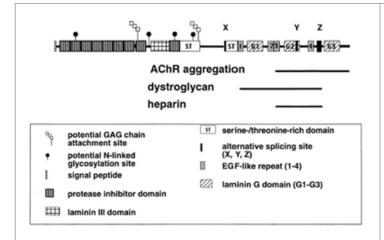
Ordering Information

Order Online »

ADI-AGR-530-D	50µg
ADI-AGR-530-F	200μg

Manuals, SDS & CofA

View Online »



The diagram indicates the domain structure and functional regions of agrin, as well as domains required for AChR aggregation and alpha-dystroglycan and heparin binding.



Immunohistochemistry analysis of rat brain tissue using Agrin, mAb (Agr-131) (prod.no ADI-AGR-530).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ICC, IHC (PS), Microarray, WB

Application Notes Detects a band of ~100kDa by Western blot.

Clone Agr-131

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09%

sodium azide.

GenBank ID M64780

Host Mouse

Immunogen Recombinant rat agrin (C-terminal fragment).

lsotype lgG2a

Purity Detail Protein G affinity purified.

Recommendation Dilutions/ConditionsWestern Blot (1:1,000)Suggested dilutions/conditions may

not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from mouse ascites.

Species Reactivity Mouse, Rat

UniProt ID P25304

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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com