

# Agrin monoclonal antibody (Agr-247)

Agrin is an essential extracellular matrix component which promotes clustering of nicotinic acetylcholine receptors (nAChRs) and other proteins during development at the neuromuscular junction. 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. The central region contains two O-linked glycosylation sites and a domain homologous to domain III of laminin. The N-terminal region anchors agrin to the extracellular matrix via other basal membrane components. This region also contains a protease inhibitor domain and glycosaminoglycan attachment sites, increasing the predicted MW from 200kDa to ~600kDa. The diagram shown indicates the domain structure and functional regions of agrin, as well as domains required for AChR aggregation and alpha-dystroglycan and heparin binding. Various agrin isoforms are generated by alternative splicing at the X, Y and Z sites, and differ in the presence or absence of small inserts. The isoforms can determine the biological activity of agrin and their expression in specific tissues and stages of development. While no difference in functional activity has been detected between splicing variants at site X, insertion of a 4 aa peptide at site Y modestly increases agrin's nAChR clustering activity. Insertion of an 8 aa peptide at splicing site Z increases the clustering activity of soluble rat agrin 10,000 fold.

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

Citations: 1

[View Online »](#)

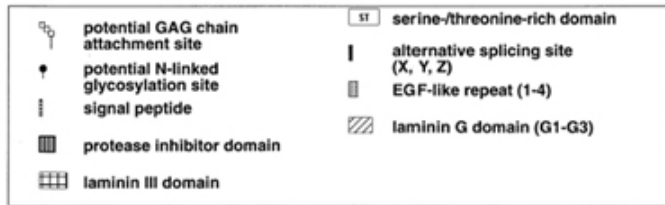
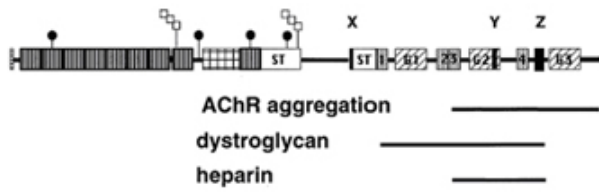
## Ordering Information

[Order Online »](#)

ADI-AGR-540-D	50µg
ADI-AGR-540-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.

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application IHC, WB

Clone Agr-247

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).

Isotype IgG1

Purity Detail Protein G affinity purified.

**Recommendation** Immunohistochemistry (10µg/ml)Suggested dilutions/conditions may not be available for  
**Dilutions/Conditions** 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](#).



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-  
fr@enzolifesciences.com](mailto:info-fr@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)