

# Nicotinic acetylcholine receptor monoclonal antibody (88B)

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

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-802-0020	100µl
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** nAChR

**Application** IHC (FS), IP, WB

**Application Notes** Detects bands of ~60kDa ( $\gamma$  subunit ) and ~65kDa ( $\delta$  subunit) by Western blot. Under non-reducing conditions detects a band of ~130kDa (dimer of  $\delta$  subunit).

**Clone** 88B

**Crossreactivity** Does not cross-react with the  $\alpha 1$  and  $\beta 1$  subunits in muscle or the neuronal nAChRs.

**Formulation** Liquid. Ascites diluted in PBS containing 0.05% sodium azide.

**Host** Mouse

**Immunogen** Purified *Torpedo californica* AChR (acetylcholine receptor).

**Isotype** IgG1

**Recommendation Dilutions/Conditions** Immunohistochemistry (frozen sections, 1:1,000)Western Blot (1:5,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Chicken, Fish, Frog, Mouse, Rat, *Torpedo californica*

**Specificity** Recognizes the intracellular domain of skeletal muscle nAChR  $\gamma$  and  $\delta$  subunits in *Torpedo californica* and the  $\delta$  subunit in mouse, rat, chicken, frog and fish.

Last modified: May 29, 2024



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)