

Bassoon monoclonal antibody (SAP7F407)

Bassoon is a 420 kDa protein that localizes at the presynaptic nerve terminals and is believed to play a role in the structural and functional organization of the synaptic vesicle cycle. Bassoon is concentrated at sites opposite to postsynaptic densities in synaptic terminals and in cultured neurons, it is found to colocalize with GABA (A) and glutamate (GluR1) receptors. These observations suggested that they serve specific functions at synaptic junctions and may be involved in organization of the cytoskeleton at the site of neurotransmitter release.

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

Citations: 2

[View Online »](#)

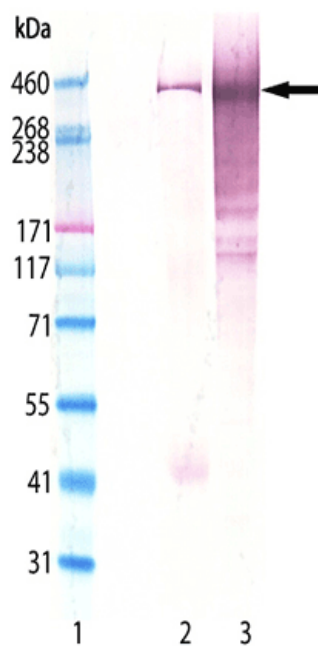
Ordering Information

[Order Online »](#)

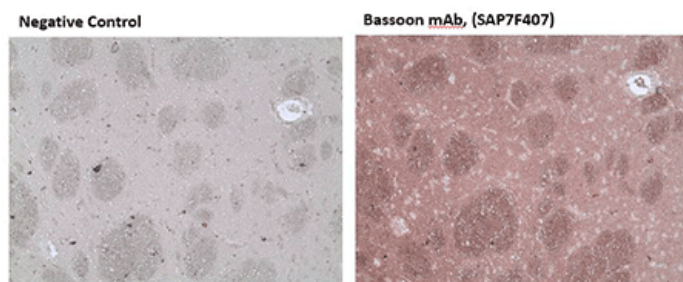
ENZ-ABS666-0050	50µg
ENZ-ABS666-0200	200µg

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of ENZ-ABS666: Lane 1: NovexTMHiMarkTM MW Marker, Lane 2: Mouse Brain lysate (Prod. No. ADI-LYT-MB100), Lane 3: Rat Brain lysate (Prod. No. ADI-LYT-RB100).



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded rat brain tissue stained with Bassoon monoclonal antibody (SAP7F407), (Prod. No. ENZ-ABS666) at 10 μ g/ml.

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ICC, IF, IHC, IP, WB
Clone	SAP7F407
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
Gene/Protein Identifier	29999 (UniGene)
Host	Mouse
Immunogen	Recombinant rat Bassoon.
Isotype	IgG2ak
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (1:400)Immunoprecipitation (2.5µg)Western Blot (1:500 – 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 hybridoma tissue culture supernatant.
Species Reactivity	Bovine, Dog, Human, Mouse, Rat
UniProt ID	O88778
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-
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)