

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: 182

[View Online »](#)

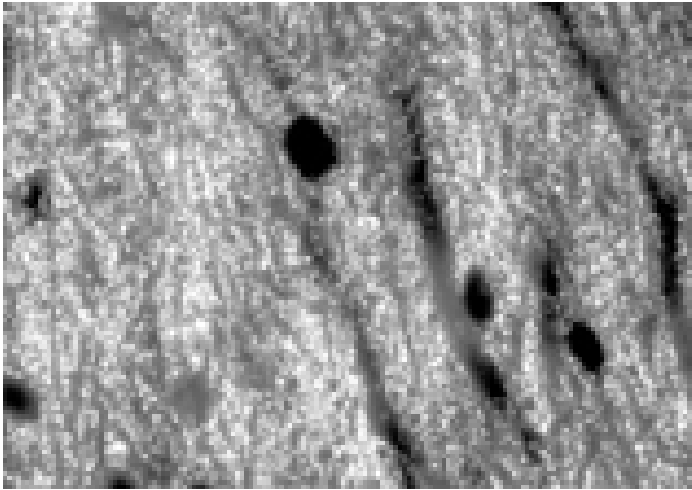
Ordering Information

[Order Online »](#)

ADI-VAM-PS003-D	50µg
ADI-VAM-PS003-F	200µg

Manuals, SDS & CofA

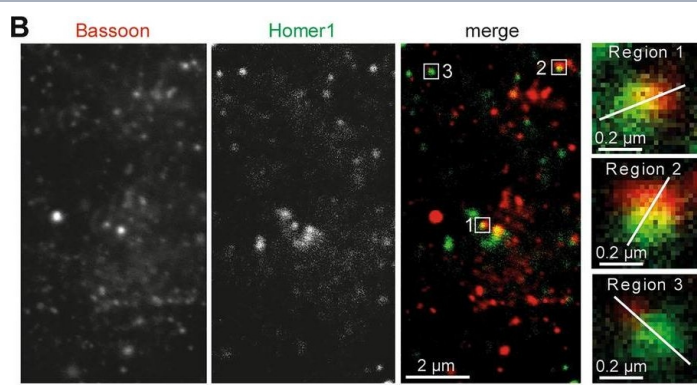
[View Online »](#)



Immunohistochemistry analysis of bassoon localization in the rat brain cerebellar molecular layer using Bassoon, mAb (SAP7F407).

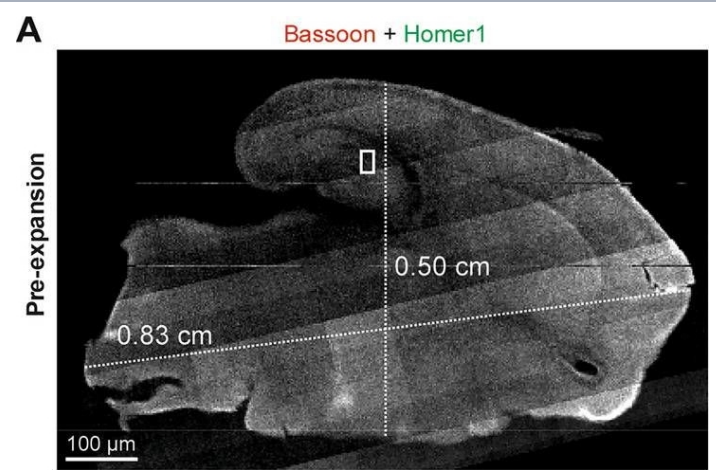


Western blot analysis of Bassoon: Lane 1: MW marker, Lane 2: Rat Brain Tissue Lysate.



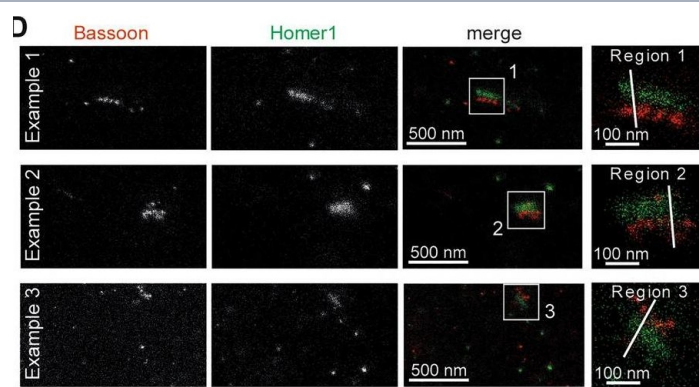
X10ht of 200 μm thick rat brain slices immunostained for the synaptic proteins Bassoon and Homer. (A) Epifluorescence tile image of merged channels displays a non-expanded 200 μm thick rat brain slice, containing the region of the hippocampal formation. (B) STED images of regions of the hippocampal formation (identified with white rectangles in A) show Bassoon (red) and Homer1 (green) in not expanded tissue. Zoom depict the marked regions (1–3). (E) Plots of the line scans over the zoomed areas. (C) Epifluorescence tile acquisition of the same brain slice after expansion labeled with NHS-Fluorescein, depicting the retention of tissue shape and the prolongation of the slice length from 0.83 to 4.98 cm and width from 0.5 to 3.05 cm. (D) Representative STED images of different regions from the C1 region of the hippocampus, showing highly resolved pre- (red) and postsynaptic (green) compartments. (F) The respective line scans are plotted, and the distance between the pre- and postsynaptic compartments is analyzed in (G). N = 20 ROI from two stained rat brain slices for not expanded STED; N = 17 ROI from 2 independent STED-X10ht experiments. (H) The graph shows an expansion factor of 6 with N = 3 independent experiments.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Heat denaturation enables multicolor X10-STED microscopy. *Sci Rep* (2023)



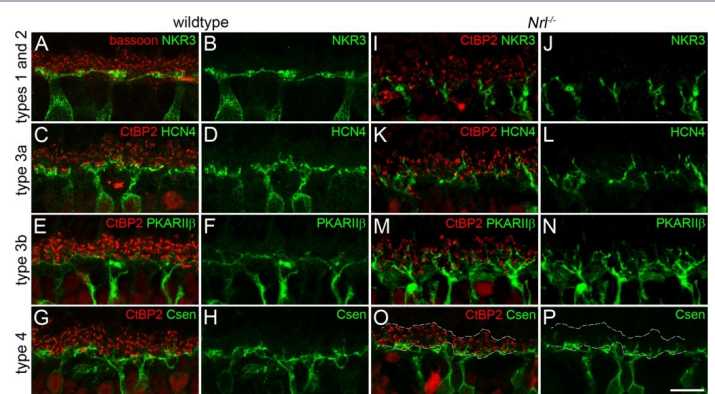
X10ht of 200 μm thick rat brain slices immunostained for the synaptic proteins Bassoon and Homer. (A) Epifluorescence tile image of merged channels displays a non-expanded 200 μm thick rat brain slice, containing the region of the hippocampal formation. (B) STED images of regions of the hippocampal formation (identified with white rectangles in A) show Bassoon (red) and Homer1 (green) in not expanded tissue. Zoom depict the marked regions (1–3). (E) Plots of the line scans over the zoomed areas. (C) Epifluorescence tile acquisition of the same brain slice after expansion labeled with NHS-Fluorescein, depicting the retention of tissue shape and the prolongation of the slice length from 0.83 to 4.98 cm and width from 0.5 to 3.05 cm. (D) Representative STED images of different regions from the C1 region of the hippocampus, showing highly resolved pre- (red) and postsynaptic (green) compartments. (F) The respective line scans are plotted, and the distance between the pre- and postsynaptic compartments is analyzed in (G). N = 20 ROI from two stained rat brain slices for not expanded STED; N = 17 ROI from 2 independent STED-X10ht experiments. (H) The graph shows an expansion factor of 6 with N = 3 independent experiments.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Heat denaturation enables multicolor X10-STED microscopy. *Sci Rep* (2023)



X10ht of 200 μm thick rat brain slices immunostained for the synaptic proteins Bassoon and Homer. (A) Epifluorescence tile image of merged channels displays a non-expanded 200 μm thick rat brain slice, containing the region of the hippocampal formation. (B) STED images of regions of the hippocampal formation (identified with white rectangles in A) show Bassoon (red) and Homer1 (green) in not expanded tissue. Zoom depict the marked regions (1–3). (E) Plots of the line scans over the zoomed areas. (C) Epifluorescence tile acquisition of the same brain slice after expansion labeled with NHS-Fluorescein, depicting the retention of tissue shape and the prolongation of the slice length from 0.83 to 4.98 cm and width from 0.5 to 3.05 cm. (D) Representative STED images of different regions from the C1 region of the hippocampus, showing highly resolved pre- (red) and postsynaptic (green) compartments. (F) The respective line scans are plotted, and the distance between the pre- and postsynaptic compartments is analyzed in (G). $N = 20$ ROI from two stained rat brain slices for not expanded STED; $N = 17$ ROI from 2 independent STED-X10ht experiments. (H) The graph shows an expansion factor of 6 with $N = 3$ independent experiments.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Heat denaturation enables multicolor X10-STED microscopy. *Sci Rep* (2023)



Dendritic stratification patterns of OFF bipolar cells. A-P: Maximum intensity projections of confocal image stacks from vertical cryosections of wild-type (A-H) and *Nrl*^{-/-} (I-P) mice double-labeled with ribbon markers bassoon (A) or CtBP2 (all other). OFF bipolar cells were labeled with antibodies against NK3R (types 1 and 2; A,B,I,J), HCN4 (type 3a; C,D,K,L), PKARII β (type 3b; E,F,M,N), and calsenilin (Csen, type 4; G,H,O,P). Dashed lines in O indicate the inner and outer border of the CtBP2 labeled area (see Fig 2A for more details). Scale bar = 10 μm for all panels.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Dendritic stratification differs among retinal OFF bipolar cell types in the absence of rod photoreceptors. *PLoS One* (2017)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application ICC, IF, IHC, IP, WB

Application Notes Detects a band of ~400kDa by Western blot.

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 ascites.

Species Reactivity Mouse, Rat

Technical Info / Product Notes **Cited samples:**
[*For an overview on cited samples please click here.*](#)

UniProt ID O88778

Worry-free Guarantee

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

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)