

Bradykinin receptor B₂ polyclonal antibody

Bradykinin (BK) is a nine-amino acid peptide produced by kallikrein cleavage of kininogen precursor proteins in tissue and plasma, and is an important regulator of vascular and pain responses to tissue injury. Two receptors (B1 and B2) have been identified as mediators of BK signaling, and are both seven-transmembrane domain-containing GPCRs. The B2 receptor is constitutively expressed in smooth muscle and neurons, whereas B1 expression is induced following injury. B2 receptors signal via Gq alpha subunits, leading to activation of phospholipase Cbeta, hydrolysis of PI, and an intracellular increase in free calcium, although coupling through Gi, Gs, and G12/13 has also been observed.

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

Citations: 1

[View Online »](#)

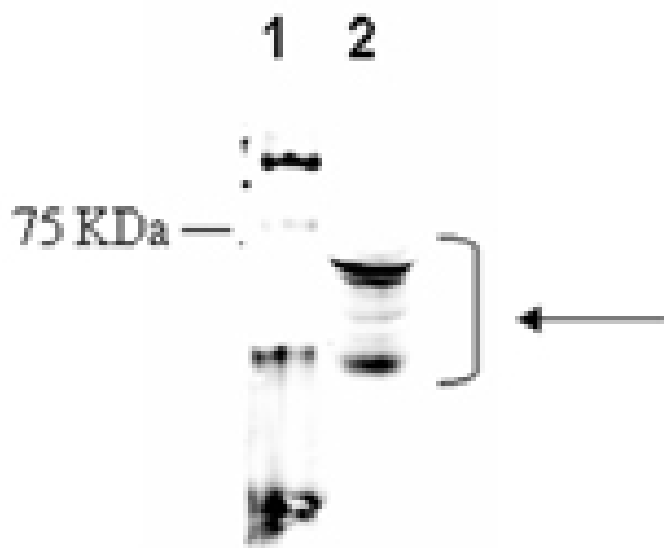
Ordering Information

[Order Online »](#)

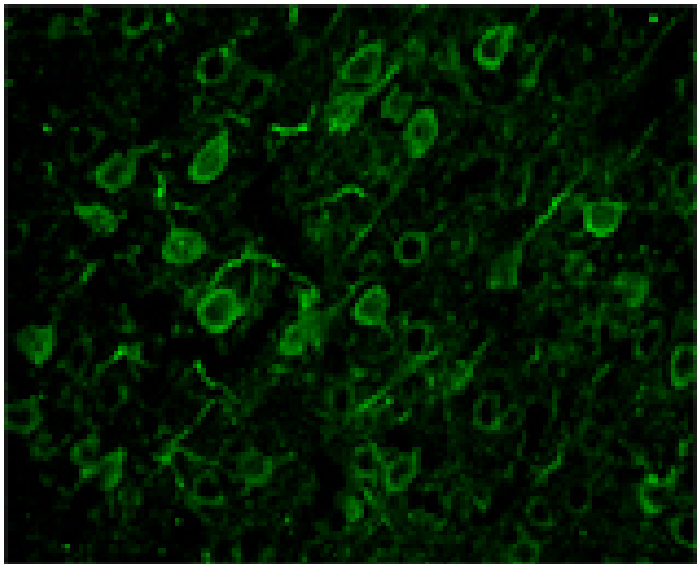
ADI-905-747-100	100µg
-----------------	-------

Manuals, SDS & CofA

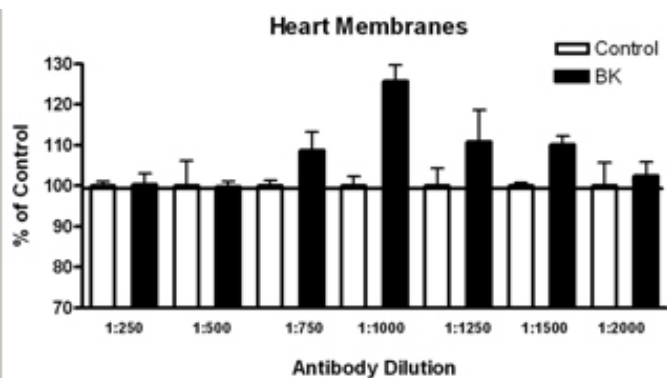
[View Online »](#)



Western blot analysis: MW marker (1) and 20µg rat brain extract (2) probed with Bradykinin Receptor (B2) pAb at 2.2µg/ml.



Immunohistochemistry analysis: Mouse brain fixed with 4% paraformaldehyde, stained with Bradykinin Receptor (B2) pAb at 1:1000, and visualized at 40x.



Membrane ELISA: Lewis rat heart membranes (5µg/well) were treated with 1 µM concentrations of agonist (BK) and probed with Bradykinin Receptor (B2) pAb (1:250 to 1:2000 of a 1µg/µL stock solution) by ELISA. Data from vehicle treated cells were taken as 100%. Results are the mean ± SEM (n=3).

Handling & Storage

Handling Keep on ice at all times.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name B2BKR

Application ELISA, IHC, WB

Application Notes Suitable for membrane-based ELISA.
Predicted MW of ~45kDa. Higher MW species (~50kDa) and intermediates also observed by Western blot which reflect post-translational modification or dimerization of the receptor.

Crossreactivity Predicted species reactivity based on epitope sequence homology: guinea pig, dog.

Formulation Liquid. In PBS containing 50% glycerol and 0.01% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the N-terminus of rat bradykinin receptor B2.

Purity Detail Peptide affinity purified.

Species Reactivity Mouse, Rat

UniProt ID P25023

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)

Last modified: May 29, 2024

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.