Adrenergic receptor β ₂ polyclonal antibody

Adrenergic receptors mediate the hormone and neurotransmitter functions of adrenaline and noradrenaline. Adrenoreceptors function in peripheral tissues and the central nervous system to regulate numerous vascular, metabolic, and neurological functions, and are the target for many therapeutically relevant drugs including those for the treatment of hypertension, asthma, benign prostatic hypertrophy, and depression. Adrenoreceptors are classified by three types and multiple subclasses (alpha1A,1B,1D; alpha2A,2B,2C; and beta1,2,3) based on pharmacological and molecular characteristics.

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

Ordering Information

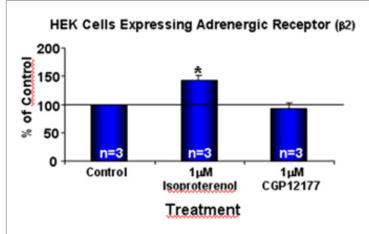
Order Online »

ADI-905-742-100

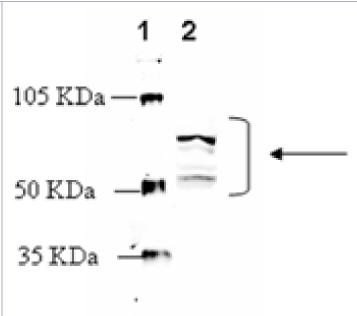
100µg

Manuals, SDS & CofA

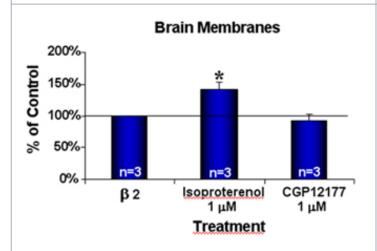
View Online »



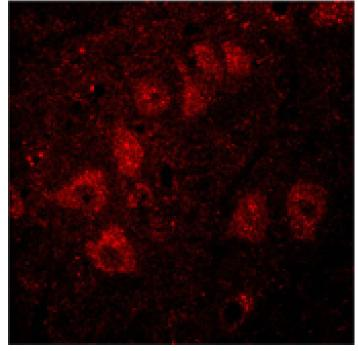
Cellular ELISA: HEK293 cells expressing Adrenergic Receptor $\beta 2$ were treated with indicated concentrations of agonist (Isoproterenol) and antagonist (CPG12177) and probed with Adrenergic Receptor ($\beta 2$) pAb (1:1000 of a stock solution of $1\mu g/\mu L$) by ELISA. Data from vehicle treated cells were taken as 100%. Results are the mean \pm SEM (n=3).



Western blot analysis: MW marker (1) and 20 μ g rat heart membrane extract (2) probed with Adrenergic Receptor (β 2) pAb at 2.5 μ g/ml.



Membrane ELISA: Lewis rat brain membranes (5 μ g/well) were treated with indicated concentrations of agonist (Isoproterenol) and antagonist (CGP12177) and probed with Adrenergic Receptor (β 2) pAb (1:1000 of a stock solution of 1 μ g/ μ L) by ELISA. Data from vehicle treated cells were taken as 100%. Results are the mean \pm SEM (n=3).



Immunohistochemistry analysis: Fluorescent confocal photomicrograph from rat brainstem. Neurons are labeled with Adrenergic Receptor (β 2) pAb and Cy3 conjugated rabbit IgG antibody (H+L).

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 ADRB2, AR β -2

Application ELISA, IHC, WB

Application NotesSuitable for cell-based and membrane-based ELISA.

Predicted MW of ~47kDa. Higher MW species (~65kDa) and intermediates also observed by Western blot which reflect post-translational modification of the receptor.

Crossreactivity Predicted species reactivity based on epitope sequence

homology: chimpanzee, orangutan, tamarin monkey, gorilla, rhesus macaque, horse, hamster, guinea pig, wild

boar, cat.

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

azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to the N-terminus of

mouse adrenergic receptor β2.

Purity Detail Peptide affinity purified.

Species Reactivity Human, Mouse, Rat

UniProt ID P18762

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com