Opioid receptor δ polyclonal antibody

The opioid receptors δ , κ , and μ are members of the seven-transmembrane domain-containing receptor family of G-protein coupled receptors. δ - and μ -receptors serve as functional receptors for endogenous enkephalin and β -endorphin peptide ligands, respectively. δ -opioid receptors are highly expressed in the olfactory bulb, neocortex, caudate putamen and nucleus accumbens, with peripheral expression in the gastro-intestinal tract and vas deferens. δ -receptors inhibit neurotransmitter release resulting in a variety of biological effects including analgesia, motor integration, gastro-intestinal motility, olfaction, respiration, and memory. Neural μ -receptor expression is prominent in the caudate putamen, neocortex, thalamus, nucleus accumbens, hippocampus, and amygdala, as well as in peripheral nervous tissues in the gut. Named for its pharmacological mediation of analgesia by morphine, the μ receptor typically functions as a depressant in physiological processes including respiration, cardiovascular function, gastrointestinal motility, and thermoregulation.

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

Ordering Information

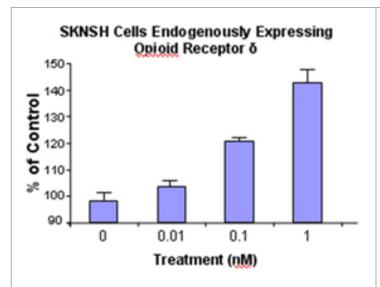
Order Online »

ADI-905-745-100

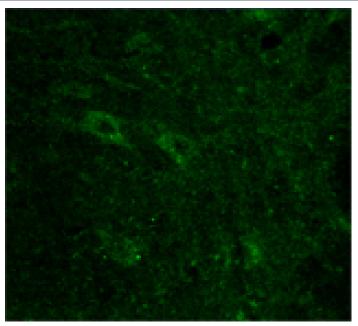
100µg

Manuals, SDS & CofA

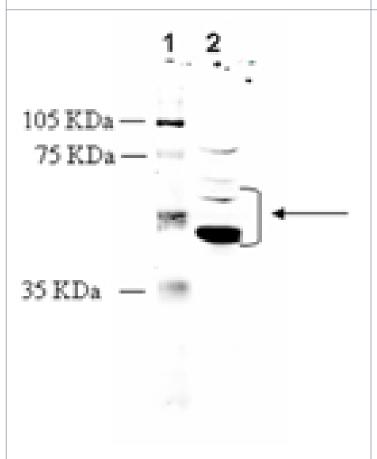
View Online »



Cellular ELISA: SKNSH cells endogenously expressing Opioid Receptor δ were treated with indicated concentrations of agonist (Deltorphin II) and probed with Opioid Receptor δ pAb (1:500 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: Mouse brain cortex, fixed with 4% paraformaldehyde, stained with Opioid Receptor δ pAb at 1:1000, and visualized at 40x.



Western blot analysis: MW marker (1) and 50 μ g rat brain extract (2) probed with Opioid Receptor δ pAb at 1.6 μ g/ml.

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 OPRD1, DOR-1

Application ELISA, IHC, WB

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

Predicted MW of ~40kDa. Higher MW species (~75kDa) and intermediates also observed by Western blot which

may reflect post-translational modification or

heterodimerization of the receptor.

Crossreactivity Predicted species reactivity based on epitope sequence

homology: dog.

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

azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to the N-terminus of

mouse opioid receptor δ .

Purity Detail Peptide affinity purified.

Species Reactivity Mouse, Rat

UniProt ID P32300

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg
Phone: +32 3 466 0420
infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com

Last modified: May 29, 2024