

Opioid receptor μ 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 »](#)

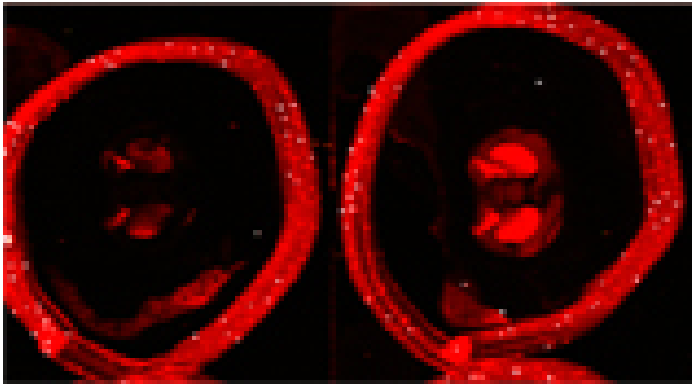
Ordering Information

[Order Online »](#)

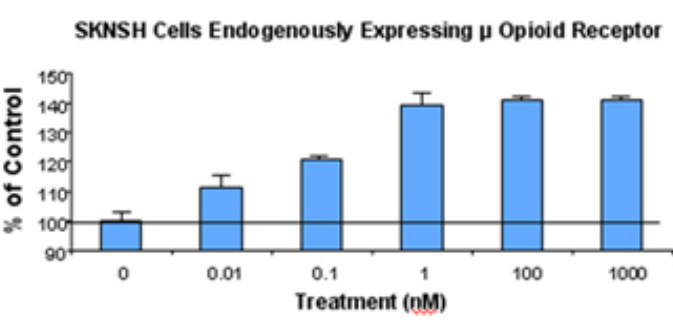
ADI-905-744-100	100 μ g
-----------------	-------------

Manuals, SDS & CofA

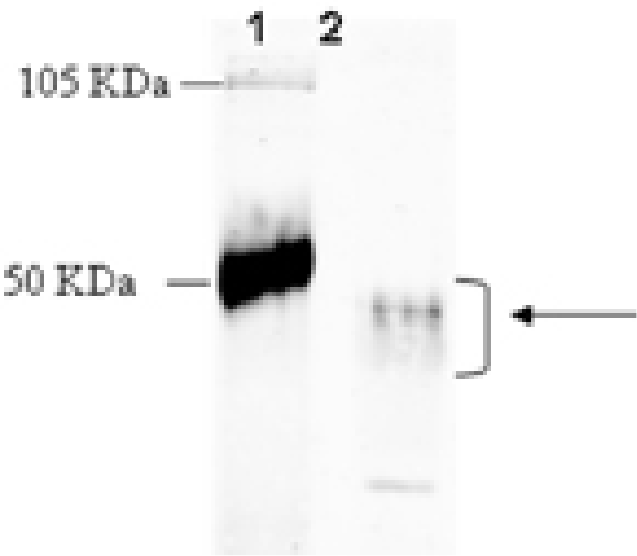
[View Online »](#)



Immunohistochemistry analysis: Mouse brain sections untreated (left) or morphine-treated (right) were stained with Opioid Receptor μ pAb.



Cellular ELISA: SKNSH cells endogenously expressing Opioid Receptor μ were treated with indicated concentrations of agonist (DAMGO) and probed with Opioid Receptor μ pAb (1:500 of a stock solution of $1\mu\text{g}/\mu\text{l}$) 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 $20\mu\text{g}$ rat brain extract (2) probed with Opioid Receptor μ pAb at $1.3\mu\text{g}/\text{ml}$.

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 OPRM1, MOR-1

Application ELISA, IHC, WB

Application Notes Suitable for cell-based and membrane-based ELISA.
Predicted MW of ~44kDa. Higher MW species (~50kDa) also observed by Western blot which may reflect post-translational modification of the receptor.

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 mouse Opioid Receptor μ .

Purity Detail Peptide affinity purified.

Species Reactivity Human, Mouse, Rat

UniProt ID P42866

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)