

PGF_{2α} receptor (FP) polyclonal antibody

Arachidonic acid is converted by cyclooxygenase (COX) to prostaglandin H₂ (PGH₂), which is further converted to the derivative prostanoids PGD₂, PGE₂, PGF_{2α}, PGI₂ (prostacyclin) and thromboxane A₂ (TXA₂). Five classes of prostanoid receptors are recognized and named DP, EP, FP, IP and TP receptors corresponding to the agonist properties of the associated prostanoids. FP receptors are found in smooth muscle where they mediate contraction, and are believed to be coupled via Gq/11 subunits to stimulate phosphatidyl inositol (PI) hydrolysis. The FP agonist Fluprostenol (a PGF_{2α} analogue) is used in animal husbandry to synchronize oestrus and induce parturition. FP receptors are known to mediate intraocular pressure, and FP agonists such as latanoprost and travoprost are drugs commonly used to treat glaucoma.

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

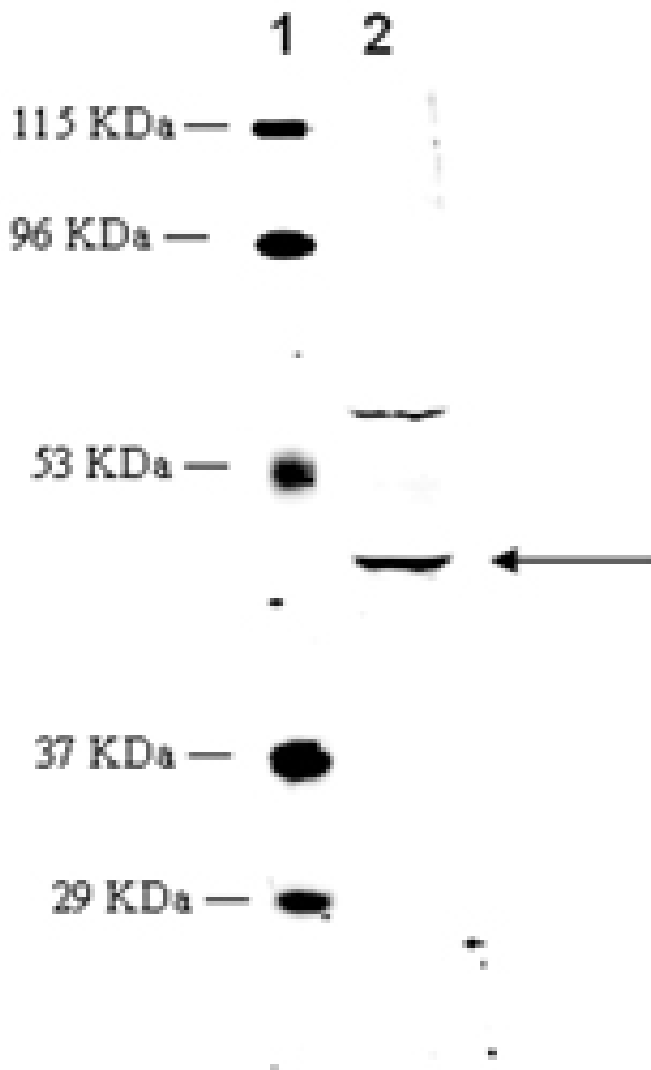
Ordering Information

[Order Online »](#)

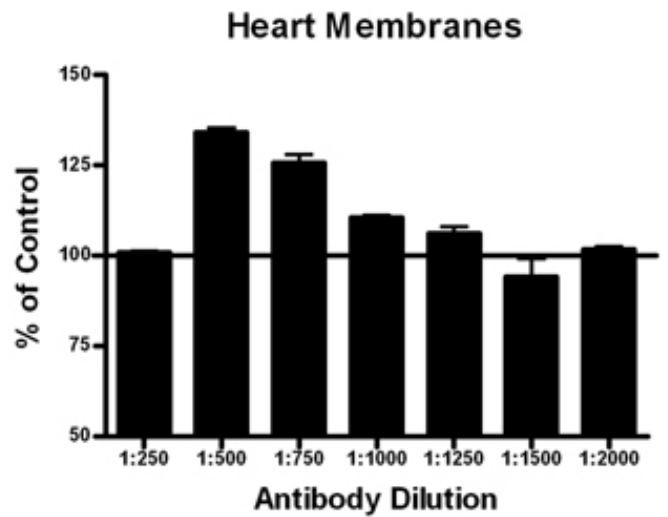
ADI-905-901-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of MW marker (1) and 25µg rat kidney extract (2), probed with PGF2α Receptor (FP) pAb.



Membrane ELISA: Lewis rat striatum membranes (5µg/well) were treated with 1 µM concentrations of agonist (cyclosin) and probed with PGF2α Receptor (FP) 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=4).

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 PTGFR, Prostaglandin F2α receptor

Application ELISA, IHC, WB

Application Notes Suitable for membrane-based ELISA.
Predicted MW of ~41kDa, observed by Western blot ~45kDa. Higher MW species (~60kDa) and intermediates also observed by Western blot which reflect post-translational modification of the receptor.

Crossreactivity Predicted species reactivity based on epitope sequence homology: wild boar, 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 PGF2 Receptor.

Purity Detail Peptide affinity purified.

Species Reactivity Mouse, Rat

UniProt ID P43118

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)