8-iso-PGF_{2α} polyclonal antibody

Prostaglandin F2alpha (PGF2alpha) is formed in a variety of cells from PGH2, which itself is synthesized from arachidonic acid by the enzyme prostaglandin synthetase. PGF2alpha is often viewed as an antagonist to PGE2 due to their opposing effects on various tissues. PGF2alpha is a potent bronchoconstrictor and has been implicated in asthma attacks. PGF2alpha is also involved in reproductive functions including corpus luteum regulation, uterine contractions, and sperm motility. This has led to its use in terminating pregnancies and inducing labor at term. High levels of PGF2alpha have also been associated with preeclampsia.

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

Ordering Information

Order Online »

ADI-905-015-100

100µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-iso-Prostaglandin F2α

Application ELISA

Formulation Liquid. Serum.

Host Rabbit

Immunogen 8-iso-Prostaglandin F2α conjugated to porcine

thyroglobulin.

Source Whole rabbit serum.

Species Reactivity Species independent

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

•

Last modified: May 29, 2024



eu@enzolifesciences.com