

Prostaglandin F2 α . tromethamine

Prostanoid receptor ligand

A major uterine luteolytic prostaglandin. Endogenous agonist of the prostanoid FP receptor ($K_d=1.0$ nM). Inhibits differentiation of 3T3-L1 preadipocytes. The product is not sterile.

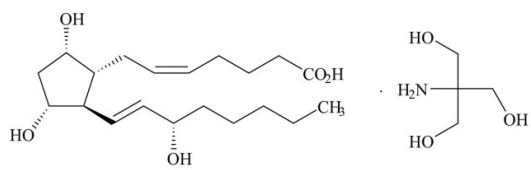
Ordering Information

[Order Online »](#)

BML-PG009-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name PGF2α . THAM

Appearance White solid.

CAS 38562-01-5

Couple Target Prostaglandin receptor, Prostanoid receptor

Couple Type Ligand

Formula $C_{20}H_{34}O_5 \cdot C_4H_{11}NO_3$

Identity Determined by NMR.

MW 475.6

Purity ≥99% (HPLC)

Solubility Soluble in water.

Source Synthetic.

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@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)