

Prostaglandin F_{2α}

Prostanoid receptor ligand

A major uterine luteolytic prostaglandin. Endogenous agonist of the prostanoid FP receptor ($K_D=1.0\text{nM}$). Inhibits differentiation of 3T3-L1 preadipocytes.

Citations: 4

[View Online »](#)

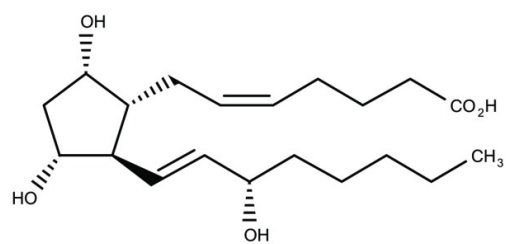
Ordering Information

[Order Online »](#)

BML-PG008-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Dinoprost, 9 α ,11 α ,15S-Trihydroxy-prosta-5Z,13E-dien-1-oic acid, PGF2 α
Appearance	Low melting white solid or colorless oil.
CAS	551-11-1
Couple Target	Prostaglandin receptor, Prostanoid receptor
Couple Type	Ligand
Formula	C ₂₀ H ₃₄ O ₅
MW	354.5
Purity	≥99% (TLC)
Solubility	Soluble in 100% ethanol (50mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com