

Prostaglandin A₁

Prostanoid receptor ligand

Biosynthesis from dihomo- γ -linolenic acid. Stimulates renin release. Antitumor activity. Inhibits NF κ B activation in several human cell lines. Induces the synthesis of heat shock proteins. Activates PPARs in human osteosarcoma cells at 10 μ M.

Citations: 7

[View Online »](#)

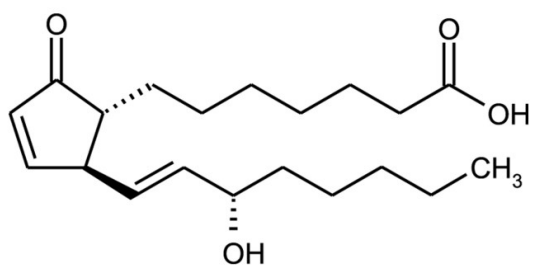
Ordering Information

[Order Online »](#)

BML-PG001-0010	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.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PGA1
Appearance	White solid.
CAS	14152-28-4
Couple Target	Prostanoid receptor
Couple Type	Ligand
Formula	$C_{20}H_{32}O_4$
MW	336.5
Purity	≥98% (HPLC)
Solubility	Soluble in 100% ethanol (50mg/ml), DMSO (50mg/ml) or PBS pH7.2 (2.4mg/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-fr@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