

U-46619

Prostaglandin H₂/thromboxane A₂ receptor agonist

U-46619 is a Prostaglandin H₂/thromboxane A₂ analog and a potent thromboxane A₂ receptor (TP) agonist ($EC_{50} = 0.035\mu M$). U-46619 induces platelet shape change and aggregation at 32nM. U-46619 also activates p38 MAPK.

Citations: 34

[View Online »](#)

Ordering Information

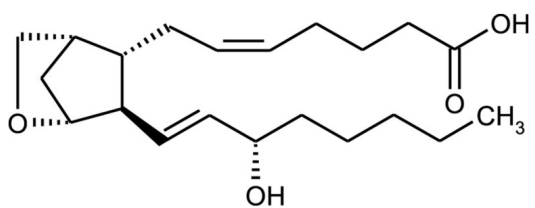
[Order Online »](#)

BML-PG023-0001	1mg
BML-PG023-0010	10mg

Manuals, SDS & CofA

[View Online »](#)

- Potent TP agonist
- Used in cardiovascular studies
- highly cited



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 years after receipt when stored at -20°C. We do not recommend storing aqueous solutions for more than one day.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	9,11-Dideoxy-9 α ,11 α -methanoepoxy prostaglandin F2 α , 9,11-Dideoxy-9 α ,11 α -methanoepoxy prosta5Z,13E-dien-1-oic acid
Appearance	Pale yellow oil.
CAS	56985-40-1
Couple Target	p38, Prostanoid receptor, Thromboxane receptor
Couple Type	Activator, Ligand
Formula	C ₂₁ H ₃₄ O ₄
Identity	Identity determined by 1H-NMR.
MW	350.5
Purity	≥98% (TLC)
Solubility	Soluble in DMF (100mg/ml), DMSO, 100% ethanol, DMSO, or PBS, pH 7.2 (2mg/ml).
Source	Synthetic.

Technical Info / Product Notes

Note: Product is not sterile.



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

Last modified: May 29, 2024