

# 13,14-Dihydro- prostaglandin E<sub>1</sub>

Prostanoid receptor ligand

Circulating metabolite of PGE<sub>1</sub> in man. Vasodilator equipotent with PGE<sub>12</sub>.  
Inhibits human platelet aggregation. Antimitotic activity.

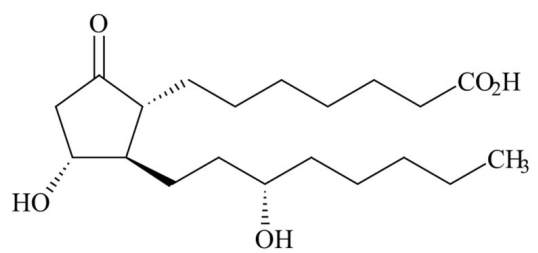
## Ordering Information

[Order Online »](#)

BML-PG048-0001	1mg
----------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -80°C for up to 1 year. Store solutions at -80°C for up to 3 months.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	13,14-Dihydro-PGE1
<b>Appearance</b>	Colorless to pale yellow oil
<b>CAS</b>	19313-28-1
<b>Couple Target</b>	Prostanoid receptor
<b>Couple Type</b>	Ligand
<b>Formula</b>	$C_{20}H_{36}O_5$
<b>MW</b>	356.5
<b>Purity</b>	≥98% (TLC)

**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-be@enzolifesciences.com](mailto:info-be@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)