

# Misoprostol, free acid

## Prostanoid receptor ligand

Free acid form of the cytoprotective PGE<sub>1</sub> analog misoprostol. Misoprostol prevents NSAID-induced gastric ulceration. The free acid form is active in cell free assays and in cell culture and receptor preparations (receptor binding assays etc.). The methyl ester form (misoprostol, Prod. No. BML-PG051) is a prodrug and must be hydrolyzed via intracellular or other esterases in order to display biological activity.

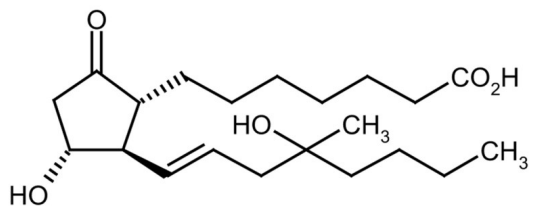
### Ordering Information

[Order Online »](#)

BML-PG052-0001	1mg
----------------	-----

### 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** (±)-15-Deoxy-(16R,S)-16-hydroxy-16-methyl-prostaglandin E1

**Appearance** Yellow viscous oil.

**CAS** 112137-89-0

**Couple Target** Prostanoid receptor

**Couple Type** Ligand

**Formula**  $C_{21}H_{36}O_5$

**MW** 368.5

**Purity** ≥98% (TLC)

**Solubility** Soluble in 100% ethanol (>15mg/ml) or DMSO (>20mg/ml).

**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)