

# Prostaglandin B<sub>2</sub>

Prostanoid receptor ligand

Most abundant prostaglandin released from osteoblasts. Inhibits glucagon-stimulated hepatocyte cAMP accumulation.

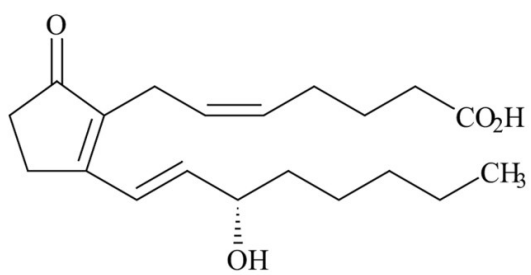
## Ordering Information

[Order Online »](#)

|                |      |
|----------------|------|
| BML-PG004-0010 | 10mg |
|----------------|------|

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |                     |
|------------------|---------------------|
| Alternative Name | PGB2                |
| Appearance       | Waxy solid.         |
| CAS              | 13367-85-6          |
| Couple Target    | Prostanoid receptor |
| Couple Type      | Ligand              |
| Formula          | $C_{20}H_{30}O_4$   |
| MW               | 334.5               |
| Purity           | ≥99%                |
| Solubility       | Soluble in ethanol. |

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)