

Benzylphosphonic acid-(AM)₂

PTP inhibitor

Benzylphosphonic acid-(AM)₂ is a bis-acetoxymethyl ester (AM) derivative of benzylphosphonic acid which is cell permeable and is converted to the active phosphonic acid via intracellular esterase-catalyzed cleavage of the AM esters.

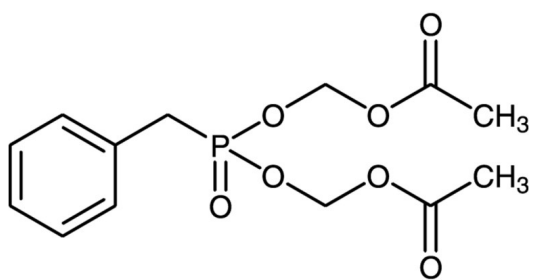
Ordering Information

[Order Online »](#)

BML-EI245-0010	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 -20C for up to 1 year. Store solutions at -20°C for up to 3 months.
Handling	Avoid alkaline media (pH>8).
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	200716-68-3
Couple Target	PTP
Couple Type	Inhibitor
Formula	$C_{13}H_{17}O_7P$
MW	316.3
MeltingPoint	38-39°C
Purity	≥98% (TLC)
Solubility	Soluble in ethanol (20 mg/ml) or DMSO (20 mg/ml).



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)