

Hydroxy-2-naphthalenylmethylphosphonic acid

Insulin receptor inhibitor

HNMPA is a membrane impermeable inhibitor of insulin receptor tyrosine kinase ($IC_{50} = 100\mu M$); in rat adipocytes ($IC_{50} = 10\mu M$). Inhibits serine and tyrosine autophosphorylation by the human insulin receptor.

Citations: 2

[View Online »](#)

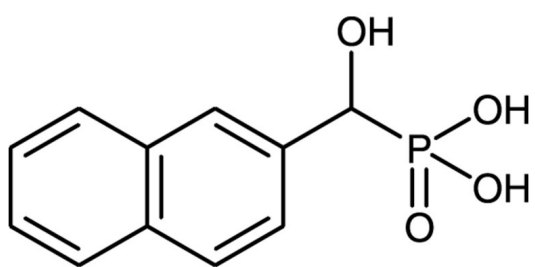
Ordering Information

[Order Online »](#)

BML-EI247-0005	5mg
BML-EI247-0025	25mg

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	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	132541-52-7
Couple Target	Insulin receptor
Couple Type	Inhibitor
Formula	$C_{11}H_{11}O_4P$
MW	238.2
MeltingPoint	≥250°C dec.
Purity	≥98% (HPLC)
Solubility	Soluble in 100% ethanol (25mg/ml), DMSO (25ml/ml) or 5% Sodium bicarbonate (25mg/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)