

HNMPA-(AM)₃

Insulin receptor inhibitor

HNMPA-(AM)₃ is a cell-permeable analog of HNMPA. It inhibits insulin receptor tyrosine kinase activity (IC_{50} =100 μ M) and insulin-stimulated glucose oxidation in isolated rat adipocytes (IC_{50} =10 μ M). It has no effect on PKA (at concs. up to 1 mM) or PKC (at concs. up to 420 μ M). Inhibits both serine and tyrosine autophosphorylation by the human insulin receptor. Does not affect protein kinase C (PKC) or cAMP-dependent protein kinase (PKA).

Citations: 7

[View Online »](#)

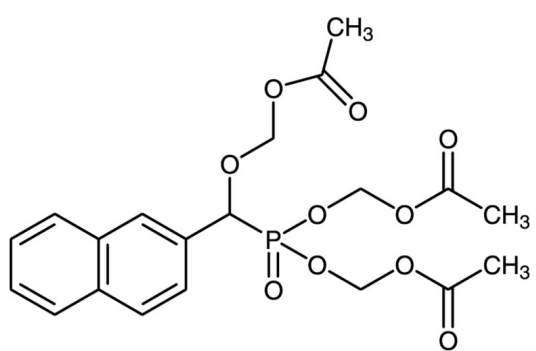
Ordering Information

[Order Online »](#)

BML-EI248-0005	5mg
BML-EI248-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions are stable for up to 3 months when stored at -20°C. Avoid alkaline medium, pH>8.0
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Hydroxy-2-naphthalenylmethylphosphonic acid tri-acetoxymethyl ester
Appearance	Colorless oil.
CAS	120944-03-8
Couple Target	Insulin receptor
Couple Type	Inhibitor
Formula	$C_{20}H_{23}O_{10}P$
Identity	Determined by NMR.
MW	454.4
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (>20mg/ml) or 100% ethanol (>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)