

Phaclofen

GABA_B receptor antagonist

Selective GABA_B receptor antagonist. Antagonizes the postsynaptic action of baclofen (Prod. No. BML-EA128).

Citations: 6

[View Online »](#)

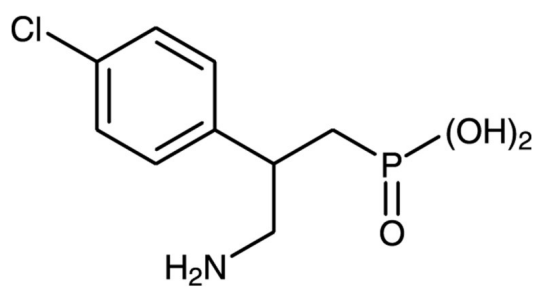
Ordering Information

[Order Online »](#)

BML-EA127-0025	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 room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(±)-3-Amino-2-(4-chlorophenyl)-propylphosphonic acid
Appearance	White solid.
CAS	114012-12-3
Couple Target	GABA receptor
Couple Type	Ligand
Formula	$C_9H_{13}ClNO_3P$
Identity	Determined by 1H -NMR and MS.
MW	249.6
Purity	≥98% (HPLC)
Solubility	Soluble in water, methanol or DMSO.

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)