

4,4-

Pentamethylenepiperidine . HCl

M2 proton channel inhibitor

A novel member of the spiro-piperidine class of M2 proton channel inhibitors similar to BL-1743. Displays increased potency ($IC_{50}=0.92\mu M$) compared with amantadine.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-CM128-0050	50mg
----------------	------

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-Azaspiro[5.5]undecane . HCl
Appearance	White solid.
CAS	180-44-9
Couple Target	M2 proton channel
Couple Type	Inhibitor
Formula	$C_{10}H_{19}N \cdot HCl$
MW	153.2 . 36.5
Purity	≥97% (HPLC)
Solubility	Soluble in water (>25mg/ml), DMSO (>25mg/ml) or 100% ethanol (>25mg/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)