

Compound 48/80 . trihydrochloride

G protein activator

Compound 48/80 is an oligomeric mixture of condensation products of N-methyl-p-methoxyphenethylamine and formaldehyde. It activates G proteins by a mechanism analogous to that of mastoparan. It inhibits calmodulin and human platelet PLC and exhibits a concentration dependent biphasic modulation of human platelet PLA_2 .

Citations: 13

[View Online »](#)

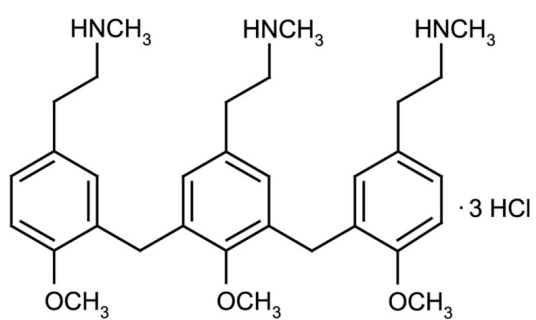
Ordering Information

[Order Online »](#)

ALX-420-008-M100	100mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 3 months when stored at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	94724-12-6
Couple Target	Calmodulin, G Protein, PLC
Couple Type	Activator, Inhibitor
Formula	$C_{32}H_{46}N_3O_3 \cdot 3HCl$
Identity	Determined by 1H-NMR.
MW	520.7 . 109.4
Solubility	Soluble in water (100mg/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-
be@enzolifesciences.com](mailto:info-be@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)