

# 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 »](#)

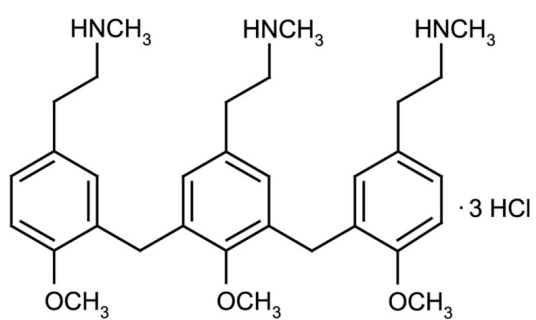
Ordering Information

[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 <sup>1</sup> H-NMR.
MW	520.7 . 109.4
Solubility	Soluble in water (100mg/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-  
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)