

WS-12

TRMP8 agonist

Cooling agent. High-affinity TRPM8 (cold menthol receptor 1; CMR1) agonist. Has an EC_{50} value ~2'000 times lower than that of (-)-menthol (Prod. No. ALX-420-042).

Citations: 2

[View Online »](#)

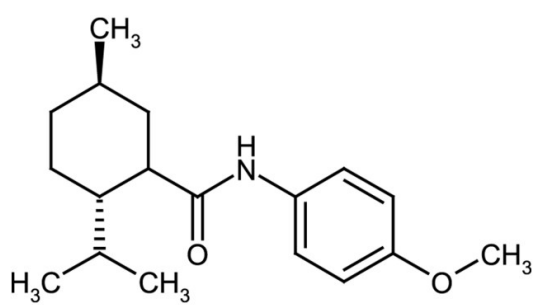
Ordering Information

[Order Online »](#)

ALX-420-041-M025	25mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (2S,5R)-2-Isopropyl-N-(4-methoxyphenyl)-5-methylcyclohexanecarboximide

Appearance White to off-white powder.

CAS 68489-09-8

Couple Target TRP channel

Couple Type Activator, Ligand

Formula $C_{18}H_{27}NO_2$

Identity Determined by 1H -NMR and MS.

MW 289.4

Purity $\geq 98\%$ (HPLC)

Solubility Soluble in 100% ethanol, methanol, ethyl acetate, , dimethyl formamide or acetonitrile; insoluble in water or DMSO.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com