

A-967079

TRPA channel blocker

Potent and selective TRPA1 channel blocker. Exhibits greater than 1000-fold selectivity for TRPA1 over other TRP channels. Exhibits analgesic actions.

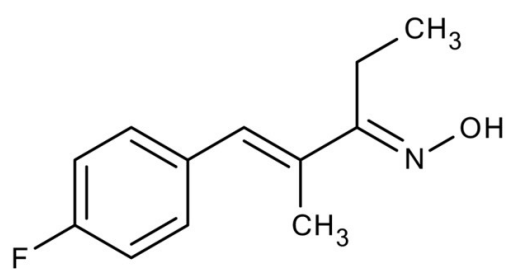
Ordering Information

[Order Online »](#)

ENZ-CHM210-1000	1g
-----------------	----

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (1E,3E)-1-(4-Fluorophenyl)-2-methyl-1-penten-3-one oxime

Appearance White solid.

CAS 1170613-55-4

Couple Target TRP channel

Couple Type Blocker

Formula $C_{12}H_{14}NOF$

MW 207.3

Purity ≥99% (TLC)

Solubility Soluble in DMSO, DMF or ethanol.

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)