

MSK-195

TRP channel agonist

Potent analgesic with $EC_{50}=0.96\mu\text{g/kg}$ in the acetic acid-induced writhing test. RTX binding affinity ($K_i=603\text{nM}$), agonism (calcium influx; $EC_{50}=240\text{nM}$).

Citations: 1

[View Online »](#)

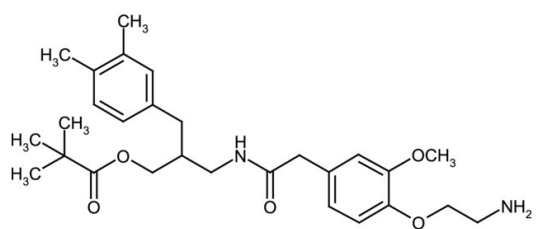
Ordering Information

[Order Online »](#)

ALX-550-396-M005	5mg
------------------	-----

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 N-[2-(3,4-Dimethylbenzyl)-3-(pivalyloxy)propyl]-2-[4-(2-aminoethoxy)-3-methoxyphenyl]acetamide

Appearance White solid.

CAS 289902-82-5

Couple Target TRP channel

Couple Type Activator

Formula $C_{28}H_{40}N_2O_5$

MW 484.6

Purity ≥98%

Solubility Soluble in methanol, 100% ethanol, DMSO or other organic solvents; insoluble in water.

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-
be@enzolifesciences.com](mailto:info-be@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)