

AP-18

TRPA1 blocker

AP-18 is a novel TRPA1 channel blocker (CHO cells, $IC_{50} = 3.1\mu M$; human and mouse cell lines, $IC_{50} = 4.5\mu M$). It reduces cinnamaldehyde-induced nociception *in vivo* and blocks cold- and mustard oil-induced activation of mouse TRPA1 but not capsaicin-induced activation demonstrating selectivity vs TRPV1. AP-18 was unable to block activation of TRPV1, TRPV2, TRPV3, TRPV4 or TRPM8 at concentrations up to $50\mu M$. It reverses CFA-induced mechanical hyperalgesia in mice.

Citations: 4

[View Online »](#)

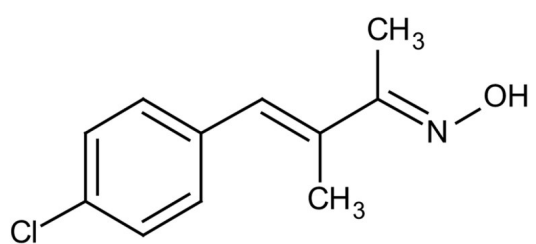
Ordering Information

[Order Online »](#)

BML-CA239-0025	25mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 4-(4-Chlorophenyl)-3-methylbut-3-en-2-oxime

Appearance White solid.

CAS 55224-94-7

Couple Target TRP channel

Couple Type Blocker

Formula $C_{11}H_{12}ClNO$

MW 209.7

Purity $\geq 98\%$ (TLC)

Solubility Soluble in DMSO (>25mg/ml); ethanol (25mg/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)