

Polygodial

TRPA1 activator

Produced from the bark of *Drymis winteri*. Naturally occurring drimane sesquiterpene found in water pepper (*Polygonum hydropiper*) which has been used in folk medicine for a variety of indications. Potent and selective activator of TRPA1 ($EC_{50} = 400\text{nM}$).

Citations: 5

[View Online »](#)

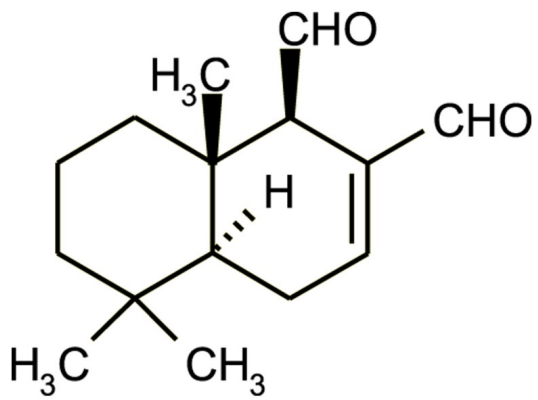
Ordering Information

[Order Online »](#)

BML-CA242-0005	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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (1R,4aS,8aS)-1,4,4a,5,6,7,8,8a-Octahydro-5,5,8a-trimethyl-1,2-naphthalenedicarboxaldehyde

Appearance White to off white solid or waxy solid.

CAS 6754-20-7

Couple Target TRP channel

Couple Type Activator

Formula $C_{15}H_{22}O_2$

MW 234.3

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (25 mg/ml) or ethanol (25 mg/ml).

Source Isolated from *Polygonum hydropiper*.

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

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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