

Kavain (+/-)

Na⁺ channel blocker

Kavain is one constituent of kava-kava, an herbal drug of the South Pacific islands, which displays anti-convulsive, anxiolytic, analgesic and muscle relaxing properties. Kavain inhibits voltage-dependent Na⁺ channels in rat CA1 hippocampal neurons. It displays protective effects in the MPTP mouse model of Parkinson's disease but potentiates NMDA toxicity in the developing hippocampus.

Citations: 3

[View Online »](#)

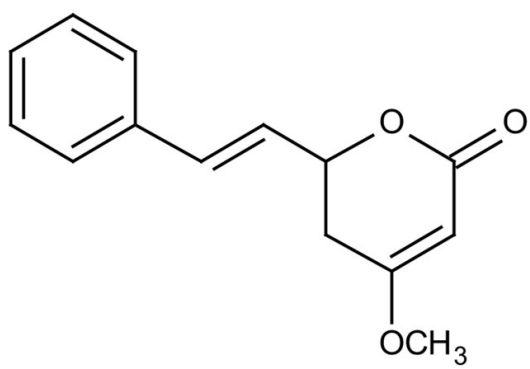
Ordering Information

[Order Online »](#)

BML-NA136-0050	50mg
BML-NA136-0250	250mg

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-Methoxy-6-styryl-5,6-dihydro- α -pyrone

CAS 500-64-1

Couple Target Sodium channel

Couple Type Blocker

Formula $C_{14}H_{14}O_3$

MW 230.3

Purity $\geq 98\%$

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)