

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

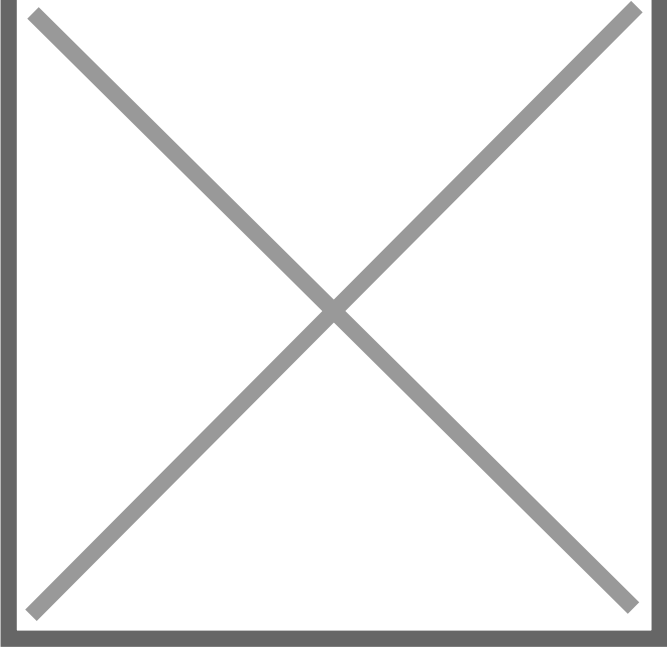
[Order Online »](#)

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

Manuals, SDS & CofA

[View Online »](#)

Image not found or type unknown



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	≥98%

Last modified: May 29, 2024

Image not found or type unknown
ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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