

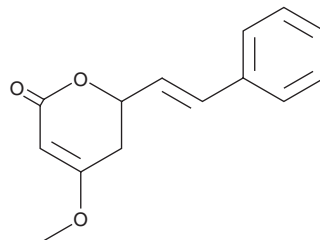
PRODUCT INFORMATION



(±)-Kavain

Item No. 39921

CAS Registry No.: 3155-48-4
Formal Name: 5,6-dihydro-4-methoxy-6-[(1E)-2-phenylethenyl]-2H-pyran-2-one
Synonym: DL-Kavain
MF: C₁₄H₁₄O₃
FW: 230.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

(±)-Kavain (Item No. 39921) is an analytical reference standard characterized as a kava pyrone.¹ (±)-Kavain is a racemic mixture of (+)-kavain (Item Nos. 33881 | 26524) and (-)-kavain. This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 26524).

Reference

- Schirmmacher, K., Büsselberg, D., Langosch, J.M., *et al.* Effects of (±)-kavain on voltage-activated inward currents of dorsal root ganglion cells from neonatal rats. *Eur. Neuropsychopharmacol.* **9(1-2)**, 171-176 (1999).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/30/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM