

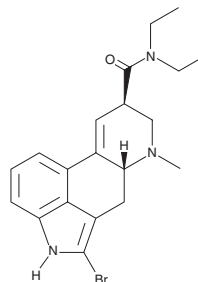
PRODUCT INFORMATION



2-bromo LSD

Item No. 40633

CAS Registry No.: 478-84-2
Formal Name: 2-bromo-9,10-didehydro-N,N-diethyl-6-methyl-ergoline-8 β -carboxamide
Synonyms: BOL-148, 2-bromo Lysergic Acid diethylamide
MF: C₂₀H₂₄BrN₃O
FW: 402.3
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 2 years



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

Description

2-bromo LSD (Item No. 40633) is an analytical reference standard categorized as a lysergamide. Unlike LSD (Item Nos. 35189 | 28289 | 37663), 2-bromo LSD does not induce myoclonic jumping activity in guinea pigs, indicating a lack of hallucinogenic potential.¹ This product is intended for research and forensic applications.

Reference

1. Carvey, P., Nausieda, P., Weertz, R., *et al.* LSD and other related hallucinogens elicit myoclonic jumping behavior in the guinea pig. *Prog. Neuropsychopharmacol. Biol. Psychiatry* **13**(1-2), 199-210 (1989).

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, 06/21/2024

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