

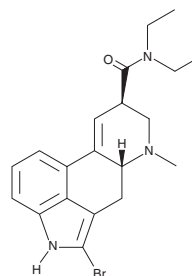
PRODUCT INFORMATION



2-bromo LSD (solution)

Item No. 36791

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, Bromolysergic Acid diethylamide, Bromolysergide
MF: C₂₀H₂₄BrN₃O
FW: 402.3
Supplied as: A 100 μ g/ml solution in acetonitrile
Storage: -20°C
Stability: \geq 3 years



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

Description

2-bromo LSD (solution) (Item No. 36791) is an analytical reference material categorized as a lysergamide. Unlike LSD (Item No. 28289), 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.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

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, 06/29/2022

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