

PRODUCT INFORMATION



GC-MS LSD Isobar Mixture

Item No. 35874

Supplied as: A solution in acetonitrile (1 mg/ml of each compound)

Fill volume: >1 ml

Storage: -20°C

Stability: ≥5 years

Special Conditions: Vortex or sonicate before use

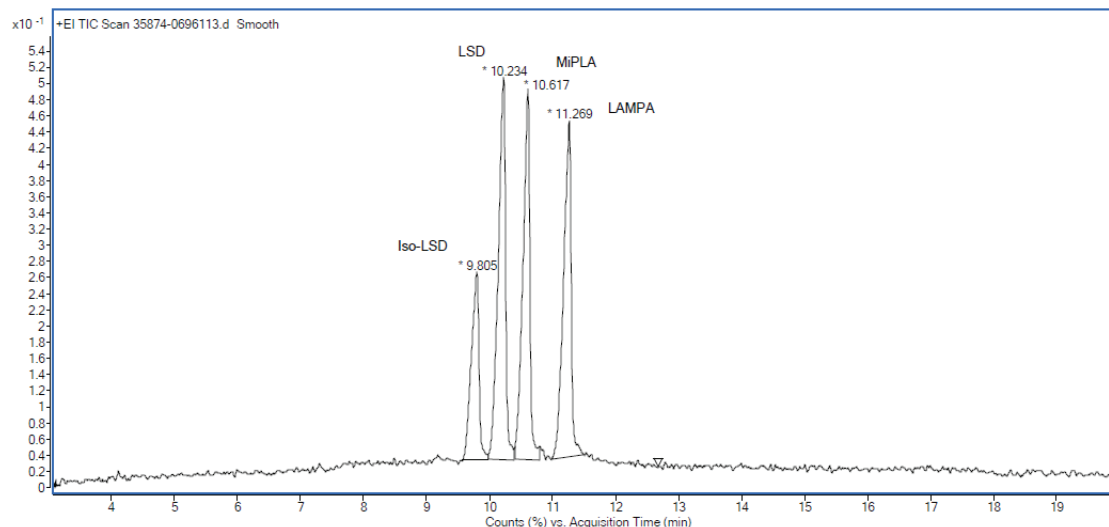
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

GC-MS LSD Isobar Mixture (Item No. 35874) is a mixture of analytical reference materials containing lysergamides. The mixture contains LSD, iso-LSD, MiPLA, and LAMPA. The mixture is supplied in an amber ampule in which the headspace has been purged with argon. It is designed for the detection and positive identification of LSD and its related isobaric compounds, as well as for use in method development and as a reference standard for these substances during routine analytical checks using GC-MS instrumentation. GC-MS LSD Isobar Mixture (Item No. 35874) contains compounds that are regulated as Schedule I compounds in the United States. This product is intended for research and forensic applications.

This mixture contains lysergamides. Included are:

Item No. 35874	GC-MS LSD Isobar Mixture	
Item Name	CAS Number	Concentration (mg/ml)
LSD	50-37-3	1.0
iso-LSD	2126-78-5	1.0
MiPLA	100768-08-9	1.0
LAMPA	40158-98-3	1.0



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, 06/16/2026

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