

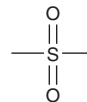
PRODUCT INFORMATION



Dimethyl sulfone

Item No. 20131

CAS Registry No.: 67-71-0
Formal Name: 1,1'-sulfonylbis-methane
Synonyms: DMSO₂, Methylsulfonylmethane, MSM, NSC 63345
MF: C₂H₆O₂S
FW: 94.1
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

Dimethyl sulfone (Item No. 20131) is an analytical reference standard that is structurally categorized as an organic sulfur derivative. It is a natural oxidized form of DMSO and has been described to have anti-inflammatory and analgesic properties.¹ Dimethyl sulfone is commonly found as an adulterant in methamphetamine samples.^{2,3} This product is intended for research and forensic applications.

References

1. Brien, S., Prescott, P., Bashir, N., *et al.* Systematic review of the nutritional supplements dimethyl sulfoxide (DMSO) and methylsulfonylmethane (MSM) in the treatment of osteoarthritis. *Osteoarthritis Cartilage* **16(11)**, 1277-1288 (2008).
2. Cole, C., Jones, L., McVeigh, J., *et al.* CUT: A guide to adulterants, bulking agents and other contaminants found in illicit drugs. (2010).
3. Vu, D.T. SPME/GC-MS characterization of volatiles associated with methamphetamine: Toward the development of a pseudomethamphetamine training material. *J. Forensic Sci.* **46(5)**, 1014-1024 (2001).

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, 06/20/2023

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