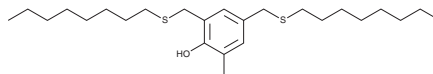


PRODUCT INFORMATION

2-Methyl-4,6-bis(octylsulfanylmethyl)phenol

Item No. 42594

CAS Registry No.: 110553-27-0
 Formal Name: 2-methyl-4,6-bis[(octylthio)methyl]-phenol
 Synonym: BIOP
 MF: $C_{25}H_{44}OS_2$
 FW: 424.7
 Purity: $\geq 95\%$
 Supplied as: A neat oil
 Storage: $-20^{\circ}C$
 Stability: ≥ 4 years



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

Laboratory Procedures

2-Methyl-4,6-bis(octylsulfanylmethyl)phenol (BIOP) is supplied as a neat oil. A stock solution may be made by dissolving the BIOP in the solvent of choice, which should be purged with an inert gas. BIOP is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

BIOP is a phenolic antioxidant.¹ Formulations containing BIOP have been used as stabilizing additives in rubber and styrene synthetic elastomers.

Reference

1. Mohanty, S., Prabhu, P., Kumar, A., *et al.* Degradation and stability studies of butadiene rubber: Correlation between experimental and computational values. *Mater. Today Commun.* **37**, 107558 (2023).

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, 05/01/2025

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