

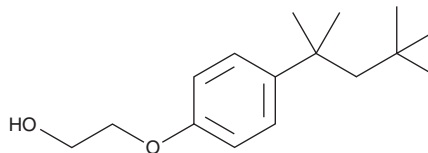
PRODUCT INFORMATION



4-*tert*-Octylphenol monoethoxylate

Item No. 37366

CAS Registry No.: 2315-67-5
Formal Name: 2-[4-(1,1,3,3-tetramethylbutyl)phenoxy]-ethanol
Synonyms: NSC 5259, OP1EO
MF: C₁₆H₂₆O₂
FW: 250.4
Purity: ≥98%
UV/Vis.: λ_{max}: 226 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

4-*tert*-Octylphenol monoethoxylate is supplied as a solid. A stock solution may be made by dissolving the 4-*tert*-octylphenol monoethoxylate in the solvent of choice, which should be purged with an inert gas. 4-*tert*-Octylphenol monoethoxylate is soluble in chloroform and methanol.

Description

4-*tert*-Octylphenol monoethoxylate is an alkylphenoxyethoxylate (APE) and a degradation product of non-ionic surfactants, such as 4-*tert*-octylphenol polyethoxylate.¹ It is also a non-steroidal estrogen.² 4-*tert*-Octylphenol monoethoxylate has been found in wastewater effluent.³

References

1. Nishio, E., Yoshikawa, H., Wakayama, M., *et al.* Isolation and identification of *Sphingomonas* sp. that yields *tert*-octylphenol monoethoxylate under aerobic conditions. *Biosci. Biotechnol. Biochem.* **69(7)**, 1226-1231 (2005).
2. Vajda, A.M., Barber, L.B., Gray, J.L., *et al.* Reproductive disruption in fish downstream from an estrogenic wastewater effluent. *Environ. Sci. Technol.* **42(9)**, 3407-3414 (2008).
3. Barber, L.B., Lee, K.E., Swackhamer, D.L., *et al.* Reproductive responses of male fathead minnows exposed to wastewater treatment plant effluent, effluent treated with XAD8 resin, and an environmentally relevant mixture of alkylphenol compounds. *Aquat. Toxicol.* **82(1)**, 36-46 (2007).

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, 11/09/2022

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