

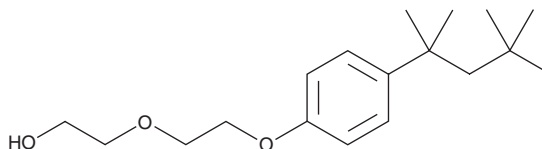
PRODUCT INFORMATION



4-*tert*-Octylphenol diethoxylate

Item No. 18655

CAS Registry No.: 2315-61-9
Formal Name: 2-[2-[4-(1,1,3,3-tetramethylbutyl)phenoxy]ethoxy]-ethanol
Synonym: OP2EO
MF: $C_{18}H_{30}O_3$
FW: 294.4
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 224, 276, 283 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥ 2 years



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

Description

4-*tert*-Octylphenol diethoxylate is a degradation product of the multi-purpose surfactant, Triton X. A known environmental contaminant, 4-*tert*-octylphenol diethoxylate is reported to be an endocrine disruptor in animal and human research studies, producing weak estrogenic effects.¹⁻⁴

References

1. Hanioka, N., Jinno, H., Chung, Y.S., *et al.* Inhibition of rat hepatic cytochrome P450 activities by biodegradation products of 4-*tert*-octylphenol ethoxylate. *Xenobiotica* **29**(9), 873-883 (1999).
2. Nishihara, T., Nishikawa, J., Kanayama, T., *et al.* Estrogenic activities of 517 chemicals by yeast two-hybrid assay. *J. Health Sci.* **46**(4), 282-298 (2000).
3. Schmieder, P.K., Aptula, A.O., Routledge, E.J., *et al.* Estrogenicity of alkylphenolic compounds: A 3-D structure-activity evaluation of gene activation. *Environ. Toxicol. Chem.* **19**(7), 1727-1740 (2000).
4. Colborn, T., vom Saal, F.S., and Soto, A.M. Developmental effects of endocrine-disrupting chemicals in wildlife and humans. *Environ. Health Perspect.* **101**(5), 378-384 (1993).

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, 12/13/2022

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