

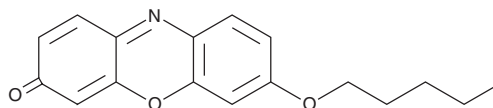
PRODUCT INFORMATION



Pentoxylresorufin

Item No. 14853

CAS Registry No.: 87687-03-4
Formal Name: 7-(pentyloxy)-3H-phenoxazin-3-one
Synonyms: 7-Pentoxylresorufin,
O-Pentylresorufin
MF: C₁₇H₁₇NO₃
FW: 283.3
Purity: ≥95%
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

Pentoxylresorufin is supplied as a solid. A stock solution may be made by dissolving the pentoxylresorufin in the solvent of choice, which should be purged with an inert gas. Pentoxylresorufin is slightly soluble in chloroform.

Description

Pentoxylresorufin is a fluorogenic substrate for cytochrome P450 (CYP) enzymes.^{1,2}

References

1. Lubet, R.A., Mayer, R.T., Cameron, J.W., *et al.* Dealkylation of pentoxylresorufin: A rapid and sensitive assay for measuring induction of cytochrome(s) P-450 by phenobarbital and other xenobiotics in the rat. *Arch. Biochem. Biophys.* **238(1)**, 43-48 (1985).
2. Uehara, S., Murayama, N., Nakanishi, Y., *et al.* Immunochemical detection of cytochrome P450 enzymes in liver microsomes of 27 cynomolgus monkeys. *J. Pharmacol. Exp. Ther.* **339(2)**, 651-661 (2011).

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, 11/08/2022

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