

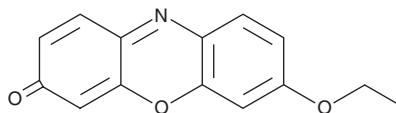
PRODUCT INFORMATION



7-Ethoxyresorufin

Item No. 16122

CAS Registry No.: 5725-91-7
Formal Name: 7-ethoxy-3H-phenoxazin-3-one
Synonyms: 7-ER, 7-Ethoxyphenoxazone, Resorufin ethyl ether
MF: C₁₄H₁₁NO₃
FW: 241.2
Purity: ≥95%
Ex./Em. Max: 570/580 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Special Conditions: Protect from light and moisture.

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

Laboratory Procedures

7-Ethoxyresorufin (7-ER) is supplied as a solid. A stock solution may be made by dissolving the 7-ER in the solvent of choice, which should be purged with an inert gas. 7-ER is soluble in organic solvents such as chloroform, DMSO, and dimethyl formamide.

Description

7-Ethoxyresorufin (7-ER) is a fluorogenic substrate for, and competitive inhibitor of, the cytochrome P450 (CYP) isoform CYP1A1 (IC₅₀ = 0.1 μM).¹⁻³ Upon enzymatic cleavage by CYP1A1, resorufin is released and its fluorescence can be used to quantify CYP1A1 activity. Resorufin displays excitation/emission maxima of 570/580 nm, respectively.⁴ 7-ER has been used in the study of arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) metabolism and the mechanisms of vasodilation.^{5,6} 7-ER is a noncompetitive inhibitor of neuronal nitric oxide synthase (nNOS; K_i = 0.76 μM).^{7,8}

References

1. Dutton, D.R., Reed, G.A., and Parkinson, A. Redox cycling of resorufin catalyzed by rat liver microsomal NADPH-cytochrome P450 reductase. *Arch. Biochem. Biophys.* **268(2)**, 605-616 (1989).
2. Bourrié, M., Meunier, V., Berger, Y., et al. Cytochrome P450 isoform inhibitors as a tool for the investigation of metabolic reactions catalyzed by human liver microsomes. *J. Pharmacol. Exp. Ther.* **277(1)**, 321-332 (1996).
3. Tassaneeyakul, W., Birkett, D.J., Veronese, M.E., et al. Specificity of substrate and inhibitor probes for human cytochromes P450 1A1 and 1A2. *J. Pharmacol. Exp. Ther.* **265(1)**, 401-407 (1993).
4. Hofmann, J. and Sernetz, M. Immobilized enzyme kinetics analyzed by flow-through microfluorimetry: Resorufin-β-D-galactopyranoside as a new fluorogenic substrate for β-galactosidase. *Analytica Chimica Acta* **163**, 67-72 (1984).
5. Labitzke, E.M., Diani-Moore, S., and Rifkind, A.B. Mitochondrial P450-dependent arachidonic acid metabolism by TCDD-induced hepatic CYP1A5; conversion of EETs to DHETs by mitochondrial soluble epoxide hydrolase. *Arch. Biochem. Biophys.* **468(1)**, 70-81 (2007).
6. Imig, J.D., Zou, A.-P., de Montellano, P.R.O., et al. Cytochrome P-450 inhibitors alter afferent arteriolar responses to elevations in pressure. *Am. J. Physiol.* **266(5Pt 2)**, H1879-H1885 (1994).
7. Li, C.G. and Rand, M.J. Inhibition of NO-mediate responses by 7-ethoxyresorufin, a substrate and competitive inhibitor of cytochrome P450. *Br. J. Pharmacol.* **118(1)**, 57-62 (1996).
8. Jiang, H.B. and Ichikawa, Y. Neuronal nitric oxide synthase catalyzes the reduction of 7-ethoxyresorufin. *Life Sci.* **65(12)**, 1257-1264 (1999).

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, 11/28/2022

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