

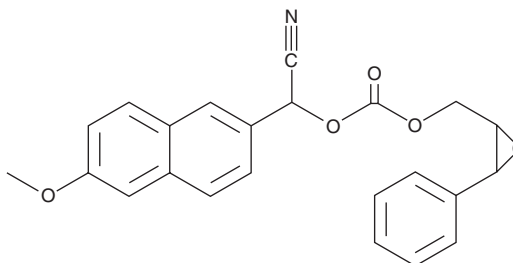
PRODUCT INFORMATION



Epoxy Fluor 7

Item No. 10008610

CAS Registry No.: 863223-43-2
Formal Name: cyano(6-methoxy-2-naphthalenyl)
methyl[(2,3)-3-phenyloxiranyl]
methyl ester, carbonic acid
Synonym: CMNPC
MF: C₂₃H₁₉NO₅
FW: 389.4
Purity: ≥98%
UV/Vis.: λ_{max}: 234 nm
Ex./Em. Max: 330/465 nm
Supplied as: A crystalline 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

Epoxy fluor 7 is supplied as a crystalline solid. A stock solution may be made by dissolving the epoxy fluor 7 in the solvent of choice. Epoxy fluor 7 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of epoxy fluor 7 is approximately 5 mg/ml in ethanol and approximately 10 mg/ml in DMSO and DMF.

Description

Epoxy fluor 7 is a fluorescent substrate for soluble epoxide hydrolase (sEH).^{1,2} Upon hydrolysis by sEH, the resulting diol undergoes cyclization and decomposition to a fluorescent aldehyde product, 6-methoxy-2-naphthaldehyde, which can be used to quantify the activity of purified human or mouse sEH. 6-Methoxy-2-naphthaldehyde displays excitation/emission maxima of 330/465 nm, respectively. Epoxy fluor 7 is intended for use with purified sEH or a preparation free of competing enzyme activities, such as esterase and glutathione S-transferase activity, and can be used for high-throughput screening.²

References

1. Jones, P.D., Wolf, N.M., Morisseau, C., *et al.* Fluorescent substrates for soluble epoxide hydrolase and application to inhibition studies. *Anal. Biochem.* **343**(1), 66-75 (2005).
2. Morisseau, C. and Hammock, B.D. Measurement of soluble epoxide hydrolase (sEH) activity. *Curr. Protoc. Toxicol.* **4.23.2(Suppl. 33)**, (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, 07/03/2024

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