

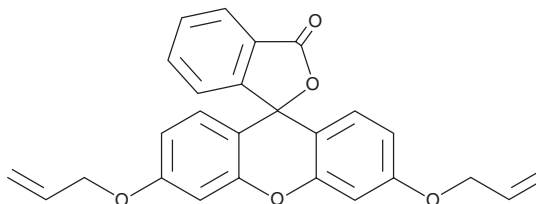
PRODUCT INFORMATION



CAY10732

Item No. 26750

CAS Registry No.: 855751-82-5
Formal Name: 3',6'-bis(2-propen-1-yloxy)-
spiro[isobenzofuran-1(3H),9'-[9H]
xanthen]-3-one
MF: C₂₆H₂₀O₅
FW: 412.4
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 276 nm
Ex./Em. Maxima: 493/527 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

CAY10732 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10732 in the solvent of choice. CAY10732 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of CAY10732 in these solvents is approximately 1, 10, and 12 mg/ml, respectively.

Description

CAY10732 is a fluorescent probe for the detection of carbon monoxide.¹ In the presence of palladium (Pd) and carbon monoxide, CAY10732 undergoes the Pd⁰-mediated Tsuji-Trost reaction to release 2,7-dichlorofluorescein, and its fluorescence can be used to detect carbon monoxide production *in vitro* and in live cells. 2,7-Dichlorofluorescein displays excitation/emission maxima of 493/527 nm, respectively.

Reference

1. Feng, S., Liu, D., Feng, W., *et al.* Allyl fluorescein ethers as promising fluorescent probes for carbon monoxide imaging in living cells. *Anal. Chem.* **89(6)**, 3754-3760 (2017).

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, 02/16/2024

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