

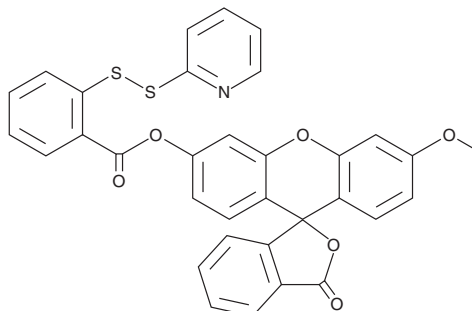
PRODUCT INFORMATION



WSP-1

Item No. 11179

CAS Registry No.: 1352750-34-5
Formal Name: 3'-methoxy-3-oxo-3H-spiro[isobenzofuran-1,9'-xanthen]-6'-yl 2-(pyridin-2-yl)disulfanylbenzoate
Synonym: Washington State Probe-1
MF: C₃₃H₂₁NO₆S₂
FW: 591.7
Purity: ≥95%
UV/Vis.: λ_{max}: 225 nm
Ex./Em. Max: 465/515 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

WSP-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the WSP-1 in the solvent of choice. WSP-1 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of WSP-1 in these solvents is approximately 2 and 20 mg/ml, respectively.

Description

Hydrogen sulfide (H₂S) is an important gaseous mediator, like nitric oxide, that has significant effects on the immunological, neurological, cardiovascular and pulmonary systems of mammals. WSP-1 is a reactive disulfide-containing, fluorescent probe designed to detect H₂S in biological samples and cells. Through a reaction-based fluorescent turn-on strategy, WSP-1 selectively and rapidly reacts with H₂S to generate benzodithiolone and a fluorophore with excitation and emission maxima of 465 and 515 nm, respectively.¹

Reference

1. Liu, C., Pan, J., Li, S., *et al.* Capture and visualization of hydrogen sulfide via a fluorescent probe. *Angew. Chem. Int. Ed. Engl.* **50(44)**, 10327-10329 (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, 01/11/2018

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