

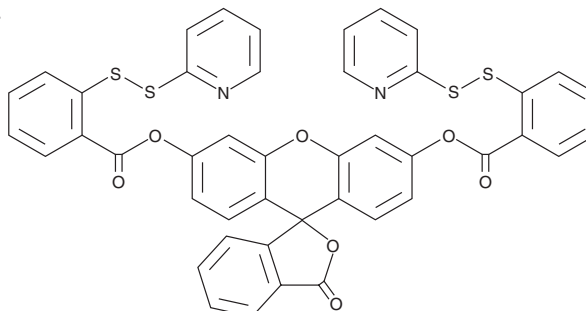
# PRODUCT INFORMATION



## WSP-5

Item No. 16929

**CAS Registry No.:** 1593024-78-2  
**Formal Name:** 3-oxo-3H-spiro[isobenzofuran-1,9'-xanthene]-3',6'-diyl bis(2-(pyridin-2-yl)disulfanyl)benzoate)  
**Synonym:** Washington State Probe-5  
**MF:** C<sub>44</sub>H<sub>26</sub>N<sub>2</sub>O<sub>7</sub>S<sub>4</sub>  
**FW:** 822.9  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 225 nm  
**Ex./Em. Max:** 502/525 nm, respectively  
**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-5 is supplied as a crystalline solid. A stock solution may be made by dissolving the WSP-5 in the solvent of choice. WSP-5 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of WSP-5 in these solvents is approximately 25 and 30 mg/ml.

### Description

WSP-5 is a turn-on fluorescent probe for hydrogen sulfide (H<sub>2</sub>S).<sup>1</sup> Upon reaction with H<sub>2</sub>S, WSP-5 releases a fluorophore that displays excitation/emission maxima of 502/525 nm, respectively. WSP-5 exhibits a faster turn-on rate and a more sensitive detection limit toward H<sub>2</sub>S than WSP-1 (Item No. 11179).

### Reference

1. Peng, B., Chen, W., Liu, C., *et al.* Fluorescent probes based on nucleophilic substitution-cyclization for hydrogen sulfide detection and bioimaging. *Chemistry* **20(4)**, 1010-1016 (2014).

#### 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, 07/09/2020

#### 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