

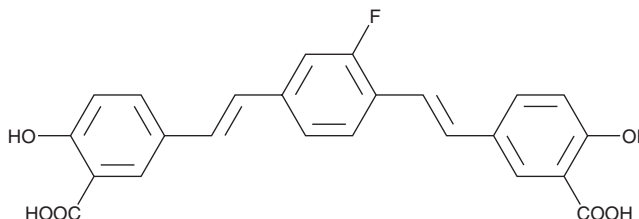
PRODUCT INFORMATION



FSB

Item No. 10651

CAS Registry No.: 760988-03-2
Formal Name: 3,3'-[(2-fluoro-1,4-phenylene)di-(1E)-2,1-ethenediyl]bis-6-hydroxy-benzoic acid
MF: C₂₄H₁₇FO₆
FW: 420.4
Purity: ≥95%
Ex./Em. Max: 390/520 nm
Supplied as: A 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

FSB is supplied as a solid. A stock solution may be made by dissolving the FSB in the solvent of choice, which should be purged with an inert gas. FSB is soluble in the organic solvent DMSO at a concentration of approximately 200 mg/ml.

Description

FSB is a fluorescent probe that has been used for the detection of amyloid plaques in human postmortem brain from patients with Alzheimer's disease and of curly amyloid fibers in *E. coli* biofilms.¹ It has also been used to label tau inclusions in mouse and human brain sections and in mouse spinal cord following intravenous injection.² FSB has excitation/emission maxima of 390/520 nm, respectively.

References

1. Sato, K., Higuchi, M., Iwata, N., *et al.* Fluoro-substituted and ¹³C-labeled styrylbenzene derivatives for detecting brain amyloid plaques. *Eur. J. Med. Chem.* **39(7)**, 573-578 (2004).
2. Velasco, A., Fraser, G., Delobel, P., *et al.* Detection of filamentous tau inclusions by the fluorescent Congo red derivative FSB [(*trans,trans*)-1-fluoro-2,5-bis(3-hydroxycarbonyl-4-hydroxy)styrylbenzene]. *FEBS Lett.* **582(6)**, 901-906 (2008).

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, 10/06/2022

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