

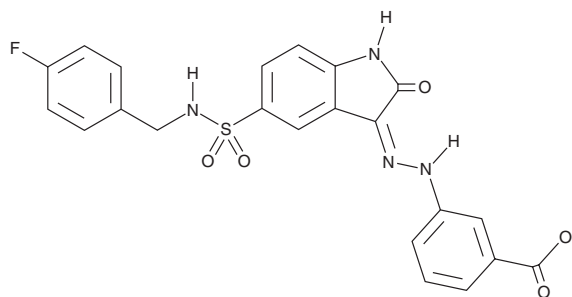
PRODUCT INFORMATION



SPI 112

Item No. 29474

CAS Registry No.: 1051387-90-6
Formal Name: 3-[[[2Z]-2-[5-[[[(4-fluorophenyl)methyl]amino]sulfonyl]-1,2-dihydro-2-oxo-3H-indol-3-ylidene]hydrazinyl]-benzoic acid
MF: C₂₂H₁₇FN₄O₅S
FW: 468.5
Purity: ≥95%
UV/Vis.: λ_{max}: 216, 408 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

SPI 112 is supplied as a crystalline solid. A stock solution may be made by dissolving the SPI 112 in the solvent of choice, which should be purged with an inert gas. SPI 112 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of SPI 112 in these solvents is approximately 2 and 1 mg/ml, respectively.

SPI 112 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SPI 112 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SPI 112 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SPI 112 is cell-impermeable inhibitor of Src homology region 2 domain-containing phosphatase 2 (SHP-2) in cell-free assays (IC₅₀ = 1 μM).¹ It selectively inhibits SHP-2 over SHP-1 and protein tyrosine phosphatase 1B (PTP1B; IC₅₀s = 18.3 and 14.5 μM, respectively).

Reference

1. Lawrence, H.R., Pireddu, R., Chen, L., *et al.* Inhibitors of Src homology-2 domain containing protein tyrosine phosphatase-2 (Shp2) based on oxindole scaffolds. *J. Med. Chem.* **51**(16), 4548-4956 (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, 12/02/2024

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