

PRODUCT INFORMATION



SPI 112Me

Item No. 36819

CAS Registry No.: 1243685-62-2

Formal Name: 3-[(2Z)-2-[5-[[[4-fluorophenyl)methyl]amino]sulfonyl]-1,2-dihydro-2-oxo-3H-indol-3-ylidene]hydrazinyl]-benzoic acid, methyl ester

MF: $C_{23}H_{19}FN_4O_5S$

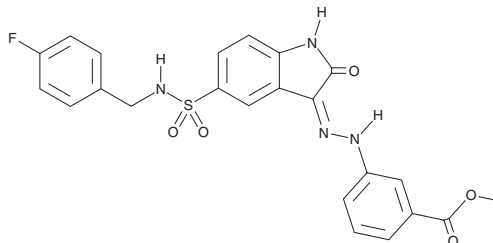
FW: 482.5

Purity: $\geq 90\%$

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

SPI 112Me is supplied as a solid. A stock solution may be made by dissolving the SPI 112Me in the solvent of choice, which should be purged with an inert gas. SPI 112Me is soluble (≥ 10 mg/ml) in acetonitrile and DMSO.

Description

SPI 112Me is a prodrug form of the Src homology region 2 domain-containing phosphatase 2 (SHP-2) inhibitor SPI 112 (Item No. 29474).¹ It is hydrolyzed by esterases intracellularly to form SPI 112. SPI 112Me inhibits the protein tyrosine phosphatase (PTP) activity of SHP-2 containing a constitutively activating glutamate-to-lysine substitution at position 76 (SHP-2^{E76K}) in TF-1 leukemia cells expressing the mutant enzyme. SPI 112Me (20 μM) inhibits Paxillin dephosphorylation, ERK1/2 activation, and cell migration in EGF-stimulated MDA-MB-468 breast cancer cells.

Reference

1. Chen, L., Pernazza, D., Scott, L.M., *et al.* Inhibition of cellular Shp2 activity by a methyl ester analog of SPI-112. *Biochem. Pharmacol.* **80**(6), 801-810 (2010).

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, 03/05/2026

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