

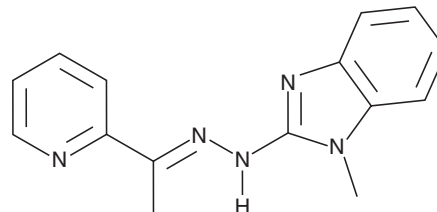
PRODUCT INFORMATION



SI-2

Item No. 19565

CAS Registry No.: 380537-35-9
Formal Name: (1E)-1-(2-pyridinyl)-ethanone-2-(1-methyl-1H-benzimidazol-2-yl)hydrazone
Synonym: EPH 116
MF: C₁₅H₁₅N₅
FW: 265.3
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 369 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

SI-2 is supplied as a crystalline solid. A stock solution may be made by dissolving the SI-2 in the solvent of choice, which should be purged with an inert gas. SI-2 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SI-2 in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

SI-2 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SI-2 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SI-2 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

SI-2 is an inhibitor of steroid receptor coactivator 3 (SRC-3). It has been shown to selectively reduce the transcriptional activities and the protein concentrations of SRC-3 in cells through direct physical interactions with SRC-3.¹ SI-2 induces death in various breast cancer cells with IC₅₀ values of 3-20 nM without affecting normal cell viability.¹ At 2 mg/kg, SI-2 can also inhibit primary tumor growth and reduce SRC-3 protein levels in an MDA-MB-468 breast cancer mouse model.

Reference

1. Song, X., Chen, J., Zhao, M., *et al.* Development of potent small-molecule inhibitors to drug the undruggable steroid receptor coactivator-3. *Proc. Natl. Acad. Sci. USA* **113(18)**, 4970-4975 (2016).

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, 02/22/2023

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