

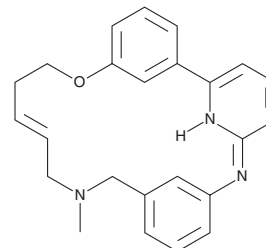
PRODUCT INFORMATION



SB-1317

Item No. 26817

CAS Registry No.: 937270-47-8
Formal Name: 14-methyl-20-oxa-5,7,14,27-tetraazatetracyclo[19.3.1.12,6.18,12]heptacos-1(25),2,4,6(27),8,10,12(26),16,21,23-decaene
Synonyms: TG02, Zotiraciclib
MF: C₂₃H₂₄N₄O
FW: 372.5
Purity: ≥98%
UV/Vis.: λ_{max}: 282 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

SB-1317 is supplied as a solid. A stock solution may be made by dissolving the SB-1317 in the solvent of choice, which should be purged with an inert gas. SB-1317 is soluble in DMSO at a concentration of approximately 30 mg/ml. SB-1317 is slightly soluble in dimethyl formamide.

Description

SB-1317 is a multi-kinase inhibitor.¹ It inhibits Cdk1, -2, -3, -5, and -9 (IC₅₀s = 3-9 nM), as well as JAK family kinases (IC₅₀s = 14-59 nM), wild-type and mutant FLT3s (IC₅₀s = 19-27 nM), and Src family kinases (IC₅₀s = 11-15 nM). SB-1317 inhibits proliferation in a panel of 12 liquid and six solid tumor cell lines (mean IC₅₀ = 0.19 μM). It induces cell cycle arrest at the G₁ phase and apoptosis in MV4-11 acute myeloid leukemia (AML) cells. SB-1317 (10, 20, and 40 mg/kg) reduces tumor volume and increases survival in an MV4-11 mouse xenograft model.

Reference

1. Goh, K.C., Novotny-Diermayr, V., Hart, S., et al. TG02, a novel oral multi-kinase inhibitor of CDKs, JAK2 and FLT3 with potent anti-leukemic properties. *Leukemia* **26(2)**, 236-243 (2012).

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, 11/18/2022

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