

PRODUCT INFORMATION

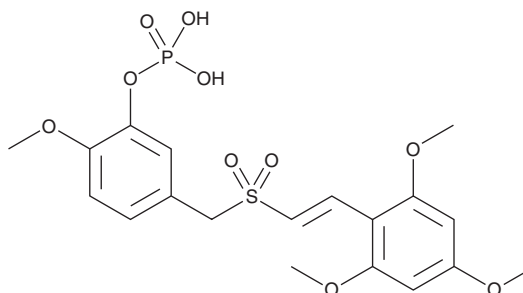


Briciclib

Item No. 22989

CAS Registry No.: 865783-99-9
Formal Name: 2-methoxy-5-[[[(1E)-2-(2,4,6-trimethoxyphenyl)ethenyl]sulfonyl]methyl]-phenol 1-(dihydrogen phosphate)

Synonym: ON-014185
MF: C₁₉H₂₃O₁₀PS
FW: 474.4
Purity: ≥98%
UV/Vis.: λ_{max}: 236, 308 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

Briciclib is supplied as a crystalline solid. A stock solution may be made by dissolving the briciclib in the solvent of choice. Briciclib is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of briciclib in these solvents is approximately 30 mg/ml.

Description

Briciclib is a water soluble derivative of ON 013100, a small molecule that binds to and inhibits eukaryotic translation initiation factor 4E (eIF4E).¹ Briciclib inhibits proliferation of JEKO-1 and Mino mantle cell leukemia (MCL), MCF-7 and MDA-MB-231 breast, AGS gastric, and OE19, OE33, and LFO-1 esophageal cancer cell lines with GI₅₀ values ranging from 9.8 to 12.2 nM. It reduces expression of cyclin D1 and c-Myc, markers of eIF4E activity, in breast and MCL cancer cell lines in a dose-dependent manner. Briciclib also induces expression of the pro-apoptotic proteins p53 and cleaved caspase-3.

Reference

1. Jasani, N., Desai, B.J., Betzu, J.M., *et al.* Abstract 1649: Potent anticancer activity of an orally bioavailable small molecule, ON 013100, and its water soluble derivative, briciclib, a clinical-stage eIF4E-targeted agent. *Cancer Res.* **75(15 Suppl)**, 1649 (2015).

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, 01/03/2023

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