

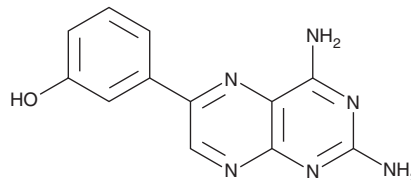
PRODUCT INFORMATION



TG100713

Item No. 30195

CAS Registry No.: 925705-73-3
Formal Name: 3-(2,4-diamino-6-pteridiny)-phenol
MF: C₁₂H₁₀N₆O
FW: 254.3
Purity: ≥98%
UV/Vis.: λ_{max}: 277, 384 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

TG100713 is supplied as a crystalline solid. A stock solution may be made by dissolving the TG100713 in the solvent of choice, which should be purged with an inert gas. TG100713 is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

TG100713 is an inhibitor of PI3K (IC₅₀s = 50, 24, 165, and 215 nM for PI3Kγ, PI3Kδ, PI3Kα, and PI3Kβ, respectively).¹ It inhibits proliferation of human umbilical vein endothelial cells (HUVECs) when used at a concentration of 10 μM. TG100713 (2 μM) inhibits reactivation of latent HIV-1 induced by panobinostat (Item No. 13280), but not prostratin (Item No. 10272) or JQ-1, in 24ST1NLESG cells.²

References

1. Doukas, J., Wrasidlo, W., Noronha, G., *et al.* Phosphoinositide 3-kinase γ/δ inhibition limits infarct size after myocardial ischemia/reperfusion injury. *Proc. Natl. Acad. Sci. USA* **103**(52), 19866-19871 (2006).
2. Vargas, B., Giacobbi, N.S., Sanyal, A., *et al.* Inhibitors of signaling pathways that block reversal of HIV-1 latency. *Antimicrob. Agents Chemother.* **63**(2), e01744-18 (2019).

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, 05/28/2020

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