

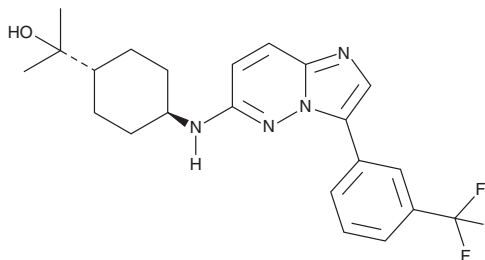
PRODUCT INFORMATION



TP-3654

Item No. 29571

CAS Registry No.: 1361951-15-6
Formal Name: *trans*- α,α -dimethyl-4-[[3-(trifluoromethyl)phenyl]imidazo[1,2-*b*]pyridazin-6-yl]amino]-cyclohexanemethanol
MF: C₂₂H₂₅F₃N₄O
FW: 418.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 217, 274, 342 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

TP-3654 is supplied as a crystalline solid. A stock solution may be made by dissolving the TP-3654 in the solvent of choice, which should be purged with an inert gas. TP-3654 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TP-3654 in ethanol is approximately 10 mg/ml and approximately 16 mg/ml in DMSO and DMF.

TP-3654 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TP-3654 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. TP-3654 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

TP-3654 is an inhibitor of Pim kinases (K_{is} = 5, 239, and 42 nM for Pim-1, Pim-2, and Pim-3, respectively).¹ It inhibits 19 additional kinases, including PI3K α , PI3K γ , and PI3K δ , with IC₅₀ values of less than 300 nM. TP-3654 inhibits phosphorylation of Bad in HEK293 cells overexpressing Bad and Pim-1 (EC₅₀ = 67 nM), as well as in UM-UC-3 cells. It reduces tumor growth in UM-UC-3 bladder carcinoma and PC3 prostate adenocarcinoma mouse xenograft models when administered at a dose of 200 mg/kg.

Reference

1. Foulks, J.M., Carpenter, K.J., Luo, B., *et al.* A small-molecule inhibitor of PIM kinases as a potential treatment for urothelial carcinomas. *Neoplasia* **16**(5), 403-412 (2014).

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, 10/17/2022

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