

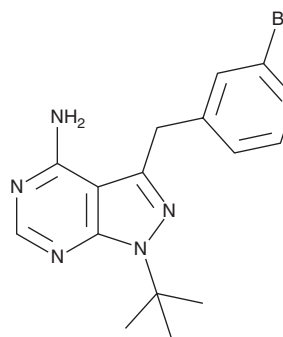
PRODUCT INFORMATION



3BrB-PP1

Item No. 45582

CAS Registry No.: 956025-99-3
Formal Name: 3-[(3-bromophenyl)methyl]1-(1,1-dimethylethyl)-1H-pyrazolo[3,4-d]pyrimidin-4-amine
MF: C₁₆H₁₈BrN₅
FW: 360.3
Purity: ≥98%
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

3BrB-PP1 is supplied as a solid. A stock solution may be made by dissolving the 3BrB-PP1 in the solvent of choice, which should be purged with an inert gas. 3BrB-PP1 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

3BrB-PP1 is a bumped kinase inhibitor (BKI) that acts as a selective, ATP-competitive inhibitor of kinases with gatekeeper mutations.¹ It increases the activity of the *T. gondii* bradyzoite-specific BAG1 promoter. 3BrB-PP1 has been used to study kinase activity using kinases containing mutations at gatekeeper residues in the ATP-binding pocket, thereby allowing 3BrB-PP1 to selectively inhibit the kinase.^{2,3}

References

1. Sugia, T., Masatania, T., Murakoshia, F., *et al.* Microplate assay for screening *Toxoplasma gondii* bradyzoite differentiation with DUAL luciferase assay. *Anal. Biochem.* **464**, 9-11 (2014).
2. Mutavchiev, D.R., Leda, M., and Sawin, K.E. Remodeling of the fission yeast Cdc42 cell-polarity module via the Sty1 p38 stress-activated protein kinase pathway. *Curr. Biol.* **26(21)**, 2921-2928 (2016).
3. Tay, Y.D., Leda, M., Spanos, C., *et al.* Fission yeast NDR/LATS kinase Orb6 regulates exocytosis via phosphorylation of the exocyst complex. *Cell Rep.* **26(6)**, 1654-1667 (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, 05/06/2026

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