

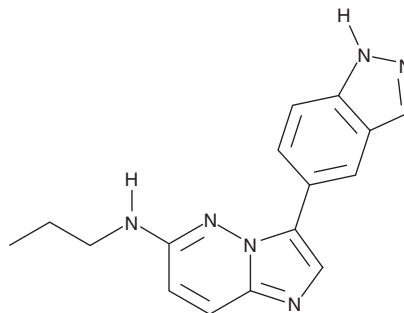
PRODUCT INFORMATION



CHR-6494

Item No. 11478

CAS Registry No.: 1333377-65-3
Formal Name: 3-(1H-indazol-5-yl)-N-propyl-imidazo[1,2-b]pyridazin-6-amine
MF: C₁₆H₁₆N₆
FW: 292.3
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 252, 335 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

CHR-6494 is supplied as a crystalline solid. A stock solution may be made by dissolving the CHR-6494 in the solvent of choice. CHR-6494 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of CHR-6494 in these solvents is approximately 0.25, 20, and 12.5 mg/ml, respectively.

CHR-6494 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CHR-6494 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CHR-6494 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CHR-6494 is a haspin protein kinase inhibitor (IC₅₀ = 2 nM) selective over 27 other protein kinases, including Aurora B kinase (IC₅₀ > 1 μM). Haspin is a kinase necessary for mitosis that selectively phosphorylates histone 3 at the threonine 3 residue (H3T3ph).¹ CHR-6494 decreases phosphorylation of H3T3 and arrests the cell cycle in the G₂/M phase, which leads to apoptosis. It showed promising results in a xenograft tumor model.

Reference

1. Huertas, D., Soler, M., Moreto, J., *et al.* Antitumor activity of a small-molecule inhibitor of the histone kinase Haspin. *Oncogene*, 1-11 (2011).

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