

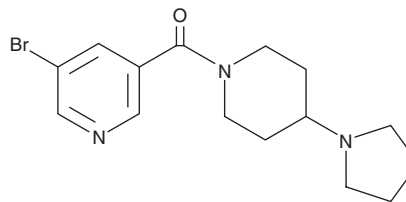
PRODUCT INFORMATION



UNC669

Item No. 10875

CAS Registry No.: 1314241-44-5
Formal Name: (5-bromo-3-pyridinyl)[4-(1-pyrrolidinyl)-1-piperidinyl]-methanone
MF: C₁₅H₂₀BrN₃O
FW: 338.2
Purity: ≥98%
UV/Vis.: λ_{max}: 204, 274 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

UNC669 is supplied as a crystalline solid. A stock solution may be made by dissolving the UNC669 in the solvent of choice. UNC669 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of UNC669 in these solvents is approximately 10, 0.5, and 20 mg/ml, respectively.

UNC669 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, UNC669 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. UNC669 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

UNC669 is an inhibitor of the MBT domain of L3MBTL1 with an IC₅₀ value of 6 μM.¹ It is 5- and 10-fold selective for L3MBTL1 over the related proteins L3MBTL3 and L3MBTL4.

Reference

1. Herold, J.M., Wigle, T.J., Norris, J.L., *et al.* Small-molecule ligands of methyl-lysine binding proteins. *J. Med. Chem.* **54**, 2504-11 (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, 11/08/2022

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