

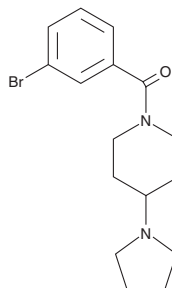
PRODUCT INFORMATION



UNC926

Item No. 27302

CAS Registry No.: 1184136-10-4
Formal Name: (3-bromophenyl)[4-(1-pyrrolidinyl)-1-piperidinyl]-methanone
MF: C₁₆H₂₁BrN₂O
FW: 337.3
Purity: ≥98%
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.

Description

UNC926 is an antagonist of the MBT domain-containing methyl-lysine reader proteins L3MBTL1, L3MBTL3, and L3MBTL4.¹ It inhibits methyl-lysine histone peptide substrate binding of L3MBTL1, L3MBTL3, and L3MBTL4 (IC₅₀s = 3.9, 3.2, and 15.6 μM, respectively) selectively over MBT domain-containing MBTD1 and SFMBT1 and chromo domain-containing CBX7 (IC₅₀s = >30 μM).

Reference

1. Herold, J.M., James, L.I., Korboukh, V.K., *et al.* Structure-activity relationships of methyl-lysine reader antagonists. *Med. Chem. Commun.* **3**(1), 45-51 (2012).

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, 11/30/2022

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