

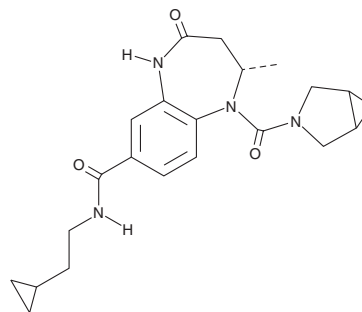
PRODUCT INFORMATION



BAY-6035

Item No. 25925

CAS Registry No.: 2247890-13-5
Formal Name: (2S)-1-(3-azabicyclo[3.1.0]hex-3-ylcarbonyl)-N-(2-cyclopropylethyl)-2,3,4,5-tetrahydro-2-methyl-4-oxo-1H-1,5-benzodiazepine-7-carboxamide
MF: C₂₂H₂₈N₄O₃
FW: 396.5
Purity: ≥95%
UV/Vis.: λ_{max}: 273 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

BAY-6035 is supplied as a crystalline solid. A stock solution may be made by dissolving the BAY-6035 in the solvent of choice, which should be purged with an inert gas. BAY-6035 is soluble in DMSO.

Description

BAY-6035 is an inhibitor of SET and MYND domain-containing protein 3 (SMYD3). It is 100-fold selective for SMYD3 over other histone methyltransferases. BAY-6035 inhibits methylation of MEKK2 peptide in a cell-free assay and in cells (IC₅₀s = 88 and 70 nM, respectively). See the Structural Genomic Consortium (SGC) website for more information.

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, 12/19/2022

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