

PRODUCT INFORMATION



AT-1

Item No. 33265

CAS Registry No.: 2098836-45-2
Formal Name: (4R)-N-acetyl-3-[[6-[[2-[[[(6S)-4-(4-chlorophenyl)-2,3,9-trimethyl-6H-thieno[3,2-f][1,2,4]triazolo[4,3-a][1,4]diazepin-6-yl]acetyl]amino]hexyl]thio]-L-valyl-4-hydroxy-N-[[4-(4-methyl-5-thiazolyl)phenyl]methyl]-L-prolinamide

MF: C₄₈H₅₈ClN₉O₅S₃

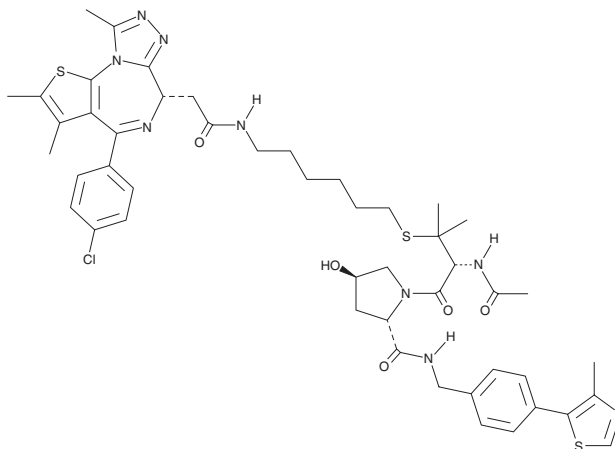
FW: 972.7

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

AT-1 is supplied as a solid. A stock solution may be made by dissolving the AT-1 in the solvent of choice, which should be purged with an inert gas. AT-1 is slightly soluble (0.1-1 mg/ml) in methanol and chloroform.

Description

AT-1 is a proteolysis-targeting chimera (PROTAC) that contains the bromodomain and extra terminal (BET) domain-containing protein inhibitor (+)-JQ1 (Item No. 11187) conjugated to the PROTAC building block VH ligand O32 (Item No. 39982).¹ It selectively degrades bromodomain-containing protein 4 (BRD4) over BRD3 and BRD2 in HeLa cells at 300 nM.

Reference

1. Löscher, W., Hönack, D., Nolting, B., *et al.* Trans-2-en-valproate: Reevaluation of its anticonvulsant efficacy in standardized seizure models in mice, rats and dogs. *Epilepsy Res.* **9(3)**, 195-210 (1991).

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, 06/14/2024

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