

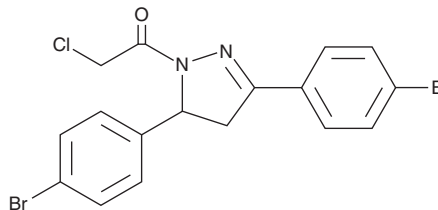
PRODUCT INFORMATION



EN219

Item No. 33957

CAS Registry No.: 380351-29-1
Formal Name: 1-[3,5-bis(4-bromophenyl)-4,5-dihydro-1H-pyrazol-1-yl]-2-chloro-ethanone
MF: C₁₇H₁₃Br₂ClN₂O
FW: 456.6
Purity: ≥95%
UV/Vis.: λ_{max}: 224, 295, 305 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

EN219 is supplied as a crystalline solid. A stock solution may be made by dissolving the EN219 in the solvent of choice, which should be purged with an inert gas. EN219 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of EN219 in these solvents is approximately 1, 5, and 10 mg/ml, respectively.

Description

EN219 is a synthetic recruiter of the E3 ubiquitin ligase RNF114.¹ It binds to cysteine 8 (C8) in the intrinsically disordered region of RNF114 (RNF114-C8; IC₅₀ = 470 nM) and inhibits RNF114-induced autoubiquitination and p21 ubiquitination in a cell-free assay when used at a concentration of 50 μM. EN219 (1 μM) also interacts with cysteine residues in the tubulin β1 chain (TUBB1), heat shock protein 60 (Hsp60), also known as Hsp family D member 1 (HspD1), and histone H3.1 (HIST1H3A) in 231MFP human breast cancer cells in a proteomic profiling assay. It has been linked to the bromodomain and extra terminal domain (BET) inhibitor ligand (+)-JQ1 for use as a proteolysis-targeting chimera (PROTAC) to degrade bromodomain-containing protein 4 (BRD4) in 231MFP cells.

Reference

1. Luo, M., Spradlin, J.N., Boike, L., *et al.* Chemoproteomics-enabled discovery of covalent RNF114-based degraders that mimic natural product function. *Cell Chem. Biol.* **28(4)**, 559-566 (2021).

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, 03/26/2024

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