

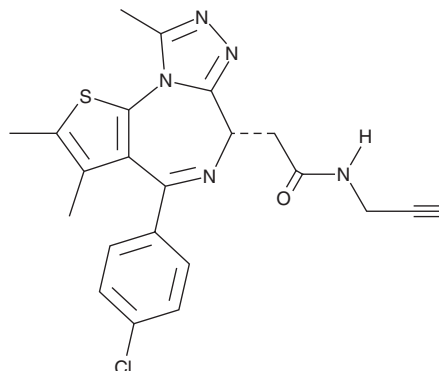
PRODUCT INFORMATION



(+)-JQ1 PA

Item No. 30414

CAS Registry No.: 2115701-93-2
Formal Name: (6S)-4-(4-chlorophenyl)-2,3,9-trimethyl-N-2-propyn-1-yl-6H-thieno[3,2-f][1,2,4]triazolo[4,3-a][1,4]diazepine-6-acetamide
MF: C₂₂H₂₀ClN₅OS
FW: 437.9
Purity: ≥98%
UV/Vis.: λ_{max}: 255 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

(+)-JQ1 PA is supplied as a crystalline solid. A stock solution may be made by dissolving the (+)-JQ1 PA in the solvent of choice, which should be purged with an inert gas. (+)-JQ1 PA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (+)-JQ1 PA in ethanol is approximately 25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

(+)-JQ1 PA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (+)-JQ1 PA should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. (+)-JQ1 PA has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(+)-JQ1 PA is a clickable form of the bromodomain and extraterminal domain (BET) inhibitor (+)-JQ1 (Item Nos. 11187 | 9002910).¹ (+)-JQ1 PA binds to bromodomain-containing protein 4 (BRD4) and displaces it from *Myc* in a ChIP-qPCR assay. It reduces proliferation of MV4-11 cells (IC₅₀ = 10.4 nM).

Reference

1. Tyler, D.S., Vappiani, J., Caneque, T., *et al.* Click chemistry enables preclinical evaluation of targeted epigenetic therapies. *Science* **356**(6345), 1397-1401 (2017).

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, 05/15/2020

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