

PRODUCT INFORMATION



A-395 (hydrochloride)

Item No. 20257

Formal Name: (3R,4S)-1-(7-fluoro-2,3-dihydro-1H-inden-1-yl)-N,N-dimethyl-4-(4-(4-(methylsulfonyl)piperazin-1-yl)phenyl)pyrrolidin-3-amine, hydrochloride

MF: C₂₆H₃₅FN₄O₂S • XHCl

FW: 486.6

Purity: ≥98%

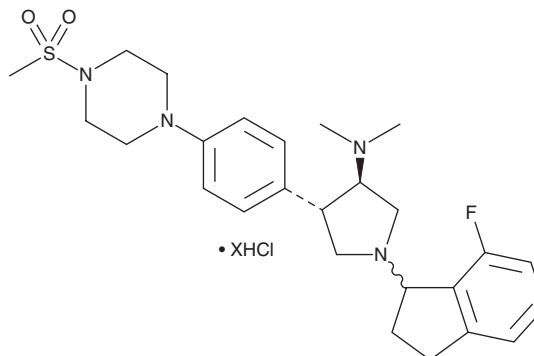
UV/Vis.: λ_{max}: 254 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years

Special Conditions: Hygroscopic



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

Laboratory Procedures

A-395 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the A-395 (hydrochloride) in water. We do not recommend storing the aqueous solution for more than one day.

Description

A-395 is an inhibitor of the interaction between embryonic ectoderm development (EED) and histone H3 lysine 27 (H3K27) that binds to the EED subunit of polycomb repressor complex 2 (PRC2) and inhibits H3K27 methylation ($K_i = 0.4$ nM).¹ It inhibits H3K27 di- and trimethylation in RD rhabdoid tumor cells (IC_{50} s = 390 and 90 nM, respectively). *In vivo*, A-395 (200 mg/kg) reduces tumor volume and intratumor H3K27 trimethylation in a Pfeiffer diffuse large B cell lymphoma (DLBCL) mouse xenograft model.

Reference

1. Hey, Y., Selvaraju, S., Curtin, M.L., *et al.* The EED protein-protein interaction inhibitor A-395 inactivates the PRC2 complex. *Nat. Chem. Biol.* **13**(4), 389-395 (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, 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