

# 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-(methylsulfonyl)piperazin-1-yl)phenyl pyrrolidin-3-amine, hydrochloride

**MF:**  $C_{26}H_{35}FN_4O_2S \cdot XHCl$

**FW:** 486.6

**Purity:**  $\geq 98\%$

**UV/Vis.:**  $\lambda_{max}$ : 254 nm

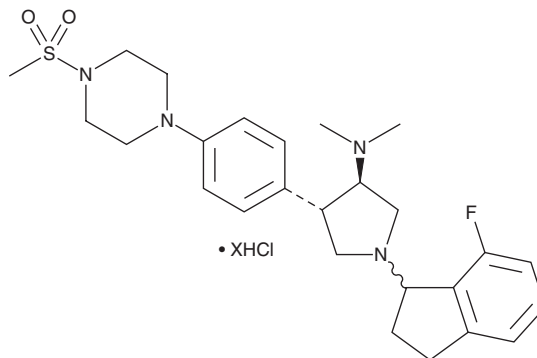
**Supplied as:** A crystalline solid

**Storage:**  $-20^{\circ}C$

**Stability:**  $\geq 4$  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).<sup>1</sup> 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. See the Structural Genomics Consortium (SGC) website for more information.

### 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, 08/05/2024

#### 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