

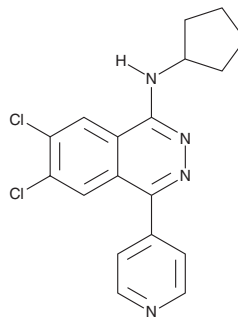
PRODUCT INFORMATION



A-196

Item No. 18317

CAS Registry No.: 1982372-88-2
Formal Name: 6,7-dichloro-N-cyclopentyl-4-(4-pyridinyl)-1-phthalazinamine
MF: C₁₈H₁₆Cl₂N₄
FW: 359.3
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 228, 273, 351 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

A-196 is supplied as a crystalline solid. A stock solution may be made by dissolving the A-196 in the solvent of choice, which should be purged with an inert gas. A-196 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of A-196 in ethanol is approximately 2 mg/ml and approximately 10 mg/ml in DMSO and DMF.

A-196 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, A-196 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. A-196 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

A-196 is a selective inhibitor of SUV420H1 and SUV420H2 (IC₅₀s = 25 and 144 nM, respectively) that is more than 100-fold selective over other histone methyltransferases and non-epigenetic targets. A-196 has been shown to inhibit the di- and tri-methylation of lysine 20 of histone H4 (H4K20me) in multiple cell lines with IC₅₀ values less than 1 μM. For more information on A-196 please visit the Structural Genomics Consortium (SGC). The negative control, SGC2043 (A-197), for A-196 is also available exclusively through the SGC.

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