

PRODUCT INFORMATION



L-Moses (hydrochloride)

Item No. 21288

Formal Name: (1S,2S)-N¹,N¹-dimethyl-N²-(3-methyl-1,2,4-triazolo[3,4-a]phthalazin-6-yl)-1-phenyl-1,2-propanediamine, dihydrochloride

Synonym: L-45

MF: C₂₁H₂₄N₆ • 2HCl

FW: 433.4

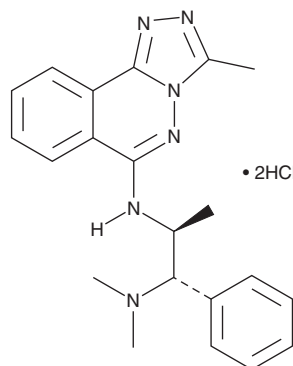
Purity: ≥98%

UV/Vis.: λ_{max}: 233, 241, 557 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

L-Moses (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the L-moses (hydrochloride) in the solvent of choice, which should be purged with an inert gas. L-Moses (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of L-moses (hydrochloride) in these solvents is approximately 10, 20, and 25 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of L-moses (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of L-moses (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

L-Moses is an inhibitor of the p300/CBP-associated factor (PCAF) bromodomain ($K_d = 126$ nM).¹ It is selective for PCAF over GCN5 ($K_d = 600$ nM) and >4,500-fold selective for PCAF over BRD4. L-Moses displaces the PCAF bromodomain from histone H3.3 in a nanoBRET target engagement assay when used at a concentration of 5 μM.

Reference

1. Moustakim, M., Clark, P.G.K., Trulli, L., *et al.* Discovery of a PCAF bromodomain chemical probe. *Angew Chem. Int. Ed. Engl.* **56**(3), 827-831 (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, 11/07/2022

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