

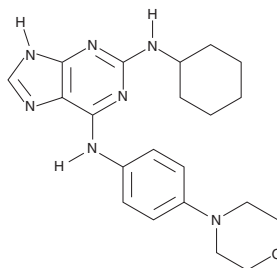
PRODUCT INFORMATION



Retreversine

Item No. 10006606

CAS Registry No.: 1028332-91-3
Formal Name: N²-cyclohexyl-N⁶-[4-(4-morpholinyl)phenyl]-1H-purine-2,6-diamin
MF: C₂₁H₂₇N₇O
FW: 393.5
Purity: ≥98%
UV/Vis.: λ_{max}: 263, 318 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

Retreversine is supplied as a crystalline solid. A stock solution may be made by dissolving the retreversine in the solvent of choice, which should be purged with an inert gas. Retreversine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of retreversine in these solvents is approximately 2.5 and 1.6 mg/ml, respectively.

Retreversine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, retreversine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Retreversine has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Terminal differentiation of embryonic progenitor cells to form distinct, adult tissues is the hallmark of organogenesis in complex organisms. However, in mammals including humans, this process is not readily reversible. Reversine is a 2,6-disubstituted purine derivative that was identified using a screening system based on the loss of terminal differentiation markers in C2C12 myoblast cells.¹ Retreversine is an inactive control molecule, created by simply interchanging the 2 and 6-substituents on the purine ring. Retreversine is inactive at 10-fold higher concentrations than reversine in the myoblast assay.¹

Reference

1. Chen, S., Zhang, Q., Wu, X., *et al.* Dedifferentiation of lineage-committed cells by a small molecule. *J. Am. Chem. Soc.* **126(2)**, 410-411 (2004).

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, 11/09/2023

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