

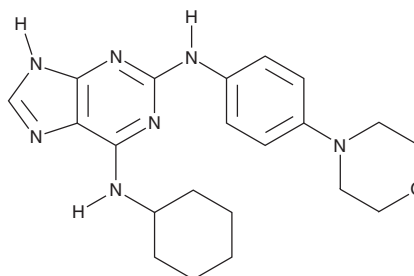
PRODUCT INFORMATION



Reversine

Item No. 10004412

CAS Registry No.: 656820-32-5
Formal Name: N⁶-cyclohexyl-N²-[4-(4-morpholinyl)phenyl]-9H-purine-2,6-diamine
MF: C₂₁H₂₇N₇O
FW: 393.5
Purity: ≥95%
UV/Vis.: λ_{max}: 283 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

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

Reversine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, reversine should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Reversine has a solubility of approximately 0.5 mg/ml in a 1:8 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Reversine is a 2,6-disubstituted purine derivative that was originally found to induce dedifferentiation of C2C12 culture myoblast cells into stem cell progenitors when used at a concentration of 5 μM for four days.¹ Depending on cell type, reversine promotes either differentiation or dedifferentiation. For example, in NT2 neuronal and HL-60 human promyelocytic leukemia cells, it induces differentiation.² It inhibits the Aurora A, B, and C kinases with IC₅₀ values of 98-876 nM and acts as an antagonist at the adenosine A₃ receptor with a K_i value of 0.66 μM.²⁻⁴ Reversine is also used for studies of chromosome segregation. It inhibits the mitotic spindle checkpoint enzyme MPS1 with IC₅₀ values of 6 and 2.8 nM for its kinase domain and full-length version, respectively.⁴ Reversine induces autophagy in WRO human follicular thyroid cancer cells and decreases Akt/mTOR signaling.⁵

References

1. Chen, S., Zhang, Q., Wu, X., *et al.* *J. Am. Chem. Soc.* **126**(2), 410-411 (2004).
2. D'Alise, A.M., Amabile, G., Iovino, M., *et al.* *Mol. Cancer Ther.* **7**(5), 1140-1149 (2008).
3. Pereira, M., Jiang, J.K., Klutz, A.M., *et al.* *J. Med. Chem.* **48**(15), 4910-4918 (2005).
4. Santaguida, S., Tighe, A., D'Alise, A.M., *et al.* *J. Cell. Biol.* **190**(1), 73-87 (2010).
5. Lu, C.-H., Liu, Y.-W., Hua, S.-C., *et al.* *Biomed. Pharmacother.* **66**(8), 642-647 (2012).

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, 12/15/2022

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