

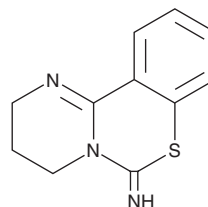
PRODUCT INFORMATION



PD 404182

Item No. 34396

CAS Registry No.: 72596-74-8
Formal Name: 3,4-dihydro-2H,6H-pyrimido[1,2-c]
[1,3]benzothiazin-6-imine
MF: C₁₁H₁₁N₃S
FW: 217.3
Purity: ≥98%
UV/Vis.: λ_{max}: 225 nm
Supplied as: A 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

PD 404182 is supplied as a solid. A stock solution may be made by dissolving the PD 404182 in the solvent of choice, which should be purged with an inert gas. PD 404182 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PD 404182 in these solvents is approximately 20, 25, and 30 mg/ml, respectively.

PD 404182 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PD 404182 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. PD 404182 has a solubility of approximately 0.09 mg/ml in a 1:10 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

PD 404182 is an antiviral agent.^{1,2} It inhibits the activity of the severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) main protease (M^{pro}), also known as 3C-like protease (3CL^{pro}; IC₅₀ = 0.081 μM), and prevents herpes simplex virus 1 (HSV-1) and HSV-2 infection in Vero cells when used at a concentration of 0.2 μM. PD 404182 also inhibits dimethylarginine dimethylaminohydrolase 1 (DDAH-1) and histone deacetylase 8 (HDAC8; IC₅₀s = 9 and 0.011 μM for the human enzymes, respectively), as well as bacterial KDO 8-P synthase (K_i = 0.026 μM), an enzyme important for LPS biosynthesis.³⁻⁵ It reduces endothelial tube formation and LPS-induced nitric oxide (NO) production in primary human dermal microvascular endothelial cells.³

References

1. Brown, A.S., Ackerley, D.F., and Calcott, M.J. High-throughput screening for inhibitors of the SARS-CoV-2 protease using a FRET-biosensor. *Molecules* **25**(20), 4666 (2020).
2. Chamoun-Emanuelli, A.M., Bobart, M., Moncla, B., et al. Evaluation of PD 404,182 as an anti-HIV and anti-herpes simplex virus microbicide. *Antimicrob. Agents Chemother.* **58**(2), 687-697 (2014).
3. Ghebremariam, Y.T., Erlanson, D.A., and Cooke, J.P. A novel and potent inhibitor of dimethylarginine dimethylaminohydrolase: a modulator of cardiovascular nitric oxide. *J. Pharmacol. Exp. Ther.* **348**(1), 69-76 (2014).
4. Kleinschek, A., Meyners, C., Digiorgio, E., et al. Potent and selective non-hydroxamate histone deacetylase 8 inhibitors. *ChemMedChem* **11**(23), 2598-2606 (2016).
5. Birck, M.R., Holler, T.P., and Woodard, R.W. Identification of a slow tight-binding inhibitor of 3-deoxy-d-manno-octulosonic acid 8-phosphate synthase. *J. Am. Chem. Soc.* **122**, 9334-9335 (2000).

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, 12/06/2022

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