

PRODUCT INFORMATION



Cilostamide

Item No. 14455

CAS Registry No.: 68550-75-4

Formal Name: N-cyclohexyl-4-[(1,2-dihydro-2-oxo-6-quinolinyloxy)-N-methylbutanamide

Synonym: OPC 3689

MF: $C_{20}H_{26}N_2O_3$

FW: 342.4

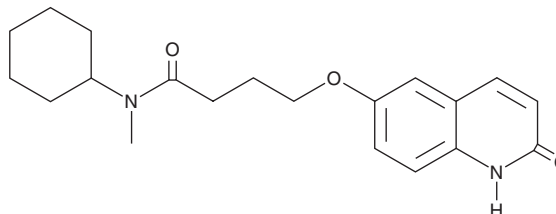
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 234, 270, 350 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

Cilostamide is supplied as a crystalline solid. A stock solution may be made by dissolving the cilostamide in the solvent of choice, which should be purged with an inert gas. Cilostamide is soluble in the organic solvent dimethyl formamide at a concentration of approximately 0.25 mg/ml.

Description

Cilostamide is a selective inhibitor of phosphodiesterase (PDE)3A and PDE3B with IC_{50} values of 27 and 50 nM, respectively.¹ It is less selective for inhibition of PDE1, PDE2, PDE4, PDE5, and PDE7 (IC_{50} s = 12 to $> 300 \mu\text{M}$).¹ Cilostamide has been shown to inhibit thrombin-induced platelet aggregation with an IC_{50} value of $1.1 \mu\text{M}$.¹

Reference

1. Sudo, T., Tachibana, K., Toga, K., *et al.* Potent effects of novel anti-platelet aggregatory cilostamide analogues on recombinant cyclic nucleotide phosphodiesterase isozyme activity. *Biochem. Pharmacol.* **59**(4), 347-356 (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, 02/18/2025

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