

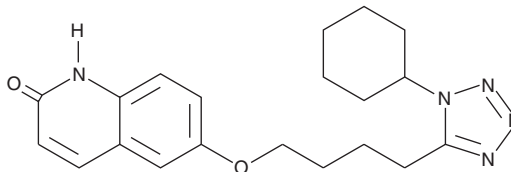
# PRODUCT INFORMATION



## 3,4-dehydro Cilostazol

Item No. 35075

**CAS Registry No.:** 73963-62-9  
**Formal Name:** 6-[4-(1-cyclohexyl-1H-tetrazol-5-yl) butoxy]-2(1H)-quinolinone  
**Synonym:** OPC 13015  
**MF:** C<sub>20</sub>H<sub>25</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 367.4  
**Purity:** ≥90%  
**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

3,4-dehydro Cilostazol is supplied as a solid. A stock solution may be made by dissolving the 3,4-dehydro cilostazol in the solvent of choice, which should be purged with an inert gas. 3,4-dehydro Cilostazol is soluble in DMSO.

### Description

3,4-dehydro Cilostazol is an active metabolite of the PDE3A inhibitor cilostazol (Item No. 15035).<sup>1,2</sup> It is formed from cilostazol by the cytochrome P450 (CYP) isoforms CYP3A4 and CYP2C19.

### References

1. Suri, A., Forbes, W.P., and Bramer, S.L. Effects of CYP3A inhibition on the metabolism of cilostazol. *Clin. Pharmacokinet.* **37**(Suppl 2), 61-68 (1999).
2. Liu, Y., Shakur, Y., Yoshitake, M., et al. Cilostazol (pletal): A dual inhibitor of cyclic nucleotide phosphodiesterase type 3 and adenosine uptake. *Cardiovasc. Drug Rev.* **19**(4), 369-386 (2001).

#### 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, 12/09/2022

#### 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](http://WWW.CAYMANCHEM.COM)