

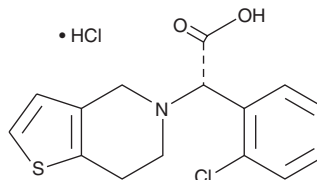
PRODUCT INFORMATION



Clopidogrel Carboxylic Acid (hydrochloride)

Item No. 17871

CAS Registry No.: 144750-42-5
Formal Name: αS-(2-chlorophenyl)-6,7-dihydro-thieno[3,2-c]pyridine-5(4H)-acetic acid, monohydrochloride
Synonym: SR26334
MF: C₁₅H₁₄ClNO₂S • HCl
FW: 344.3
Purity: ≥95%
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

Clopidogrel carboxylic acid (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the clopidogrel carboxylic acid (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Clopidogrel carboxylic acid (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of clopidogrel carboxylic acid (hydrochloride) in these solvents is approximately 50 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of clopidogrel carboxylic acid (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of clopidogrel carboxylic acid (hydrochloride) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Clopidogrel carboxylic acid is an inactive metabolite of (±)-clopidogrel (Item No. 13657) and (S)-(+)-clopidogrel (Item No. 21002).¹ It is formed from (±)-clopidogrel and (S)-(+)-clopidogrel by carboxylesterase 1 (CES1).

Reference

1. Zhu, H.-J., Wang, X., Gawronski, B.E., *et al.* Carboxylesterase 1 as a determinant of clopidogrel metabolism and activation. *J. Pharmacol. Exp. Ther.* **344**(3), 665-672 (2013).

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, 05/30/2023

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