

PRODUCT INFORMATION



LY3000328

Item No. 29729

CAS Registry No.: 1373215-15-6
Formal Name: N-[(3R,4S)-3,4-dihydro-3-[[[(methylamino)carbonyl]oxy]-6-[4-(3-oxetanyl)-1-piperazinyl]-2H-1-benzopyran-4-yl]-4-fluoro-benzamide

MF: C₂₅H₂₉FN₄O₅

FW: 484.5

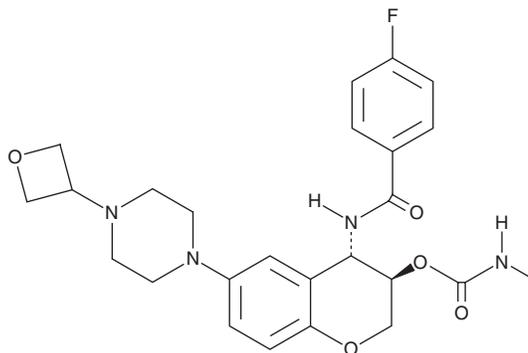
Purity: ≥98%

UV/Vis.: λ_{max}: 235 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

LY3000328 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY3000328 in the solvent of choice, which should be purged with an inert gas. LY3000328 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of LY3000328 in these solvents is approximately 30 mg/ml.

LY3000328 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LY3000328 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. LY3000328 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

LY3000328 is a potent cathepsin S inhibitor (IC₅₀ = 7.7 nM for the human enzyme).¹ It is selective for cathepsin S over cathepsin L, K, B, and V (IC₅₀s = >10,000, >10,000, 4,390, and 21,500 nM, respectively). LY3000328 (1-30 mg/kg) reduces aortic diameter in a mouse model of calcium chloride-induced abdominal aortic aneurysm.

Reference

1. Jadhav, P.K., Schiffler, M.A., Gavardinas, K., *et al.* Discovery of cathepsin S inhibitor LY3000328 for the treatment of abdominal aortic aneurysm. *ACS Med. Chem. Lett.* **5**(10), 1138-1142 (2014).

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, 11/01/2022

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