

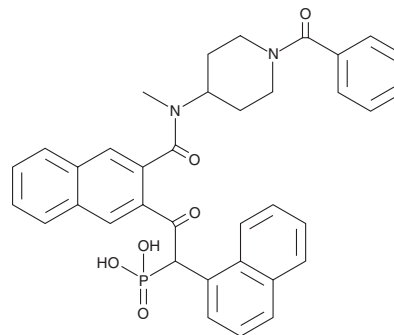
PRODUCT INFORMATION



Cathepsin G Inhibitor I

Item No. 14928

CAS Registry No.: 429676-93-7
Formal Name: P-[2-[3-[[1-benzoyl-4-piperidiny]methylamino]carbonyl]-2-naphthalenyl]-1-(1-naphthalenyl)-2-oxoethyl]-phosphonic acid
MF: C₃₆H₃₃N₂O₆P
FW: 620.6
Purity: ≥95%
UV/Vis.: λ_{max}: 224, 248 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

Cathepsin G inhibitor I is supplied as a crystalline solid. A stock solution may be made by dissolving the cathepsin G inhibitor I in the solvent of choice, which should be purged with an inert gas. Cathepsin G inhibitor I is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of cathepsin G inhibitor I in these solvents is approximately 10 and 5 mg/ml, respectively.

Cathepsin G inhibitor I is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Cathepsin G inhibitor I should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Cathepsin G inhibitor I 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

Cathepsin G inhibitor I is an inhibitor of cathepsin G (IC₅₀ = 53 nM).¹ It is selective for cathepsin G over various serine proteases, including thrombin, Factor Xa, Factor IXa, plasmin, trypsin, and leukocyte proteinase 3 (PR3; IC₅₀s = >100 μM for all), but also weakly inhibits chymotrypsin. It inhibits cathepsin G-induced activation of IL-36 in neutrophil degranulates when used at a concentration of 10 μM.² Cathepsin G inhibitor I (10 μM) reduces CD4⁺ T cell-induced secretion of IFN-γ and IL-17 in B cells incubated with tetanus toxin C-fragments (TTC) and also reduces TTC presentation in B cells.³ It decreases the size of neuromyelitis optica lesions and reduces the number of perivascular neutrophils, indicating decreased neutrophil brain entry, when used in combination with the neutrophil elastase inhibitor sivelestat (Item No. 17779) in a mouse model of IgG-induced brain injury when administered intracerebrally at a dose of 5 μg/animal.⁴

Reference

1. Greco, M.N., Hawkins, M.J., Powell, E.T., *et al.* Nonpeptide inhibitors of cathepsin G: Optimization of a novel β-ketophosphonic acid lead by structure-based drug design. *J. Am. Chem. Soc.* **124**(15), 3810-3811 (2002).
2. Henry, C.M., Sullivan, G.P., Clancy, D.M., *et al.* Neutrophil-derived proteases escalate inflammation through Aactivation of IL-36 family cytokines. *Cell Rep.* **14**(4), 708-722 (2016).
3. Reich, M., Lesner, A., Legowska, A., *et al.* Application of specific cell permeable cathepsin G inhibitors resulted in reduced antigen processing in primary dendritic cells. *Mol. Immunol.* **46**(15), 2994-2999 (2009).
4. Saadoun, S., Waters, P., MacDonald, C., *et al.* Neutrophil protease inhibition reduces neuromyelitis optica-immunoglobulin G-induced damage in mouse brain. *Ann. Neurol.* **71**(3), 323-333 (2012).

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, 11/29/2022

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