

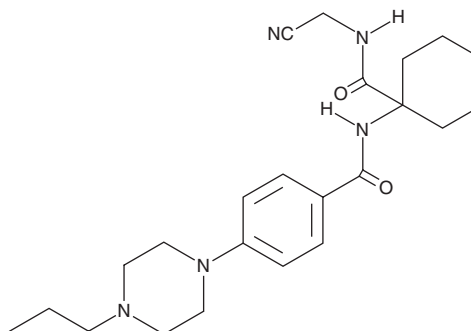
PRODUCT INFORMATION



Balicatib

Item No. 17995

CAS Registry No.: 354813-19-7
Formal Name: N-[1-[[[(cyanomethyl)amino]carbonyl]cyclohexyl]-4-(4-propyl-1-piperazinyl)-benzamide
Synonym: AAE581
MF: C₂₃H₃₃N₅O₂
FW: 411.5
Purity: ≥98%
UV/Vis.: λ_{max}: 202, 297 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

Balicatib is supplied as a crystalline solid. A stock solution may be made by dissolving the balicatib in the solvent of choice, which should be purged with an inert gas. Balicatib is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of balicatib in these solvents is approximately 0.5, 10, and 25 mg/ml, respectively.

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

Description

Balicatib is a potent inhibitor of cathepsin K (IC₅₀s = 1.4, 56, and 480 nM for human, rat, and mouse forms, respectively).^{1,2} It is at least 100-fold selective for cathepsin K over cathepsins B, L, and S.² Balicatib has a basic lipophilic nature, resulting in lysosomal trapping and increased selectivity for lysosomal cathepsin K.^{1,2}

References

- Desmarais, S., Black, W.C., Oballa, R., *et al.* Effect of cathepsin k inhibitor basicity on *in vivo* off-target activities. *Mol. Pharmacol.* **73**(1), 147-156 (2008).
- Falgueyret, J.-P., Desmarais, S., Oballa, R., *et al.* Lysosomotropism of basic cathepsin K inhibitors contributes to increased cellular potencies against off-target cathepsins and reduced functional selectivity. *J. Med. Chem.* **48**(24), 7535-7543 (2005).

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, 01/11/2022

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