

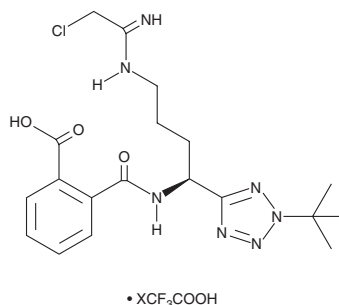
PRODUCT INFORMATION



CAY10729 (trifluoroacetate salt) (technical grade)

Item No. 26546

Formal Name: 2-[[[(1S)-4-[(2-chloro-1-iminoethyl)amino]-1-[2-(1,1-dimethylethyl)-2H-tetrazol-5-yl]butyl]amino]carbonyl]-benzoic acid, trifluoroacetate salt
MF: C₁₉H₂₆ClN₇O₃ • XCF₃COOH
FW: 435.9
Purity: ≥70%
UV/Vis.: λ_{max}: 273 nm
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

CAY10729 is a tetrazole derivative of the protein arginine deiminase (PAD) inhibitor Cl-amidine (Item No. 10599) that selectively inhibits PAD1 over PAD2, PAD3, and PAD4 ($k_{\text{inact}}/K_i = 162,200, 60,000, 17,500, \text{ and } 25,400 \text{ M}^{-1}\text{min}^{-1}$ for PAD1-4, respectively).¹

Reference

1. Subramanian, V., Knight, J.S., Parelkar, S., *et al.* Design, synthesis, and biological evaluation of tetrazole analogs of cl-amidine as protein arginine deiminase inhibitors. *J. Med. Chem.* **58**(3), 1337-1344 (2015).

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, 08/08/2019

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