

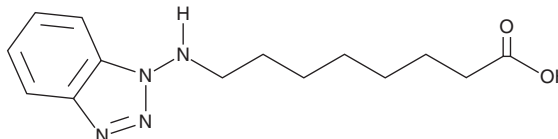
PRODUCT INFORMATION



CAY10770

Item No. 31216

CAS Registry No.: 2428734-45-4
Formal Name: 8-(1H-benzotriazol-1-ylamino)-
octanoic acid
MF: C₁₄H₂₀N₄O₂
FW: 276.3
Purity: ≥98%
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.

Laboratory Procedures

CAY10770 is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol and chloroform purged with an inert gas can be used. The solubility of CAY10770 in these solvents is approximately 1 and 10 mg/ml, respectively.

Description

CAY10770 is an inhibitor of the cytochrome P450 (CYP) isoform CYP4Z1 (IC₅₀ = 5.9 μM).¹ It is selective for CYP4Z1 over CYP4A11, CYP4F2, CYP4F3a, CYP4F3b (IC₅₀s = 187-282 μM) but does inhibit CYP4F8 and CYP4F12 (IC₅₀s = 167 and 91 μM, respectively). CAY10770 (3 μM) inhibits the production of 14(15)-EET, 19-HETE, and 14(15)-DiHET by 83, 86, and 80%, respectively, in T47D breast cancer cells expressing CYP4Z1.

Reference

1. Kowalski, J.P., McDonald, M.G., Pelletier, R.D., *et al.* Design and characterization of the first selective and potent mechanism-based inhibitor of cytochrome P450 4Z1. *J. Med. Chem.* **63(9)**, 4824-4836 (2020).

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, 08/19/2020

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