

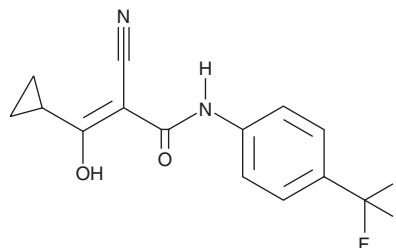
PRODUCT INFORMATION



PfDHODH-IN-1

Item No. 37728

CAS Registry No.: 183945-55-3
Formal Name: 2-cyano-3-cyclopropyl-3-hydroxy-N-[4-(trifluoromethyl)phenyl]-2-propenamide
Synonyms: CCHTFP, PfDHODH Inhibitor 1
MF: C₁₄H₁₁F₃N₂O₂
FW: 296.2
Purity: ≥98%
UV/Vis.: λ_{max}: 254, 292 nm
Supplied as: A 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

PfDHODH-IN-1 is supplied as a solid. A stock solution may be made by dissolving the PfDHODH-IN-1 in the solvent of choice, which should be purged with an inert gas. PfDHODH-IN-1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PfDHODH-IN-1 in these solvents is approximately 1, 10, and 25 mg/ml, respectively.

Description

PfDHODH-IN-1 is an inhibitor of dihydroorotate dehydrogenase (DHODH; IC₅₀s = 190, 47, and 21 nM for the human, mouse, and rat enzymes, respectively) and a derivative of the DHODH inhibitor A-771726 (Item No. 14404).^{1,2} It inhibits delayed-type hypersensitivity reactions in mice and rats when administered at doses of 30 and 10 mg/kg, respectively.²

References

1. Baldwin, J., Farajallah, A.M., Malmquist, N.A., *et al.* Malarial dihydroorotate dehydrogenase. Substrate and inhibitor specificity. *J. Biol. Chem.* **277**(44), 41827-41834 (2002).
2. Kuo, E.A., Hambleton, P.T., Kay, D.P., *et al.* Synthesis, structure-activity relationships, and pharmacokinetic properties of dihydroorotate dehydrogenase inhibitors: 2-Cyano-3-cyclopropyl-3-hydroxy-N-[3'-methyl-4'-(trifluoromethyl)phenyl]propenamide and related compounds. *J. Med. Chem.* **39**(23), 4608-4621 (1996).

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, 06/02/2023

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