

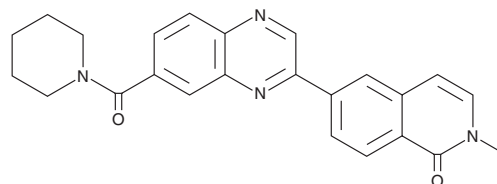
PRODUCT INFORMATION



CAY10795

Item No. 37981

CAS Registry No.: 2241676-74-2
Formal Name: 2-methyl-6-[7-(1-piperidinylcarbonyl)-2-quinoxaliny]-1(2H)-isoquinolinone
Synonyms: 15-PGDH Inhibitor,
15-hydroxy Prostaglandin Dehydrogenase Inhibitor
MF: C₂₄H₂₂N₄O₂
FW: 398.5
Purity: ≥98%
UV/Vis.: λ_{max}: 269 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

CAY10795 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10795 in the solvent of choice, which should be purged with an inert gas. CAY10795 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10795 in ethanol is approximately 25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

CAY10795 is a 15-hydroxy prostaglandin dehydrogenase (15-PGDH) inhibitor (IC₅₀ = 3 nM).¹ It increases prostaglandin E₂ (PGE₂; Item No. 14010) levels greater than 3-fold in IL-1β-stimulated A549 human lung cells when used at a concentration of 500 nM. CAY10795 reduces 15-PGDH activity in the colon of mice when administered at a dose of 20 mg/kg and limits decreases in body weight and colon length and reduces colon ulceration in a mouse model of ulcerative colitis induced by dextran sodium sulfate (DSS; Item No. 23250) when administered at a dose of 40 mg/kg. It also accelerates recovery of circulating neutrophils following whole-body radiation and bone marrow transplant in mice when administered at a dose of 10 mg/kg twice per day.

Reference

1. Hu, B., Toda, K., Wang, X., *et al.* Orally bioavailable quinoxaline inhibitors of 15-prostaglandin dehydrogenase (15-PGDH) promote tissue repair and regeneration. *J. Med. Chem.* **65(22)**, 15327-15343 (2022).

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, 02/15/2023

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