

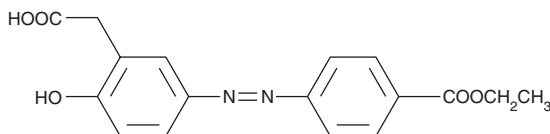
PRODUCT INFORMATION



CAY10397

Item No. 70130

CAS Registry No.: 78028-01-0
Formal Name: 5-[[4-(ethoxycarbonyl)phenyl]azo]-2-hydroxy-benzeneacetic acid
Synonym: CK47A
MF: C₁₇H₁₆N₂O₅
FW: 328.3
Purity: ≥98%
UV/Vis.: λ_{max}: 261, 370 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

CAY10397 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10397 in the solvent of choice, which should be purged with an inert gas. CAY10397 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CAY10397 in these solvents is approximately 10, 25, and 35 mg/ml, respectively. The solubility of CAY10397 in 0.1 M Na₂CO₃ is approximately 200 µg/ml.

Description

Prostaglandins are rapidly inactivated *in vivo* by the action of 15-hydroxy prostaglandin dehydrogenase (15-hydroxy PGDH). This enzyme oxidizes the 15-hydroxyl group and hydrogenates the 13,14-double bond, producing 13,14-dihydro-15 keto metabolites with greatly reduced biological activity. CAY10397 is a selective inhibitor of 15-hydroxy PGDH with an IC₅₀ of approximately 10 µM.¹ CAY10397 acts to prolong the lifetime and activity of endogenously produced prostaglandins both in cell culture and *in vivo*.

Reference

1. Berry, C.N., Hoult, J.R., Peers, S.H., *et al.* Inhibition of prostaglandin 15-hydroxydehydrogenase by sulphasalazine and a novel series of potent analogues. *Biochem. Pharmacol.* **32(19)**, 2863-2871 (1983).

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, 04/01/2024

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