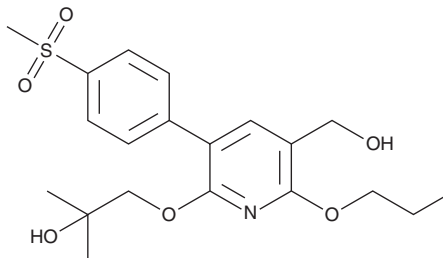


PRODUCT INFORMATION



COX-2-IN-6 Item No. 45306

CAS Registry No.: 2756347-91-6
Formal Name: 6-(2-hydroxy-2-methylpropoxy)-5-[4-(methylsulfonyl)phenyl]-2-propoxy-3-pyridinemethanol
Synonyms: COX-2 Inhibitor 6, Cyclooxygenase 2-IN-6, Cyclooxygenase 2 Inhibitor 6
MF: C₂₀H₂₇NO₆S
FW: 409.5
Purity: ≥98%
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

COX-2-IN-6 is supplied as a solid. A stock solution may be made by dissolving the COX-2-IN-6 in the solvent of choice, which should be purged with an inert gas. COX-2-IN-6 is soluble (≥10 mg/ml) in ethanol. COX-2-IN-6 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

COX-2-IN-6 is a gut-restricted COX-2 inhibitor (IC₅₀ = 0.84 μM).¹ It is selective for COX-2 over COX-1 (IC₅₀ = >50 μM). COX-2-IN-6 inhibits the production of PGE₂ (Item No. 14010) in HEK293 cells expressing COX-2 but not those expressing COX-1 (IC₅₀s = 0.6 and >50 μM, respectively). *In vivo*, COX-2-IN-6 (300 mg/kg) selectively accumulates in the colon versus the plasma at a ratio of 492:1 and reduces total ileum polyp area in an *Apc*^{Min/+} mouse model of familial adenomatous polyposis (FAP) and colorectal cancer.

Reference

1. Zhang, Z., Ghosh, A., Connolly, P. J., *et al.* Gut-restricted selective cyclooxygenase-2 (COX-2) inhibitors for chemoprevention of colorectal cancer. *J Med. Chem.* **64**(15), 11570-11596 (2021).

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, 03/17/2026

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