

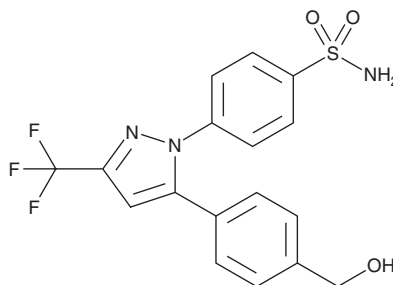
PRODUCT INFORMATION



Hydroxy Celecoxib

Item No. 33994

CAS Registry No.: 170571-00-3
Formal Name: 4-[5-[4-(hydroxymethyl)phenyl]-3-(trifluoromethyl)-1H-pyrazol-1-yl]-benzenesulfonamide
MF: C₁₇H₁₄F₃N₃O₃S
FW: 397.4
Purity: ≥95%
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

Hydroxy celecoxib is supplied as a solid. A stock solution may be made by dissolving the hydroxy celecoxib in the solvent of choice, which should be purged with an inert gas. Hydroxy celecoxib is slightly soluble in DMSO and methanol.

Description

Hydroxy celecoxib is an inactive metabolite of celecoxib (Item No. 10008672).¹ It is formed from celecoxib by the cytochrome P450 (CYP) isoforms CYP2C9, CYP3A4, and CYP2D6 in human liver microsomes.²

References

1. Penning, T.D., Talley, J.J., Bertenshaw, S.R., *et al.* Synthesis and biological evaluation of the 1,5-diarylpyrazole class of cyclooxygenase-2 inhibitors: Identification of 4-[5-(4-methylphenyl)-3-(trifluoromethyl)-1H-pyrazol-1-yl]benzenesulfonamide (SC-58635, celecoxib). *J. Med. Chem.* **40**(9), 1347-1365 (1997).
2. Siu, Y.A., Hao, M.-H., Dixit, V., *et al.* Celecoxib is a substrate of CYP2D6: Impact on celecoxib metabolism in individuals with CYP2C9*3 variants. *Drug Metab. Pharmacokinet.* **33**(5), 219-227 (2018).

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, 12/08/2022

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