

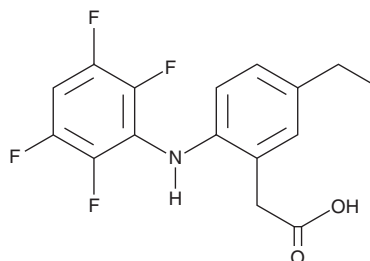
PRODUCT INFORMATION



Robenacoxib

Item No. 42218

CAS Registry No.: 220991-32-2
Formal Name: 5-ethyl-2-[(2,3,5,6-tetrafluorophenyl)amino]-benzeneacetic acid
Synonym: CGS 34975
MF: C₁₆H₁₃F₄NO₂
FW: 327.3
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

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

Description

Robenacoxib is a non-steroidal anti-inflammatory drug (NSAID) and COX-2 inhibitor ($K_i = 30$ nM for the human enzyme).¹ It is selective for COX-2 over COX-1 ($K_i = 800$ nM for the ovine enzyme). Robenacoxib inhibits IL-1 β -induced prostaglandin E₂ (PGE₂; Item No. 14010) production in human dermal fibroblasts ($IC_{50} = 31$ nM). It reduces carrageenan-induced paw edema ($ID_{50} = 0.4$ mg/kg), has antinociceptive activity in the Randall-Selitto test, and inhibits LPS-induced fever in rats. Formulations containing robenacoxib have been used in the treatment of post-operative pain and inflammation in cats.

Reference

1. King, J.N., Dawson, J., Esser, R.E., *et al.* Preclinical pharmacology of robenacoxib: A novel selective inhibitor of cyclooxygenase-2. *J. Vet. Pharmacol. Ther.* **32**(1), 1-17 (2009).

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, 10/29/2024

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