

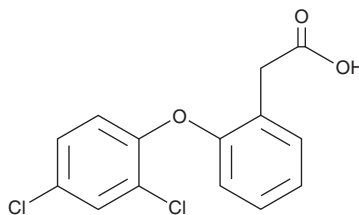
PRODUCT INFORMATION



Fenclofenac

Item No. 41709

CAS Registry No.: 34645-84-6
Formal Name: 2-(2,4-dichlorophenoxy)-
benzeneacetic acid
Synonym: RX-67408
MF: $C_{14}H_{10}Cl_2O_3$
FW: 297.1
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Fenclofenac is supplied as a solid. A stock solution may be made by dissolving the fenclofenac in the solvent of choice, which should be purged with an inert gas. Fenclofenac is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

Fenclofenac is a non-steroidal anti-inflammatory drug (NSAID) and COX inhibitor ($IC_{50}s = 7$ and $4 \mu M$ for COX-1 and COX-2, respectively).^{1,2} It reduces paw volume in a rat model of adjuvant-induced polyarthritis when administered at doses of 3, 15, and 75 mg/kg and reverses yeast-induced motor impairments, a measure of nociception, in rats ($ED_{50} = 54.3$ mg/kg).¹ Fenclofenac (20-180 mg/kg) reduces carrageenan-induced paw edema in rats. Formulations containing fenclofenac have previously been used in the treatment of rheumatoid arthritis.

References

1. Atkinson, D.C. and Leach, E.C. Anti-inflammatory and related properties of 2-(2,4-dichlorophenoxy) phenylacetic acid (fenclofenac). *Agents Actions* **6(5)**, 657-666 (1976).
2. Barnett, J., Chow, J., Ives, D., *et al.* Purification, characterization and selective inhibition of human prostaglandin G/H synthase 1 and 2 expressed in the baculovirus system. *Biochim. Biophys. Acta* **1209(1)**, 130-139 (1994).

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, 09/27/2024

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