

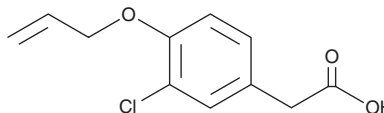
PRODUCT INFORMATION



Alclofenac

Item No. 29877

CAS Registry No.: 22131-79-9
Formal Name: 3-chloro-4-(2-propen-1-yloxy)-benzeneacetic acid
MF: C₁₁H₁₁ClO₃
FW: 226.7
Purity: ≥95%
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

Alclofenac is supplied as a crystalline solid. A stock solution may be made by dissolving the alclofenac in the solvent of choice, which should be purged with an inert gas. Alclofenac is soluble in chloroform and methanol.

Description

Alclofenac is a non-steroidal anti-inflammatory drug (NSAID) with analgesic, anti-inflammatory, and anti-pyretic properties.¹⁻³ It inhibits contractions induced by serotonin (5-HT) in isolated guinea pig ileum by 28% when used at a concentration of 200 µg/ml.⁴ Alclofenac (480 mg/kg per day) decreases implant neutrophil and macrophage infiltration in a rat model of irritant-induced inflammation induced by implantation of heat-killed *M. tuberculosis*-impregnated polyurethane cubes.⁵

References

1. Delbeke, F.T., Landuyt, J., and Debackere, M. Disposition of human drug preparations in the horse. III. Orally administered alclofenac. *J. Vet. Pharmacol. Ther.* **17(5)**, 353-358 (1994).
2. Staquet, M., Luyckx, A., and Van Cauwenberge, H. A double-blind comparison of alclofenac, pentazocine, and codeine with placebo control in pathologic pain. *J. Clin. Pharmacol. New Drugs* **11(6)**, 450-455 (1971).
3. Berry, H., Fernandes, L., Ford-Hutchinson, A.W., et al. Alclofenac and D-penicillamine. Comparative trial in rheumatoid arthritis. *Ann. Rheum. Dis.* **37(1)**, 93-97 (1978).
4. Famaey, J.P., Fontaine, J., Seaman, I., et al. A possible role of prostaglandins in guinea-pig isolated ileum contractions to serotonin. *Prostaglandins* **14(1)**, 119-124 (1977).
5. Woodland, J., Vernon-Roberts, B., Swettenham, K.V., et al. Comparison of the effects of alclofenac, flurbiprofen, and prednisolone on acute inflammatory response in the rat. *Ann. Rheum. Dis.* **36(2)**, 160-165 (1977).

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, 11/14/2022

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