

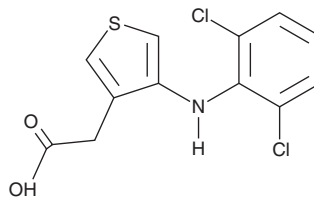
# PRODUCT INFORMATION



## Eltenac

Item No. 27673

**CAS Registry No.:** 72895-88-6  
**Formal Name:** 4-[(2,6-dichlorophenyl)amino]-3-thiopheneacetic acid  
**MF:** C<sub>12</sub>H<sub>9</sub>Cl<sub>2</sub>NO<sub>2</sub>S  
**FW:** 302.2  
**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

Eltenac is supplied as a solid. A stock solution may be made by dissolving the eltenac in the solvent of choice, which should be purged with an inert gas. Eltenac is soluble in methanol, DMSO, and dimethyl formamide.

### Description

Eltenac is a non-selective COX inhibitor (IC<sub>50</sub> = 0.03 μM for both COX-1 and COX-2 in isolated human whole blood) and a non-steroidal anti-inflammatory drug (NSAID).<sup>1</sup> *In vivo*, eltenac (1 mg/kg) reduces lameness in horses with tendinitis, pododermatitis, navicular disease, non-infectious arthritis, degenerative joint disease, phalanx fracture, and osteochondrosis.<sup>2</sup> It also reduces castration-induced edema in horses. Formulations containing eltenac have been used in the treatment of postoperative swelling in horses.

### References

1. Klein, T., Dullweber, F., Brehm, C., *et al.* Characterization of eltenac and novel COX-2 selective thiopheneacetic acid analogues in vitro and in vivo. *Biochem. Pharmacol.* **76(6)**, 717-725 (2008).
2. Prügner, W., Huber, R., and Lühmann, R. Eltenac, a new anti-inflammatory and analgesic drug for horses: clinical aspects. *J. Vet. Pharmacol. Ther.* **14(2)**, 193-199 (1991).

#### 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, 10/21/2022

#### 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