

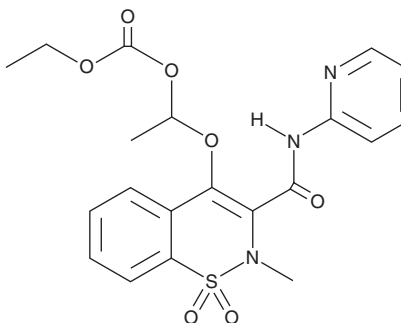
# PRODUCT INFORMATION



## Ampiroxicam

Item No. 23026

**CAS Registry No.:** 99464-64-9  
**Formal Name:** carbonic acid, ethyl 1-[[2-methyl-1,1-dioxido-3-[(2-pyridinylamino)carbonyl]-2H-1,2-benzothiazin-4-yl]oxy]ethyl ester  
**Synonym:** CP 65,703  
**MF:** C<sub>20</sub>H<sub>21</sub>N<sub>3</sub>O<sub>7</sub>S  
**FW:** 447.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 315 nm  
**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

Ampiroxicam is supplied as a crystalline solid. A stock solution may be made by dissolving the ampiroxicam in the solvent of choice, which should be purged with an inert gas. Ampiroxicam is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of ampiroxicam in these solvents is approximately 10 and 20 mg/ml, respectively.

Ampiroxicam is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ampiroxicam should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Ampiroxicam has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Ampiroxicam is a prodrug of piroxicam (Item No. 13368), a non-steroidal anti-inflammatory drug (NSAID) that non-selectively and reversibly inhibits COX-1 and COX-2 (IC<sub>50</sub>s = 1.57 and 1.69 μM, respectively).<sup>1,2</sup> Ampiroxicam is converted to piroxicam at a rate of 100, 90, 70, and 50% in humans, rats, dogs, and monkeys, respectively.<sup>1</sup>

### References

1. Carty, T.J., Marfat, A., Moore, P.F., *et al.* Ampiroxicam, an anti-inflammatory agent which is a prodrug of piroxicam. *Agents Actions* **39(3-4)**, 157-165 (1993).
2. Pairet, M., van Ryn, J., Schierok, H., *et al.* Differential inhibition of cyclooxygenases-1 and -2 by meloxicam and its 4'-isomer. *Inflamm. Res.* **47(6)**, 270-276 (1998).

#### 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