

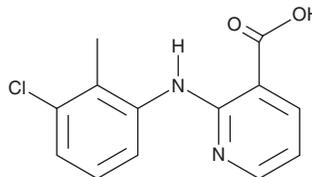
PRODUCT INFORMATION



Clonixin

Item No. 22284

CAS Registry No.: 17737-65-4
Formal Name: 2-[[3-chloro-2-methylphenyl]amino]-3-pyridinecarboxylic acid
Synonyms: CBA 93626, Clonixic Acid, NSC 335505
MF: C₁₃H₁₁ClN₂O₂
FW: 262.7
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 290, 343 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

Clonixin is supplied as a crystalline solid. A stock solution may be made by dissolving the clonixin in the solvent of choice. Clonixin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of clonixin in these solvents is approximately 3.3, 25, and 20 mg/ml, respectively.

Description

Clonixin is a non-steroidal anti-inflammatory drug (NSAID). It induces analgesia in rats in the formalin test when administered at doses of 80 and 120 mg/kg.¹ It also inhibits paw edema induced by carrageenan in rats, as well as decreases vascular permeability and acts as an antipyretic in mice.²

References

1. Bustamante, D., Miranda, H.F., Pelissier, T., *et al.* Analgesic action of clonixin, nifedipine and morphine using the formalin test. *Gen. Pharmacol.* **20(3)**, 319-322 (1989).
2. Takagi, K., Takayanagi, I., Kayaoka, S., *et al.* Pharmacological studies on clonixin. I. Antiinflammatory and analgesic actions, and effects on the gastrointestinal tract. *Nihon Yakurigaku Zasshi.* **46(5)**, 514-519 (1971).

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, 11/02/2022

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