

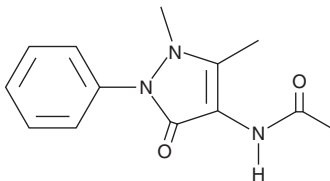
PRODUCT INFORMATION



4-Acetamido Antipyrine

Item No. 40308

CAS Registry No.: 83-15-8
Formal Name: N-(2,3-dihydro-1,5-dimethyl-3-oxo-2-phenyl-1H-pyrazol-4-yl)-acetamide
Synonyms: 4-AAAP, NSC 331807
MF: C₁₃H₁₅N₃O₂
FW: 245.3
Purity: ≥95%
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

4-Acetamido antipyrine (4-AAAP) is supplied as a solid. A stock solution may be made by dissolving the 4-AAAP in the solvent of choice, which should be purged with an inert gas. 4-AAAP is soluble in methanol and DMSO.

Description

4-AAAP is an active metabolite of the non-opioid analgesic and antipyretic prodrug metamizole (Item No. 15771).¹ It inhibits β -galactosidase activity, a marker of liver X receptor α (LXR α) activity, in a yeast two-hybrid assay (IC₂₀ = 79.9 μ M).² 4-AAAP has been found as a micropollutant in reclaimed wastewater.

References

1. Kipka, H., Tomasi, R., Hübner, M., *et al.* Simultaneous LC-ESI-MS/MS quantification of levosimendan and its metabolites for therapeutic drug monitoring of cardiac surgery patients. *Pharmaceutics* **14**(7), 1454 (2022).
2. Zhang, H., Jia, Y., Tang, Z., *et al.* Screening of chemicals with binding activities of liver X receptors from reclaimed waters. *Sci. Total Environ.* **713**:136570, (2020).

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, 02/14/2024

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