

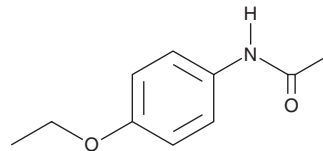
PRODUCT INFORMATION



Phenacetin

Item No. 20082

CAS Registry No.: 62-44-2
Formal Name: N-(4-ethoxyphenyl)-acetamide
Synonyms: *p*-Acetophenetidide, 4'-Ethoxyacetanilide, NSC 7651
MF: C₁₀H₁₃NO₂
FW: 179.2
Purity: ≥98%
UV/Vis.: λ_{max}: 250 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Phenacetin (Item No. 20082) is an analytical reference standard categorized as an analgesic.¹ Phenacetin has been found as an adulterant in cocaine seized by law enforcement.^{2,3} It has potential for abuse.¹ This product is intended for research and forensic applications.

References

1. Abbott, F.V. and Hellemans, K.G. Phenacetin, acetaminophen and dipyrone: Analgesic and rewarding effects. *Behav. Brain Res.* **112(1-2)**, 177-186 (2000).
2. Broséus, J., Gentile, N., and Esseiva, P. The cutting of cocaine and heroin: A critical review. *Forensic Sci. Int.* **262**, 73-83 (2016).
3. Maldaner, A.O., Botelho, É.D., Zacca, J.J., et al. Brazilian Federal District cocaine chemical profiling - mass balance approach and new adulterant routinely quantified (aminopyrine). *J. Braz. Chem. Soc.* **26(6)**, 1227-1231 (2015).

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, 06/20/2023

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