

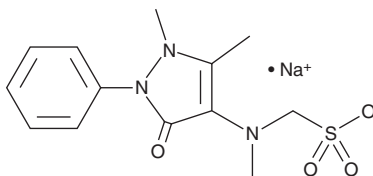
PRODUCT INFORMATION



Metamizole (sodium salt)

Item No. 32996

CAS Registry No.: 68-89-3
Formal Name: 1-[(2,3-dihydro-1,5-dimethyl-3-oxo-2-phenyl-1H-pyrazol-4-yl)methylamino]-methanesulfonic acid, monosodium salt
Synonym: Dipyrone
MF: $C_{13}H_{16}N_3O_4S \cdot Na$
FW: 333.3
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Description

Metamizole (sodium salt) (Item No. 32996) is an analytical reference standard categorized as an analgesic and anti-inflammatory agent.¹ It has been found as an adulterant in samples seized by law enforcement.^{1,2} This product is intended for research and forensic applications. This product is also available as a general research tool (Item No. 15771).

References

1. Isaacs, R.C.A., Harper, M.M., and Miller, E.C. Analytical challenges in the confirmative identification of dipyrone as an adulterant in illicit drug samples. *Forensic Sci. Int.* **270**, 185-192 (2017).
2. Fucci, N. and De Giovanni, N. Adulterants encountered in the illicit cocaine market. *Forensic Sci. Int.* **95(3)**, 247-252 (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, 02/14/2025

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