

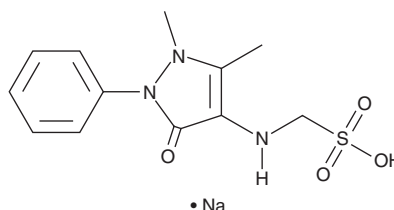
PRODUCT INFORMATION



Sulfamidopyrine (sodium salt)

Item No. 36409

CAS Registry No.: 129-89-5
Formal Name: 1-[(2,3-dihydro-1,5-dimethyl-3-oxo-2-phenyl-1H-pyrazol-4-yl)amino]-methanesulfonic acid, monosodium salt
Synonyms: 4-N-Demethylanalgin, Melaminsulfone, Sulfamidopyrine
MF: $C_{12}H_{15}N_3O_4S \cdot Na$
FW: 320.3
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 254 nm
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.

Laboratory Procedures

Sulfamidopyrine (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the sulfamidopyrine (sodium salt) in the solvent of choice, which should be purged with an inert gas. Sulfamidopyrine (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of sulfamidopyrine (sodium salt) in these solvents is approximately 0.5, 20, and 5 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of sulfamidopyrine (sodium salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of sulfamidopyrine (sodium salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Sulfamidopyrine is a derivative of the nonopioid 4-methylaminoantipyrine prodrug metamizole (Item No. 15771).

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, 08/01/2023

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