

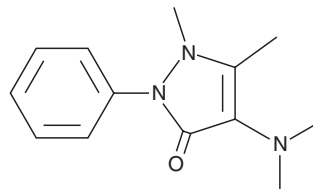
PRODUCT INFORMATION



Aminophenazone

Item No. 15770

CAS Registry No.: 58-15-1
Formal Name: 4-(dimethylamino)-1,2-dihydro-1,5-dimethyl-2-phenyl-3H-pyrazol-3-one
Synonyms: 4-Dimethylaminoantipyrine, Aminopyrine, NSC 4993
MF: C₁₃H₁₇N₃O
FW: 231.3
Purity: ≥98%
UV/Vis.: λ_{max}: 238, 271 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

Aminophenazone is supplied as a crystalline solid. A stock solution may be made by dissolving the aminophenazone in the solvent of choice. Aminophenazone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of aminophenazone in ethanol is approximately 25 mg/ml and approximately 5 mg/ml in DMSO and DMF.

Aminophenazone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, aminophenazone should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Aminophenazone has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Aminophenazone is an analgesic and anti-inflammatory compound.¹ It decreases acetic acid-induced writhing in mice (ED₅₀ = 46 mg/kg, s.c.) and reduces inflammation in a rat model of paw edema when administered at a dose of 48.2 mg/kg. Formulations containing aminophenazone have previously been used in the treatment of pain.

Reference

1. Takaya, M., Sato, M., Terashima, K., *et al.* A new nonsteroidal analgesic-antiinflammatory agent. Synthesis and activity of 4-ethoxy-2-methyl-5-morpholino-3(2H)-pyridazinone and related compounds. *J. Med. Chem.* **22**(1), 53-58 (1979).

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, 11/28/2022

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