

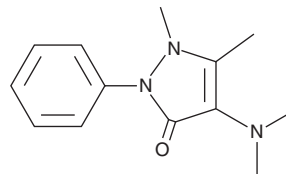
# PRODUCT INFORMATION



## Aminophenazone

Item No. 32992

**CAS Registry No.:** 58-15-1  
**Formal Name:** 4-(dimethylamino)-1,2-dihydro-1,5-dimethyl-2-phenyl-3H-pyrazol-3-one  
**Synonyms:** Aminopyrine, 4-Dimethylaminoantipyrine, NSC 4993  
**MF:** C<sub>13</sub>H<sub>17</sub>N<sub>3</sub>O  
**FW:** 231.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 241, 273 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

Aminophenazone (Item No. 32992) is an analytical reference standard categorized as an analgesic and anti-inflammatory agent.<sup>1</sup> Aminophenazone has been found as an adulterant in seized cocaine samples.<sup>2</sup> It is also a thermal degradation product of the adulterant dipyrone and has been found in seized illicit fentanyl samples.<sup>3</sup> This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 15770).

### References

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).
2. 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).
3. 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).

#### 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, 06/22/2023

#### 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