

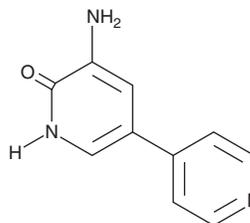
PRODUCT INFORMATION



Amrinone

Item No. 23898

CAS Registry No.: 60719-84-8
Formal Name: 5-amino-[3,4'-bipyridin]-6(1H)-one
Synonym: WIN 40,680
MF: C₁₀H₉N₃O
FW: 187.2
Purity: ≥98%
UV/Vis.: λ_{max}: 202, 246, 326 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

Amrinone is supplied as a crystalline solid. A stock solution may be made by dissolving the amrinone in the solvent of choice, which should be purged with an inert gas. Amrinone is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of amrinone in these solvents is approximately 0.5 and 0.3 mg/ml, respectively.

Description

Amrinone is an inhibitor of phosphodiesterase 3 (PDE3; IC₅₀ = 19.5 μM).¹ It increases developed tension and contractile force in isolated cat papillary muscle.¹ Amrinone (1-10 mg/kg) has positive inotropic effects, increasing the rate and force of heart contraction in anesthetized and unanesthetized dogs.

Reference

1. Alousi, A.A., Farah, A.E., Leshar, G.Y., *et al.* Cardiotoxic activity of amrinone--Win 40680 [5-amino-3,4'-bipyridine-6(1H)-one]. *Circ. Res.* **45(5)**, 666-677 (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, 11/01/2022

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