

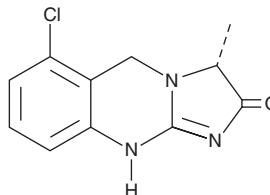
# PRODUCT INFORMATION



## Quazinone

Item No. 29151

**CAS Registry No.:** 70018-51-8  
**Formal Name:** (3R)-6-chloro-1,5-dihydro-3-methyl-imidazo[2,1-b]quinazolin-2(3H)-one  
**Synonym:** Ro 13-6438  
**MF:** C<sub>11</sub>H<sub>10</sub>ClN<sub>3</sub>O  
**FW:** 235.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 254 nm  
**Supplied as:** A 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

Quazinone is supplied as a solid. A stock solution may be made by dissolving the quazinone in the solvent of choice, which should be purged with an inert gas. Quazinone is soluble in organic solvents such as ethanol and DMSO. The solubility of quazinone in these solvents is approximately 1 and 5 mg/ml, respectively.

### Description

Quazinone is a selective inhibitor of phosphodiesterase 3 (PDE3) with positive inotropic and vasodilating properties.<sup>1,2</sup> It induces relaxation of precontracted isolated human cavernous smooth muscle (IC<sub>50</sub> = 4.2 μM).<sup>1</sup> Quazinone (10-300 μg/kg) increases myocardial contractile force in anesthetized open-chest dogs in a dose-dependent manner, as well as decreases systolic and diastolic blood pressure.<sup>2</sup> It also inhibits DNA synthesis induced by the PDGF isoform PDGF-BB in bovine coronary artery smooth muscle cells in a concentration-dependent manner.<sup>3</sup>

### References

1. Taher, A., Meyer, M., Stief, C.G., *et al.* Cyclic nucleotide phosphodiesterase in human cavernous smooth muscle. *World J. Urol.* **15(1)**, 32-35 (1997).
2. Eigenmann, R., Gerold, M., and Holck, M. Cardiovascular profile of Ro 13-6438, a novel positive inotropic agent with vasodilating properties. *J. Cardiovasc. Pharmacol.* **6(3)**, 511-519 (1984).
3. Osinski, M.T. and Schrör, K. Inhibition of platelet-derived growth factor-induced mitogenesis by phosphodiesterase 3 inhibitors: Role of protein kinase A in vascular smooth muscle cell mitogenesis. *Biochem. Pharmacol.* **60(3)**, 381-387 (2000).

#### 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, 09/30/2022

#### 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