

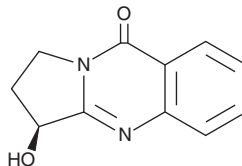
PRODUCT INFORMATION



Vasicinone

Item No. 11761

CAS Registry No.: 486-64-6
Formal Name: (3S)-2,3-dihydro-3-hydroxy-pyrrolo[2,1-b]quinazolin-9(1H)-one
Synonym: (-)-Vasicinone
MF: C₁₁H₁₀N₂O₂
FW: 202.2
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 267, 301, 313 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

Vasicinone is supplied as a crystalline solid. A stock solution may be made by dissolving the vasicinone in the solvent of choice. Vasicinone is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of vasicinone in these solvents is approximately 20 mg/ml.

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

Description

Vasicinone is a quinazoline alkaloid that can be isolated from *A. vasica* leaves and demonstrates bronchodilatory activity.¹ It also is reported to possess antioxidant activity in nitric oxide and ABTS radical scavenging assays.²

References

1. Amin, A.H. and Mehta, D.R. A bronchodilator alkaloid (vasicinone) from *Adhatoda vasica* Nees. *Nature* **184**(Suppl 17), 1317 (1959).
2. Nilani, P., Kasthuribai, N., Duraisamy, B., *et al.* *In vitro* antioxidant activity of selected antiasthmatic herbal constituents. *Anc. Sci. Life* **28**(4), 3-6 (2009).

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, 11/14/2022

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