

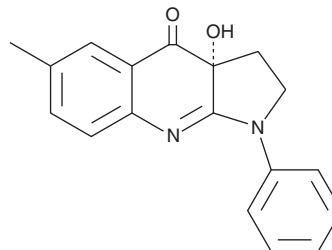
PRODUCT INFORMATION



(+)-Blebbistatin

Item No. 13165

CAS Registry No.: 1177356-70-5
Formal Name: 1,2,3,3a-tetrahydro-3aR-hydroxy-6-methyl-1-phenyl-4H-pyrrolo[2,3-b]quinolin-4-one
Synonym: (R)-Blebbistatin
MF: C₁₈H₁₆N₂O₂
FW: 292.3
Purity: ≥98%
UV/Vis.: λ_{max}: 236, 270, 298, 418 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

(+)-Blebbistatin is supplied as a crystalline solid. A stock solution may be made by dissolving the (+)-blebbistatin in an organic solvent purged with an inert gas, which should be purged with an inert gas. (+)-Blebbistatin is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of (+)-blebbistatin in these solvents is approximately 12.5 and 20 mg/ml, respectively.

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

Description

(+)-Blebbistatin is the inactive enantiomer of (-)-blebbistatin (Item No. 13013), which is a selective inhibitor of non-muscle myosin II ATPases.^{1,2} Prolonged exposure to blue light (450-490 nm) results in degradation of blebbistatin to an inactive product via cytotoxic intermediates, which may be problematic for its use in fluorescent live cell imaging applications.^{3,4} (R)-nitro-Blebbistatin (Item No. 9001935) is a more stable form of (+)-blebbistatin.⁵ The addition of a nitro group stabilizes the molecule to circumvent its degradation by prolonged blue light exposure. (R)-nitro-Blebbistatin has the same stereochemistry as the inactive (+)-blebbistatin enantiomer.

References

1. Wang, X., Grammatikakis, N., Siganou, A., *et al.* Regulation of molecular chaperone gene transcription involves the serine phosphorylation, 14-3-3e binding, and cytoplasmic sequestration of heat shock factor 1. *Mol. Cell. Biol.* **23**(17), 6013-6026 (2003).
2. Kovács, M., Tóth, J., Hetényi, C., *et al.* Mechanism of blebbistatin inhibition of myosin II. *J. Biol. Chem.* **279**(34), 35557-35563 (2004).
3. Kolega, J. Phototoxicity and photoinactivation of blebbistatin in UV and visible light. *Biochem. Biophys. Res. Commun.* **320**(3), 1020-1025 (2004).
4. Sakamoto, T., Limouze, J., Combs, C.A., *et al.* Blebbistatin, a myosin II inhibitor, is photoinactivated by blue light. *Biochemistry* **44**(2), 584-588 (2005).
5. Lucas-Lopez, C., Patterson, S., Blum, T., *et al.* Absolute stereochemical assignment and fluorescence tuning of the small molecule tool, (-)-blebbistatin. *Eur. J. Org. Chem.* 1736-1740 (2005).

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, 03/11/2024

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