

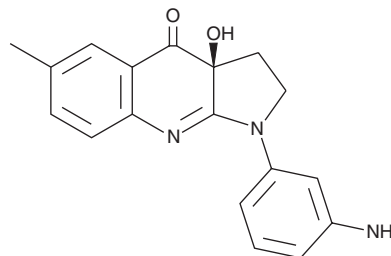
# PRODUCT INFORMATION



## (S)-3'-amino Blebbistatin

Item No. 24170

**CAS Registry No.:** 2097141-18-7  
**Formal Name:** (3aS)-1-(3-aminophenyl)-1,2,3,3a-tetrahydro-3a-hydroxy-6-methyl-4H-pyrrolo[2,3-b]quinolin-4-one  
**Synonyms:** (-)-3'-amino Blebbistatin, *m*-amino Blebbistatin, *meta*-amino Blebbistatin  
**MF:** C<sub>18</sub>H<sub>17</sub>N<sub>3</sub>O<sub>2</sub>  
**FW:** 307.3  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 244, 264, 298, 419 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

(S)-3'-amino Blebbistatin is supplied as a crystalline solid. A stock solution may be made by dissolving the (S)-3'-amino blebbistatin in the solvent of choice. (S)-3'-amino Blebbistatin is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of (S)-3'-amino blebbistatin in these solvents is approximately 12.5 and 20 mg/ml, respectively.

(S)-3'-amino Blebbistatin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (S)-3'-amino blebbistatin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. (S)-3'-amino 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

(S)-3'-amino Blebbistatin is a more stable and less phototoxic form of (-)-blebbistatin (Item No. 13013), which is a selective cell-permeable inhibitor of non-muscle myosin II ATPases.<sup>1,2</sup> (-)-Blebbistatin rapidly and reversibly inhibits Mg-ATPase activity and *in vitro* motility of non-muscle myosin IIA and IIB for several species (IC<sub>50</sub>s = 0.5-5 μM), while poorly inhibiting smooth muscle myosin (IC<sub>50</sub> = 80 μM).<sup>3</sup> Through these effects, it blocks apoptosis-related bleb formation, directed cell migration, and cytokinesis in vertebrate cells. However, 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.<sup>4,5</sup> The addition of a 3'-amino group decreases the inherent fluorescence while retaining the activity of (-)-blebbistatin.<sup>6</sup> (S)-3'-amino Blebbistatin has the same stereochemistry as the active (-)-blebbistatin enantiomer.

### References

1. Straight, A.F., Cheung, A., Limouze, J., *et al.* *Science* **299**(5613), 1743-1747 (2003).
2. Kovács, M., Tóth, J., Hetényi, C., *et al.* *J. Biol. Chem.* **279**(34), 35557-35563 (2004).
3. Limouze, J., Straight, A.F., Mitchison, T., *et al.* *J. Muscle Res. Cell Motil.* **25**(4-5), 337-341 (2004).
4. Kolega, J. *Biochem. Biophys. Res. Commun.* **320**(3), 1020-1025 (2004).
5. Sakamoto, T., Limouze, J., Combs, C.A., *et al.* *Biochemistry* **44**(2), 584-588 (2005).
6. Verhasselt, S., Roman, B.I., De Wever, O., *et al.* *Org. Biomol. Chem.* **15**(9), 2104-2118 (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, 07/31/2018

#### 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