

# PRODUCT INFORMATION



## Blasticidin S Item No. 35252

CAS Registry No.: 2079-00-7

Formal Name: 4-[[[(3S)-3-amino-5-[(aminoiminomethyl)methylamino]-1-oxopentyl]amino]-1-(4-amino-2-oxo-1(2H)-pyrimidinyl)-1,2,3,4-tetrahydro-β-D-erythro-hex-2-enopyranuronic acid

MF: C<sub>17</sub>H<sub>26</sub>N<sub>8</sub>O<sub>5</sub>

FW: 422.4

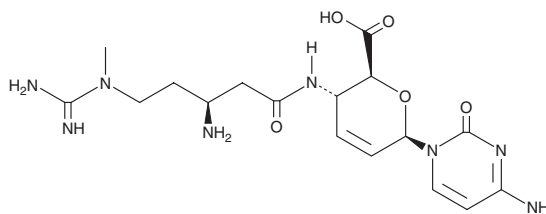
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Streptomyces* sp.



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

### Laboratory Procedures

Blasticidin S is supplied as a solid. A stock solution may be made by dissolving the blasticidin S in the solvent of choice, which should be purged with an inert gas. Blasticidin S is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Blasticidin S is a bacterial metabolite originally isolated from *S. griseochromogenes* that has antibacterial and fungicidal activities.<sup>1</sup> It is active against *B. subtilis*, *S. lutea*, *E. coli*, *P. fluorescens*, and *M. tuberculosis*, as well as the plant pathogenic fungi *P. oryzae*, *S. mali*, *G. kaki*, and *G. lacticolor* in a broth dilution assay when used at concentrations ranging from 5 to 50 µg/ml.

### Reference

1. Takeuchi, S., Hirayama, K., Ueda, K., *et al.* Blasticidin S, a new antibiotic. *J. Antibiot. (Tokyo)* **11(1)**, 1-5 (1958).

#### 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, 01/03/2023

#### 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