

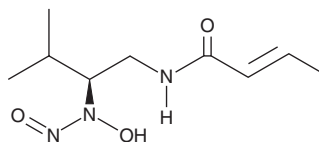
# PRODUCT INFORMATION



## Dopastin

Item No. 32896

**CAS Registry No.:** 37134-80-8  
**Formal Name:** (2E)-N-[(2S)-2-(hydroxynitrosoamino)-3-methylbutyl]-2-butenamide  
**Synonym:** NSC 252927  
**MF:** C<sub>9</sub>H<sub>17</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 215.3  
**Purity:** ≥90%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Bacterium/*Pseudomonas*



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

### Laboratory Procedures

Dopastin is supplied as a solid. A stock solution may be made by dissolving the dopastin in the solvent of choice, which should be purged with an inert gas. Dopastin is soluble in methanol, butanol, acetone, and chloroform.

### Description

Dopastin is a bacterial metabolite that has been found in *Pseudomonas* and an inhibitor of dopamine β-hydroxylase ( $K_i = 4.7 \mu\text{M}$ ).<sup>1,2</sup> *In vivo*, dopastin (20 mg/kg) reduces mean systolic blood pressure in spontaneously hypertensive rats. Dopastin also inhibits the germination of barley seeds ( $\text{IC}_{50} = 47 \mu\text{M}$ ).<sup>2</sup>

### References

1. Iinuma, H., Takeuchi, T., Kondo, S., *et al.* Dopastin, an inhibitor of dopamine β-hydroxylase. *J. Antibiot. (Tokyo)* **25(8)**, 497-500 (1972).
2. Iinuma, H., Matsuzaki, M., Nagatsu, T., *et al.* Biochemical and biological studies on dopastin, an inhibitor of dopamine β-hydroxylase. *Agric. Biol. Chem.* **38(11)**, 2107-2111 (1974).

#### 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, 11/10/2022

#### 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](http://WWW.CAYMANCHEM.COM)