

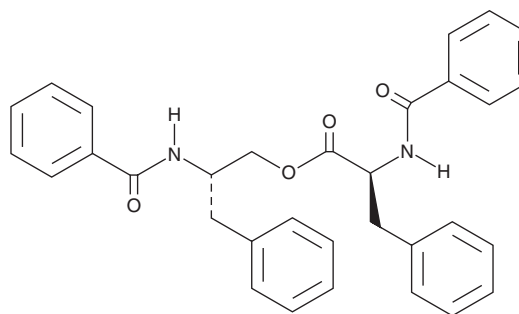
# PRODUCT INFORMATION



## Asperphenamate

Item No. 25086

**CAS Registry No.:** 63631-36-7  
**Formal Name:** N-benzoyl-L-phenylalanine,2S-(benzoylamino)-3-phenylpropyl ester  
**Synonyms:** Anabellamide, Auranamide, NSC 306231  
**MF:** C<sub>32</sub>H<sub>30</sub>N<sub>2</sub>O<sub>4</sub>  
**FW:** 506.6  
**Purity:** ≥95%  
**Supplied as:** A 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

Asperphenamate is supplied as a solid. A stock solution may be made by dissolving the asperphenamate in the solvent of choice, which should be purged with an inert gas. Asperphenamate is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

### Description

Asperphenamate is a fungal secondary metabolite originally isolated from *A. flavipes*.<sup>1</sup> It inhibits proliferation of T47D and MDA-MB-231 breast and HL-60 leukemia cancer cell lines (IC<sub>50</sub>s = 92.3, 96.5, and 97.9 μM, respectively) and inhibits yeast α-glucosidase activity (IC<sub>50</sub> = 8.3 μM).<sup>2,3</sup>

### References

1. Clark, A.M. and Hufford, C.D. Synthesis of asperphenamate, a novel fungal metabolite. *Phytochem.* **17(3)**, 552-553 (1978).
2. Yuan, L., Wang, J.H., and Sun, T.M. Total synthesis and anticancer activity studies of the stereoisomers of asperphenamate and patriscabratine. *Chinese Chem. Lett.* **21(2)**, 155-158 (2010).
3. Del Valle, P., Martinez, A.L., Figueroa, M., et al. Alkaloids from the fungus *Penicillium spathulatum* as α-glucosidase inhibitors. *Planta Med.* **82(14)**, 1286-1294 (2016).

#### 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, 12/06/2022

#### 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