

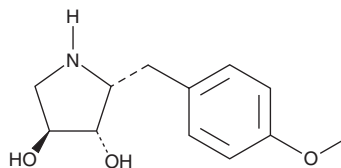
# PRODUCT INFORMATION



## Deacetylanisomycin

Item No. 28067

**CAS Registry No.:** 27958-06-1  
**Formal Name:** (2R,3S,4S)-2-[(4-methoxyphenyl)methyl]-3,4-pyrrolidinediol  
**Synonyms:** (-)-Deacetylanisomycin, SA 3097D1  
**MF:** C<sub>12</sub>H<sub>17</sub>NO<sub>3</sub>  
**FW:** 223.3  
**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

Deacetylanisomycin is supplied as a solid. A stock solution may be made by dissolving the deacetylanisomycin in the solvent of choice. Deacetylanisomycin is soluble in the organic solvent methanol.

### Description

Deacetylanisomycin is a derivative of the protein and DNA synthesis inhibitor anisomycin (Item No. 11308) with antimalarial and anticancer activity.<sup>1,2</sup> It inhibits the growth of the *P. falciparum* strains K1 and T9-96 *in vitro* (IC<sub>50</sub>s = 40.29 and 33.57 μM, respectively).<sup>1</sup> Deacetylanisomycin is cytotoxic to LU99 lung carcinoma and MCF-7 breast cancer cells (IC<sub>50</sub>s = 23 and 34 μM, respectively). It has been used as a negative control for assessment of protein synthesis inhibitor activity in molluscan neurons.<sup>3</sup>

### References

1. Ekong, R.M., Kirby, G.C., Patel, G., *et al.* Comparison of the *in vitro* activities of quassinoids with activity against *Plasmodium falciparum*, anisomycin and some other inhibitors of eukaryotic protein synthesis. *Biochem. Pharmacol.* **40(2)**, 297-301 (1990).
2. Hosoya, Y., Kameyama, T., Naganawa, H., *et al.* Anisomycin and new congeners active against human tumor cell lines. *J. Antibiot. (Tokyo)* **46(8)**, 1300-1302 (1993).
3. Alkon, D.L., Bank, B., Naito, S., *et al.* Inhibition of protein synthesis prolongs Ca<sup>2+</sup>-mediated reduction of K<sup>+</sup> currents in molluscan neurons. *Proc. Nat. Acad. Sci. USA* **84(19)**, 6948-6952 (1987).

#### 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, 12/09/2022

#### 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