

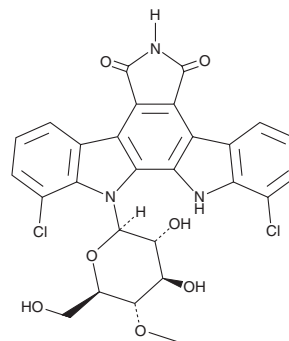
# PRODUCT INFORMATION



## Rebeccamycin

Item No. 17914

**CAS Registry No.:** 93908-02-2  
**Formal Name:** 1,11-dichloro-12,13-dihydro-12-(4-O-methyl-β-D-glucopyranosyl)-5H-indolo[2,3-a]pyrrolo[3,4-c]carbazole-5,7(6H)-dione  
**Synonyms:** BRN 4732638, NSC 359079  
**MF:** C<sub>27</sub>H<sub>21</sub>Cl<sub>2</sub>N<sub>3</sub>O<sub>7</sub>  
**FW:** 570.4  
**Purity:** ≥99%  
**Supplied as:** A lyophilisate  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

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

### Description

Rebeccamycin is an indolocarbazole antibiotic produced by *S. aerocolonigenes* ATCC39243 that weakly inhibits DNA topoisomerase I.<sup>1,2</sup> It is cytotoxic to human lung adenocarcinoma, colon carcinoma, and nasopharyngeal carcinoma cell lines (IC<sub>50</sub>s range from 0.4-98 μM), producing single-strand breaks in the DNA of these tumor cells.<sup>3</sup>

### References

1. Long, B.H., Rose, W.C., Vyas, D.M., *et al.* Discovery of antitumor indolocarbazoles: Rebeccamycin, NSC 655649, and fluoroindolocarbazoles. *Curr. Med. Chem. Anticancer Agents* **2(2)**, 255-266 (2002).
2. Sánchez, C., Butovich, I.A., Braña, A.F., *et al.* The biosynthetic gene cluster for the antitumor rebeccamycin: Characterization and generation of indolocarbazole derivatives. *Chem. Biol.* **9(4)**, 519-531 (2002).
3. Bush, J.A., Long, B.H., Catino, J.J., *et al.* Production and biological activity of rebeccamycin, a novel antitumor agent. *J. Antibiot. (Tokyo)* **40(5)**, 668-678 (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, 11/02/2022

#### 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