

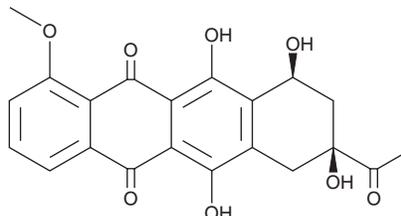
# PRODUCT INFORMATION



## Daunorubicinone

Item No. 32912

**CAS Registry No.:** 21794-55-8  
**Formal Name:** (8S,10S)-8-acetyl-7,8,9,10-tetrahydro-6,8,10,11-tetrahydroxy-1-methoxy-5,12-naphthacenedione  
**Synonyms:** (+)-Daunomycinone, NSC 109531  
**MF:** C<sub>21</sub>H<sub>18</sub>O<sub>8</sub>  
**FW:** 398.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 233, 250 nm  
**Supplied as:** A crystalline 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

Daunorubicinone is supplied as a crystalline solid. A stock solution may be made by dissolving the daunorubicinone in the solvent of choice, which should be purged with an inert gas. Daunorubicinone is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of daunorubicinone in DMSO and DMF is approximately 30 mg/ml. Daunorubicinone is slightly soluble in ethanol.

Daunorubicinone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, daunorubicinone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Daunorubicinone has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Daunorubicinone is an aglycone form of the antitumor antibiotic daunorubicin (Item No. 14159).<sup>1</sup> It has been used as a starting material in the synthesis of anticancer agents.<sup>2,3</sup>

### References

1. Arcamone, F., Franceschi, G., Orezzi, P., et al. Daunomycin. I. The structure of daunomycinone. *J. Am. Chem. Soc.* **86(23)**, 5334-5335 (1964).
2. Aligiannis, N., Pouli, N., Marakos, P., et al. Synthesis and cytotoxic activity of a new potent daunomycinone derivative. *Bioorg. Med. Chem.* **12(24)**, 3505-3507 (2002).
3. Baer, H.H. and Siemsen, L. Synthesis and biological activity of (S)-2'-fluorodaunorubicin. *Can. J. Chem.* **66(1)**, 187-190 (1988).

#### 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, 12/13/2022

#### 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