

# PRODUCT INFORMATION



**NSC 15520**

Item No. 26020

**CAS Registry No.:** 730960-98-2

**Formal Name:** 4,4a,4b,5,6,7,8,8a,9,10-decahydro-4b,8-dimethyl-2-(1-methylethyl)-3H-3,10a-ethanophenanthrene-8,11,12-tricarboxylic acid

**MF:** C<sub>24</sub>H<sub>34</sub>O<sub>6</sub>

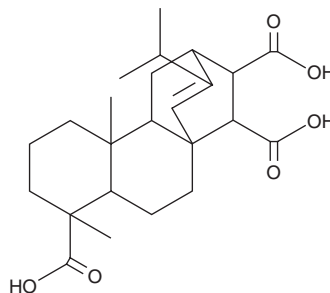
**FW:** 418.5

**Purity:** ≥98%

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥2 years



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

## Laboratory Procedures

NSC 15520 is supplied as a crystalline solid. A stock solution may be made by dissolving the NSC 15520 in the solvent of choice, which should be purged with an inert gas. NSC 15520 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of NSC 15520 in these solvents is approximately 30 mg/ml.

## Description

NSC 15520 is an inhibitor of replication protein A 70 (RPA70) dsDNA binding.<sup>1</sup> It selectively inhibits dsDNA over ssDNA binding to RPA70 when used at concentrations ranging from 25 to 400 μM. NSC 15520 also inhibits the protein-protein interaction between the cell cycle checkpoint control proteins p53 or Rad9 and RPA70.

## Reference

1. Glanzer, J.G., Liu, S., and Oakley, G.G. Small molecule inhibitor of the RPA70 N-terminal protein interaction domain discovered using in silico and in vitro methods. *Bioorg. Med. Chem.* **19(8)**, 2589-2595 (2011).

### 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, 02/21/2020

## 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)