

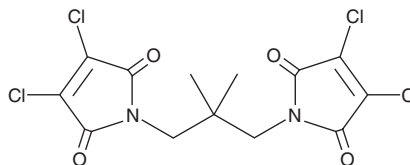
# PRODUCT INFORMATION



**NSC 617145**

Item No. 21973

**CAS Registry No.:** 203115-63-3  
**Formal Name:** 1,1'-(2,2-dimethyl-1,3-propanediyl)bis[3,4-dichloro-1H-pyrrole-2,5-dione]  
**MF:** C<sub>13</sub>H<sub>10</sub>Cl<sub>4</sub>N<sub>2</sub>O<sub>4</sub>  
**FW:** 400.0  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 240 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

NSC 617145 is supplied as a crystalline solid. A stock solution may be made by dissolving the NSC 617145 in the solvent of choice, which should be purged with an inert gas. NSC 617145 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of NSC 617145 in these solvents is approximately 1 and 2 mg/ml, respectively.

NSC 617145 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NSC 617145 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. NSC 617145 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

NSC 617145 is an inhibitor of WRN helicase (IC<sub>50</sub> = 230 nM) that inhibits the ATPase, but not exonuclease, activity of WRN helicase in a concentration-dependent manner.<sup>1</sup> It is selective for WRN over BLM, FANCF, ChlR1, RecQ, and UvrD helicases, only exhibiting 7% inhibition of RECQ1 at a concentration of 5 μM. NSC 617145 inhibits growth of HeLa cells via formation of DNA double strand breaks (DSBs) and induction of apoptosis in a WRN helicase-dependent manner. NSC 617145 also acts synergistically with mitomycin C (Item No. 11435) to inhibit growth as well as induce DSBs and chromosomal abnormalities in Fanconi-anemia deficient (FA-D2<sup>-/-</sup>) cells.

## Reference

1. Aggarwal, M., Banerjee, T., Sommers, J.A., *et al.* Werner syndrome helicase has a critical role in DNA damage responses in the absence of a functional fanconi anemia pathway. *Cancer Res.* **73(17)**, 5497-5507 (2013).

### 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, 11/28/2022

## 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