

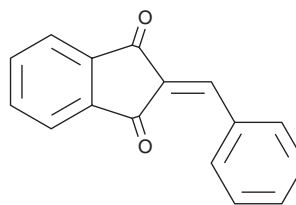
# PRODUCT INFORMATION



## PRT4165

Item No. 19093

**CAS Registry No.:** 31083-55-3  
**Formal Name:** 2-(3-pyridinylmethylene)-1H-indene-1,3(2H)-dione  
**Synonym:** NSC 600157  
**MF:** C<sub>15</sub>H<sub>9</sub>NO<sub>2</sub>  
**FW:** 235.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 245, 264, 330 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

PRT4165 is supplied as a crystalline solid. A stock solution may be made by dissolving the PRT4165 in the solvent of choice, which should be purged with an inert gas. PRT4165 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PRT4165 in these solvents is approximately 1, 2, and 10 mg/ml, respectively.

PRT4165 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PRT4165 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. PRT4165 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

BMI1/RING1A and BMI1/RNF2 are E3 ubiquitin ligase complexes that mediate the monoubiquitination of histone 2A (H2A). This is an essential function of polycomb repressive complex 1 (PRC1).<sup>1</sup> PRT4165 is an inhibitor of the ubiquitin ligase activity of PRC1.<sup>2,3</sup> It blocks BMI1/RING1A self-ubiquitination (IC<sub>50</sub> = 3.9 μM), as well as polyubiquitination of topoisomerase 2α in cells.<sup>2</sup> PRT4165 inhibits ubiquitination of H2A by either RNF2 or RING1. Through these actions, PRT4165 prevents the accumulation of detectable ubiquitin at DNA double-strand breaks, the retention of response proteins around breaks, and the repair of breaks.<sup>3</sup>

### References

1. Cao, J. and Yan, Q. Histone ubiquitination and deubiquitination in transcription, DNA damage response, and cancer. *Frontiers in Oncology* **2(26)**, 1-9 (2012).
2. Alchanati, I., Teicher, C., Cohen, G., et al. The E<sub>3</sub> ubiquitin-ligase Bmi1/Ring1A controls the proteasomal degradation of Top2α cleavage complex - a potentially new drug target. *PLoS One* **4(12)**, (2009).
3. Ismail, I.H., McDonald, D., Strickfaden, H., et al. A small molecule inhibitor of polycomb repressive complex 1 inhibits ubiquitin signaling at DNA double-strand breaks. *J. Biol. Chem.* **288(37)**, 26944-26954 (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