

PRT-4165 (2-pyridine-3-yl-methylene-indan-1,3-dione)

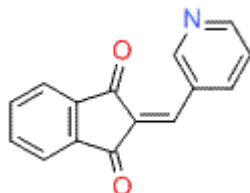
Cat. # SI9885

Background: A potent inhibitor of histone H2A ubiquitylation. Inhibits RNF2 and Ring1A, both of which are found in polycomb repressive complex 1 (PRC1). Does not inhibit RNF8 or RNF168. Inhibits accumulation of ubiquitin around DNA double-strand breaks (DSBs) as well as the repair of DSBs (1).

Application: Studying chromatin remodeling

Product Information

CAS No.	31083-55-3
Purity:	≥ 98% (TLC); NMR (conforms)
Molecular Weight:	235.2
Physical State:	Dry yellow powder
Quantity:	5mg
Solubility:	25mg/mL in DMSO
Storage:	Store powder dessicated at ambient temperature, solutions at -20°C

**Formula:** C₁₅H₉NO₂**References**

1. Ismail, H., McDonald, D. Strickfaden, H., Xu, Z., and Hendzei, M.J. 2013. A small molecule inhibitor of polycomb repressive complex 1 inhibits ubiquitin signaling at DNA double-strand breaks. *J. Biol. Chem.* **288**:26944-26954.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2015 LifeSensors, Inc. All Rights Reserved