

TCID

4,5,6,7-tetrachloro-1H-indene-1,3(2H)-dione

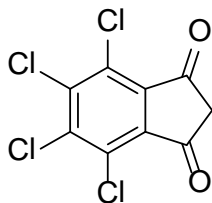
Cat. # SI9679

Background: Inhibits ubiquitin C-terminal hydrolase L3 (UCH-L3). IC₅₀=0.6 μM, 125-fold selective over UCH-L1, IC₅₀=75 μM. Cell permeable.

Application: Cell-based and in vitro studies of UCH-L3 function

Product Information:

CAS No.	30675-13-9
Purity:	>97% (TLC); NMR (Conforms)
Molecular Weight:	283.9
Physical State:	White Powder
Quantity:	10 mg, 50 mg
Solubility:	DMSO (25mg/ml)
Storage:	Store desiccated as supplied at ambient temperature for up to 1 year. Store solutions at -20°C for up to 3 months.



Formula: C₉H₂Cl₄O₂

References

1. Love K.R. *et al.* Mechanisms, biology and inhibitors of deubiquitinating enzymes. *Nat. Chem. Biol.* 2007 **3**:697
2. Mtango, N.R. *et al.* Essential role of ubiquitin C-terminal hydrolases UCHL1 and UCHL3 in mammalian oocyte maturation. *J. Cell Physiol.* 2012 **227**:2022

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2011 LifeSensors, Inc. All Rights Reserved