

## LDN-57444

(5-Chloro-1-[(2,5-dichlorophenyl)methyl]-1H-indole-2,3-dione 3-(O-acetyloxime)

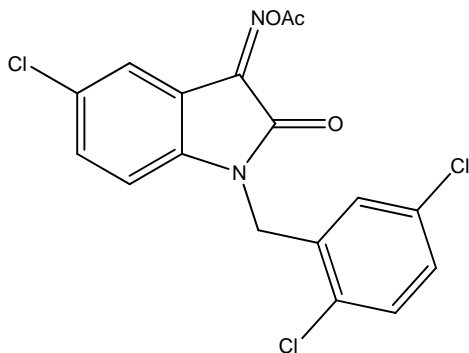
Cat. # SI9639

**Background:** LDN-57444 is a cell permeable ubiquitin C-terminal hydrolase (UCH-L1) inhibitor ( $K_i=0.4 \mu\text{M}$ ) with approximately 28-fold greater selectivity over UCH-L3. Decreases proteasome activity and increases levels of ubiquitinated proteins. Induces apoptosis and causes dramatic alterations in synaptic protein distribution and spine morphology in vivo. LDN-57444 is a useful tool to study UCH-L1 in different pathological conditions such as cancer and neurodegeneration.

**Application:** Inhibition of ubiquitin C-terminal hydrolase.

### Product Information

<b>CAS No.</b>	668467-91-2
<b>Purity:</b>	> 98% (TLC); NMR (Conforms)
<b>Molecular Weight:</b>	397.6
<b>Physical State:</b>	Yellow powder
<b>Quantity:</b>	10 mg, 50 mg
<b>Solubility:</b>	DMSO (20mg/ml)
<b>Storage:</b>	Store desiccated as supplied at 4°C for up to 2 years. Store solutions at -20°C for up to one week. Protect from light.



**Formula:**  $\text{C}_{17}\text{H}_{11}\text{Cl}_3\text{N}_2\text{O}_3$

### References

1. Tan et al Endoplasmic reticulum stress contributes to the cell death induced by UCH-L1 inhibitor. *Mol.Cell.Biochem.* **318**: 109 (2008).
2. Susor, A., et al., Proteomic analysis of porcine oocytes during in vitro maturation reveals essential role for the ubiquitin C-terminal hydrolase-L1. *Reprod.* **134**: 559 - 568, (2007)

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