

NSC-632839

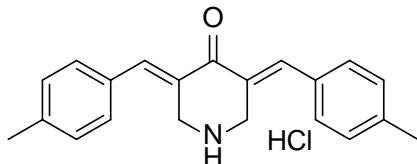
3,5-Bis[(4-methylphenyl)methylene]-4-piperidone hydrochloride

Cat. # SI9689

Background: A pan-specific ubiquitin & ubiquitin-like isopeptidase inhibitor. Inhibits USP2, USP7, and SENP2 with EC₅₀ values of 45 ± 4 μM, 37 ± 1 μM, and 9.8 ± 1.8 μM, respectively. Does not affect proteasome proteolytic activity. Induces apoptosis via a Bcl-2-dependent but apoptosome-independent mitochondrial pathway. Cell permeable.

Product Information:

CAS No.	157654-67-6
Purity:	>98% (HPLC); NMR (Conforms)
Molecular Weight:	339.9
Physical State:	Yellow Powder
Quantity:	10 mg, 50 mg
Solubility:	May be dissolved in: DMSO (8 mg/ml)
Storage:	Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month.



Formula: C₂₁H₂₁NO · HCl

References

1. Aleo E. *et al.* Identification of New Compounds That Trigger Apoptosome-Independent Caspase Activation and Apoptosis. *Cancer Res.* 2006 **66**:9235
2. Nicholson B. *et al.* Characterization of ubiquitin and ubiquitin-like-protein isopeptidase activities. *Protein Sci.* 2008 **17**:1035

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