

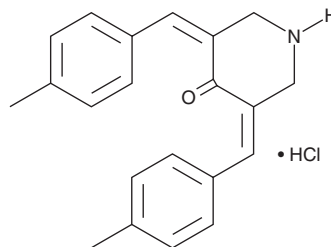
PRODUCT INFORMATION



NSC 632839

Item No. 16893

CAS Registry No.: 157654-67-6
Formal Name: 3,5-bis[4-(4-methylphenyl)methylene]-4-piperidinone, monohydrochloride
Synonyms: F6, Ubiquitin Isopeptidase Inhibitor II
MF: C₂₁H₂₁NO • HCl
FW: 339.9
Purity: ≥98%
UV/Vis.: λ_{max}: 240, 345 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

NSC 632839 is supplied as a crystalline solid. A stock solution may be made by dissolving the NSC 632839 in the solvent of choice, which should be purged with an inert gas. NSC 632839 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of NSC 632839 in ethanol is approximately 0.25 mg/ml and approximately 10 mg/ml.

Description

Protein ubiquitination is a dynamic process that can be reversed by deubiquitinating enzymes (DUBs) that remove ubiquitin from proteins, sparing them from degradation by the proteasome. The DUBs have been divided into various groups, the largest being the ubiquitin-specific proteases (USPs). NSC 632839 inhibits USP2 and USP7 as well as the ubiquitin-like SUMO peptidase SENP2 with EC₅₀ values of 45, 37, and 9.8 μM, respectively.¹ It is reported to inhibit apoptosis through a Bcl-2-dependent and apoptosome-independent pathway of caspase activation (IC₅₀ = 15.7 μM in cells sensitized by E1A, an adenoviral oncogene).²

References

1. Nicholson, B., Leach, C.A., Goldenberg, S.J., *et al.* Characterization of ubiquitin and ubiquitin-like-protein isopeptidase activities. *Protein Sci.* **17(6)**, 1035-1043 (2008).
2. Aleo, E., Henderson, C.J., Fontanini, A., *et al.* Identification of new compounds that trigger apoptosome-independent caspase activation and apoptosis. *Cancer Res.* **66(18)**, 9235-9244 (2006).

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, 10/05/2022

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