

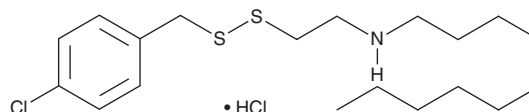
# PRODUCT INFORMATION



**NSC 624206**

Item No. 20569

**CAS Registry No.:** 13116-77-3  
**Formal Name:** N-[2-[[[4-chlorophenyl)methyl]dithio]ethyl]-1-decanamine, monohydrochloride  
**MF:** C<sub>19</sub>H<sub>32</sub>ClNS<sub>2</sub> • HCl  
**FW:** 410.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 225 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 624206 is supplied as a crystalline solid. A stock solution may be made by dissolving the NSC 624206 in the solvent of choice. NSC 624206 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of NSC 624206 in ethanol and DMSO is approximately 1 mg/ml and approximately 0.16 mg/ml in DMF.

## Description

NSC 624206 is an inhibitor of ubiquitin-activating enzyme (E1) that specifically blocks ubiquitin-thioester formation (IC<sub>50</sub> = 13 μM) but has no effect on ubiquitin adenylation.<sup>1</sup> Through this action, NSC 624206 prevents the ubiquitination and degradation of the tumor suppressor protein p27 *in vitro*.<sup>1</sup>

## Reference

1. Ungermannova, D., Parker, P.J., Nasveschuk, C.G., *et al.* Identification and mechanistic studies of a novel ubiquitin E1 inhibitor. *J. Biomol. Screen.* **17**(4), 421-434 (2012).

**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, 11/03/2022

## 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