

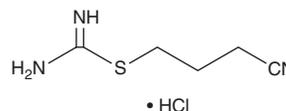
# PRODUCT INFORMATION



## Kevetrin (hydrochloride)

Item No. 28288

**CAS Registry No.:** 66592-89-0  
**Formal Name:** carbamimidothioic acid, 3-cyanopropyl ester, monohydrochloride  
**Synonyms:** 4-Isothioureidobutyronitrile, NSC 525990  
**MF:** C<sub>5</sub>H<sub>9</sub>N<sub>3</sub>S • HCl  
**FW:** 179.7  
**Purity:** ≥95%  
**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

Kevetrin (hydrochloride) is supplied as a crystalline solid. Aqueous solutions of kevetrin (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of kevetrin (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Kevetrin is an activator of p53.<sup>1,2</sup> It increases levels of activated p53 and expression of p21 in A549 cells and decreases expression of the transcription factor E2F1 in a variety of cell lines.<sup>1,2</sup> Kevetrin decreases cell viability in a panel of cancer cell lines with a mean IC<sub>50</sub> value of 0.49 μM.<sup>3</sup> It reduces tumor growth in MDA-MB-231, HT-29, PC3, HCT15, A549, NCI-H1975, CRL-1619, LNCaP, MIA PaCa, and SCC-5 mouse xenograft models when administered at a dose of 200 mg/kg.

### References

1. Kumar, A., Hiran, T., Holden, S.A., *et al.* Abstract 4470: Kevetrin<sup>TM</sup>, a novel small molecule, activates p53, enhances expression of p21, induces cell cycle arrest and apoptosis in a human cancer cell line. *Cancer Res.* **71(8 Supp)**, 4470 (2011).
2. Kumar, A., Holden, S.A., Chafai-Fadela, K., *et al.* Abstract 2874: Kevetrin targets both MDM2-p53 and Rb-E2F pathways in tumor suppression. *Cancer Res.* **72(8 Supp)**, 2874 (2012).
3. Menon, K. Nitrile derivatives and their pharmaceutical use and compositions. *Cellceutix Corporation.* **US 2012/0189537A I** (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, 12/05/2022

#### 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](http://WWW.CAYMANCHEM.COM)