

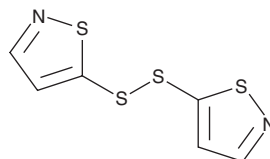
# PRODUCT INFORMATION



**NU 9056**

Item No. 18118

**CAS Registry No.:** 1450644-28-6  
**Formal Name:** 5,5'-dithiobis-isothiazole  
**MF:** C<sub>6</sub>H<sub>4</sub>N<sub>2</sub>S<sub>4</sub>  
**FW:** 232.4  
**Purity:** ≥85%  
**Supplied as:** A liquid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

## Laboratory Procedures

NU 9056 is supplied as a liquid. A stock solution may be made by dissolving the NU 9056 in the solvent of choice, which should be purged with an inert gas. NU 9056 is soluble (≥10 mg/ml) in ethanol and DMSO.

## Description

NU 9056 was originally identified as an inhibitor of TIP60, also known as lysine acetyltransferase 5 (KAT5; IC<sub>50</sub> = 2 μM).<sup>1</sup> However, it has been shown that NU 9056 is thiol-reactive in a cell-free assay and inhibits multiple targets in biochemical and cell-based assays, indicating it is a non-selective interference compound.<sup>2</sup> NU 9056 inhibits the growth of LNCaP, CWR22Ru1, and PC3 prostate cancer cells (GI<sub>50</sub>s = 24, 7.5, and 27 μM, respectively) and induces apoptosis in LNCaP cells.<sup>1</sup> It inhibits activation of the NOD-like receptor protein 3 (NLRP3) inflammasome in LPS-primed nigericin-induced primary mouse peritoneal macrophages.<sup>3</sup> NU 9056 (10 mg/kg) decreases lung injury in a mouse model of LPS-induced endotoxemia. It increases freezing time seven days after fear conditioning, indicating enhanced remote memory, in mice when administered at a dose of 2 μg/g.<sup>4</sup>

## References

1. Coffey, K., Blackburn, T.J., Cook, S., *et al.* Characterisation of a Tip60 specific inhibitor, NU9056, in prostate cancer. *PLoS One* **7(10)**, e45539 (2012).
2. Dahlin, J.L., Nelson, K.M., Strasser, J.M., *et al.* Assay interference and off-target liabilities of reported histone acetyltransferase inhibitors. *Nat. Commun.* **8(1)**, 1527 (2017).
3. Zhang, Y., Luo, L., Xu, X., *et al.* Acetylation is required for full activation of the NLRP3 inflammasome. *Nat. Commun.* **14(1)**, 8396 (2023).
4. Narkaj, K., Stefanelli, G., Wahdan, M., *et al.* Blocking H2A.Z incorporation via Tip60 inhibition promotes systems consolidation of fear memory in mice. *eNeuro* **5(5)**, (2018).

### 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, 01/05/2026

## 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