

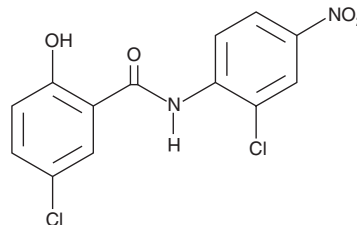
# PRODUCT INFORMATION



## Niclosamide

Item No. 10649

**CAS Registry No.:** 50-65-7  
**Formal Name:** 5-chloro-N-(2-chloro-4-nitrophenyl)-2-hydroxybenzamide  
**Synonym:** NSC 178296  
**MF:** C<sub>13</sub>H<sub>8</sub>Cl<sub>2</sub>N<sub>2</sub>O<sub>4</sub>  
**FW:** 327.1  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 333 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

Niclosamide is supplied as a crystalline solid. A stock solution may be made by dissolving the niclosamide in the solvent of choice, which should be purged with an inert gas. Niclosamide is soluble in the organic solvent ethanol at a concentration of approximately 0.5 mg/ml.

### Description

Niclosamide is an anthelmintic that disrupts mitochondrial metabolism in parasitic worms and animal models.<sup>1-3</sup> It inhibits STAT3 transcriptional activity (IC<sub>50</sub> = 0.25 μM in a reporter assay) and stimulates autophagy by reversibly inhibiting mammalian target of rapamycin complex 1 (mTORC1) signaling in MCF-7 cells expressing EGFP-LC3.<sup>2,4</sup> It completely inhibits phosphorylation of ribosomal S6 kinases (S6K) when used at a concentration of 3 μM.<sup>2</sup> Niclosamide (1 μM) induces mitochondrial uncoupling, decreasing cellular ATP concentrations and increasing the ratio of ADP to ATP in HepG2 cells, and activates AMP-activated protein kinase (AMPK) in NIH3T3 cells.<sup>3</sup> It also increases lipid oxidation by 5-fold in HepG2 cells. Niclosamide (5 and 10 μM) inhibits replication of severe acute respiratory coronavirus 2 (SARS-CoV-2) and Middle East respiratory syndrome CoV (MERS-CoV).<sup>5</sup> It prevents high-fat diet-induced hepatic steatosis and insulin resistance and improves glycemic control in *db/db* mice when administered at a dose of 125 mg/kg per day.<sup>3</sup> Niclosamide also inhibits proliferation and colony formation of DU145 prostate cancer cells, which have constitutively active STAT3 (IC<sub>50</sub>s = 0.7 and 0.1 μM, respectively).<sup>4</sup>

### References

1. Pampori, N.A., Singh, G., and Srivastava, V.M. Energy metabolism in *Cotugnia digonopora* and the effect of anthelmintics. *Mol. Biochem. Parasitol.* **11**, 205-213 (1984).
2. Balgi, A.D., Fonseca, B.D., Donohue, E., et al. Screen for chemical modulators of autophagy reveals novel therapeutic inhibitors of mTORC1 signaling. *PLoS One* **4**(9), e7124 (2009).
3. Tao, H., Zhang, Y., Zeng, X., et al. Niclosamide ethanalamine-induced mild mitochondrial uncoupling improves diabetic symptoms in mice. *Nat. Med.* **20**(11), 1263-1269 (2014).
4. Ren, X., Duan, L., He, Q., et al. Identification of niclosamide as a new small-molecule inhibitor of the STAT3 signaling pathway. *ACS Med. Chem. Lett.* **1**(9), 454-459 (2010).
5. Mostafa, A., Kandeil, A., Elshaier, Y.A.M.M., et al. FDA-approved drugs with potent in vitro antiviral activity against severe acute respiratory syndrome coronavirus 2. *Pharmaceuticals (Basel)* **13**(12), 443 (2020).

#### 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, 11/04/2022

#### 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