

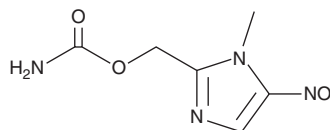
# PRODUCT INFORMATION



## Ronidazole

Item No. 33071

**CAS Registry No.:** 7681-76-7  
**Formal Name:** 1-methyl-5-nitro-1H-imidazole-2-methanol,  
2-carbamate  
**Synonym:** Ridzol  
**MF:** C<sub>6</sub>H<sub>8</sub>N<sub>4</sub>O<sub>4</sub>  
**FW:** 200.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 301 nm  
**Supplied as:** A 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

Ronidazole is supplied as a solid. A stock solution may be made by dissolving the ronidazole in the solvent of choice, which should be purged with an inert gas. Ronidazole is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ronidazole in these solvents is approximately 30 mg/ml.

Ronidazole is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ronidazole should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Ronidazole has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Ronidazole is a carbamate ester with antiprotozoal and antibiotic activities.<sup>1</sup> It is active against a panel of *T. foetus* isolates with minimum lethal concentrations (MLCs) ranging from 0.625 to 1.25 µg/ml. It is also active against *C. difficile* bacteria (MIC<sub>90</sub> = 0.125 µg/ml).<sup>2</sup> Ronidazole (50 mg/kg) inhibits infection relapse in a cat model of experimental *T. foetus* infection.<sup>3</sup> It also inhibits *T. muris* infection in mice when administered at a concentration of 400 mg/L in drinking water.<sup>4</sup> Ronidazole (10 mg/kg) increases survival in a mouse model of *C. difficile* infection.<sup>2</sup> Formulations containing ronidazole have been used in the treatment of *T. foetus* infection in cats.

### References

1. Kather, E.J., Marks, S.L., and Kass, P.H. Determination of the in vitro susceptibility of feline *Tritrichomonas foetus* to 5 antimicrobial agents. *J. Vet. Intern. Med.* **21(5)**, 966-970 (2007).
2. AbdelKhalek, A. and Seleem, M.N. Repurposing the veterinary antiprotozoal drug ronidazole for the treatment of *Clostridioides difficile* infection. *Int. J. Antimicrob. Agents* **56(6)**, 106188 (2020).
3. Gookin, J.L., Copple, C.N., Papich, M.G., et al. Efficacy of ronidazole for treatment of feline *Tritrichomonas foetus* infection. *J. Vet. Intern. Med.* **20(3)**, 536-543 (2006).
4. Steiner, J.M., Schwamberger, S., Pantchev, N., et al. Use of ronidazole and limited culling to eliminate *Tritrichomonas muris* from laboratory mice. *J. Am. Assoc. Lab. Anim. Sci.* **55(4)**, 480-483 (2016).

#### 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, 10/04/2022

#### 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