

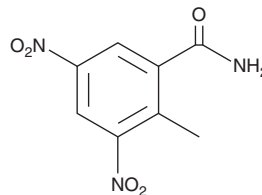
# PRODUCT INFORMATION



## Dinitolmide

Item No. 31138

CAS Registry No.: 148-01-6  
Formal Name: 2-methyl-3,5-dinitro-benzamide  
Synonyms: Dinitrotoluamide, Zoalene  
MF:  $C_8H_7N_3O_5$   
FW: 225.2  
Purity: 98%  
UV/Vis.:  $\lambda_{max}$ : 238 nm  
Supplied as: A solid  
Storage: -20°C  
Stability:  $\geq 4$  years



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

### Laboratory Procedures

Dinitolmide is supplied as a solid. A stock solution may be made by dissolving the dinitolmide in the solvent of choice, which should be purged with an inert gas. Dinitolmide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of dinitolmide in these solvents is approximately 0.5 mg/ml in ethanol and 15 mg/ml in DMSO and DMF.

Dinitolmide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, dinitolmide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Dinitolmide has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Dinitolmide is an anticoccidial agent.<sup>1</sup> Dietary administration of dinitolmide (125-500 ppm) reduces mortality and decreases the number of oocysts and oocyst sporulation in chick models of *E. maxima*, *E. brunetti*, or *E. acervulina* infection. Formulations containing dinitolmide have been used in the prevention and treatment of coccidiosis in poultry.

### Reference

1. Joyner, L.P. and Norton, C.C. The anticoccidial effects of amprolium, dinitolmide and monensin against *Eimeria maxima*, *E. brunetti* and *E. acervulina* with particular reference to oocyst sporulation. *Parasitology* 75(2), 155-164 (1977).

#### 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, 11/09/2022

#### 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