

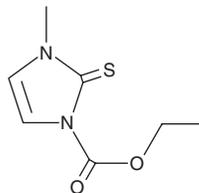
PRODUCT INFORMATION



Carbimazole

Item No. 28280

CAS Registry No.: 22232-54-8
Formal Name: 2,3-dihydro-3-methyl-2-thioxo-1H-imidazole-1-carboxylic acid, ethyl ester
MF: C₇H₁₀N₂O₂S
FW: 186.2
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 301, 352, 371 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

Carbimazole is supplied as a crystalline solid. A stock solution may be made by dissolving the carbimazole in the solvent of choice, which should be purged with an inert gas. Carbimazole is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of carbimazole in these solvents is approximately 1 mg/ml.

Carbimazole is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, carbimazole should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Carbimazole has a solubility of approximately 0.1 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Carbimazole is a prodrug form of the thyroid hormone synthesis inhibitor methimazole (Item No. 23718).¹ Carbimazole (15 mg/animal) reduces total thyroxine concentration, thyroid enlargement, polyphagia, anorexia, and alopecia in a cat model of hyperthyroidism.

Reference

1. Frénais, R., Rosenberg, D., Burgaud, S., *et al.* Clinical efficacy and safety of a once-daily formulation of carbimazole in cats with hyperthyroidism. *J. Small Anim. Pract.* **50(10)**, 510-515 (2009).

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