

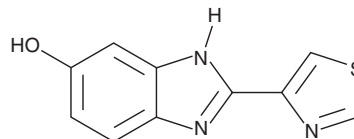
# PRODUCT INFORMATION



## 5-hydroxy Thiabendazole

Item No. 25021

CAS Registry No.: 948-71-0  
Formal Name: 2-(4-thiazolyl)-1H-benzimidazol-6-ol  
Synonym: 5-OH TBZ  
MF:  $C_{10}H_7N_3OS$   
FW: 217.2  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{\max}$ : 250, 318 nm  
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

5-hydroxy Thiabendazole (5-OH TBZ) is supplied as a crystalline solid. A stock solution may be made by dissolving the 5-OH TBZ in the solvent of choice. 5-OH TBZ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 5-OH TBZ is approximately 0.5 mg/ml in ethanol and approximately 20 mg/ml in DMSO and DMF.

5-OH TBZ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 5-OH TBZ should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 5-OH TBZ has a solubility of approximately 0.05 mg/ml in a 1:20 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

5-OH TBZ is a major metabolite of the anthelmintic thiabendazole (Item No. 23391).<sup>1</sup> Unlike thiabendazole, 5-OH TBZ has no effect on the growth of third-stage *A. caninum* larvae.

### Reference

1. Stone, O.J., Mullins, J.F., and Willis, C.J. Comparison of thiabendazole and 5-hydroxy-thiabendazole (for anthelmintic effect). *J. Invest. Dermatol.* **45**(2), 132-133 (1965).

#### 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, 12/05/2022

#### 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