

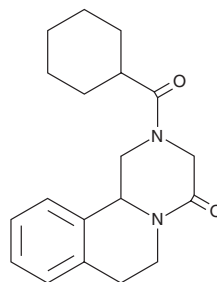
# PRODUCT INFORMATION



## Praziquantel

Item No. 20508

**CAS Registry No.:** 55268-74-1  
**Formal Name:** 2-(cyclohexylcarbonyl)-1,2,3,6,7,11b-hexahydro-4H-pyrazino[2,1-a]isoquinolin-4-one  
**Synonyms:** Embay 8440, (±)-Praziquantel  
**MF:** C<sub>19</sub>H<sub>24</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 312.4  
**Purity:** ≥98%  
**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

Praziquantel is supplied as a crystalline solid. A stock solution may be made by dissolving the praziquantel in the solvent of choice, which should be purged with an inert gas. Praziquantel is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of praziquantel in these solvents is approximately 10, 20, and 30 mg/ml, respectively.

Praziquantel is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, praziquantel should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Praziquantel 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

Praziquantel is an anthelmintic agent.<sup>1,2</sup> It is lethal to *S. mansoni* schistosomula and adults (LC<sub>50</sub>s = 0.68 and 0.03 mg/ml, respectively).<sup>2</sup> Praziquantel (172 and 592 mg/kg) reduces worm burden in a mouse model of *S. mansoni* infection. Formulations containing praziquantel have been used in the treatment of schistosomiasis.

### References

1. Cook, G.C. Anthelmintic agents: Some recent developments and their clinical application. *Postgrad. Med. J.* **67(783)**, 16-22 (1991).
2. Olliaro, P., Delgado-Romero, P. and Keiser, J. The little we know about the pharmacokinetics and pharmacodynamics of praziquantel (racemate and R-enantiomer). *J. Antimicrob. Chemother.* **69(4)**, 863-870 (2014).

#### 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, 11/30/2022

#### 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