

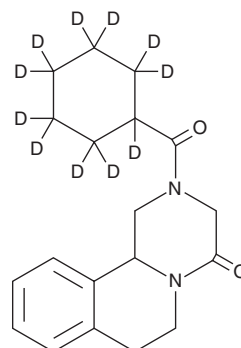
PRODUCT INFORMATION



Praziquantel-d₁₁

Item No. 30237

CAS Registry No.: 1246343-36-1
Formal Name: 2-(cyclohexyl-1,2,2,3,3,4,4,5,5,6,6-d₁₁-carbonyl)-1,2,3,6,7,11b-hexahydro-4H-pyrazino[2,1-a]isoquinolin-4-one
Synonym: (±)-Praziquantel-d₁₁
MF: C₁₉H₁₃D₁₁N₂O₂
FW: 323.5
Chemical Purity: ≥95% (Praziquantel)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₁₁); ≤1% d₀
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

Praziquantel-d₁₁ is intended for use as an internal standard for the quantification of praziquantel (Item No. 20508) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Praziquantel-d₁₁ is supplied as a solid. A stock solution may be made by dissolving the praziquantel-d₁₁ in the solvent of choice, which should be purged with an inert gas. Praziquantel-d₁₁ is soluble in chloroform and methanol.

Description

Praziquantel is an anthelmintic agent.^{1,2} It is lethal to *S. mansoni* schistosomula and adults (LC₅₀s = 0.68 and 0.03 mg/ml, respectively).² 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, 05/07/2024

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