

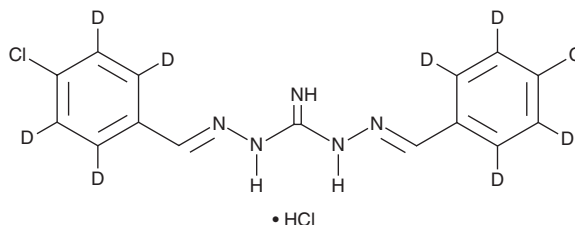
PRODUCT INFORMATION



Robenidine-d₈ (hydrochloride)

Item No. 43810

CAS Registry No.: 1173097-77-2
Formal Name: 2,2'-bis[(4-chlorophenyl-2,3,5,6-d₄)-methylene]-carbonimidic dihydrazide, monohydrochloride
MF: C₁₅H₅Cl₂D₈N₅ • HCl
FW: 378.7
Chemical Purity: ≥98% (Robenidine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₈); ≤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

Robenidine-d₈ is intended for use as an internal standard for the quantification of robenidine (Item no. 43792) 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).

Robenidine-d₈ is supplied as a solid. A stock solution may be made by dissolving the robenidine-d₈ in the solvent of choice, which should be purged with an inert gas. Robenidine-d₈ is soluble (≥10 mg/ml) in DMSO.

Description

Robenidine is an anticoccidial agent.¹ It is also active against methicillin-resistant *S. aureus* (MRSA) and vancomycin-resistant enterococci (VRE; MICs = 6.1 and 7.4 μM, respectively).² Dietary administration of robenidine reduces weight gain, decreases intestinal lesion and oocyst numbers, and increases survival in *E. acervuline*-, *E. brunetti*-, *E. mivati*-, *E. necatrix*-, or *E. tenella*-infected chickens. Formulations containing robenidine have been used in animal feed for the prevention of coccidiosis or control of necrotic enteritis in chickens.

References

1. Kennett, R.L., Kantor, S., and Gallo, A. Efficacy studies with robenidine, a new type of anticoccidial, in the diet. *Poult. Sci.* **53**(3), 978-986 (1974).
2. Abraham, R.J., Stevens, A.J., Young, K.A., et al. Robenidine analogues as gram-positive antibacterial agents. *J. Med. Chem.* **59**(5), 2126-2138 (2016).

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, 07/29/2025

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