

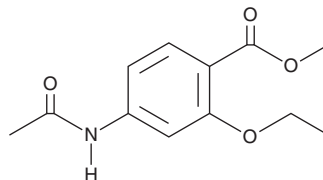
PRODUCT INFORMATION



Ethopabate

Item No. 36079

CAS Registry No.: 59-06-3
Formal Name: 4-(acetylamino)-2-ethoxy-benzoic acid, methyl ester
Synonym: Ethyl Pabate
MF: C₁₂H₁₅NO₄
FW: 237.3
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 272 nm
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

Ethopabate is supplied as a solid. A stock solution may be made by dissolving the ethopabate in the solvent of choice, which should be purged with an inert gas. Ethopabate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ethopabate in ethanol is approximately 2 mg/ml and approximately 5 mg/ml in DMSO and DMF.

Description

Ethopabate is a coccidiostat.^{1,2} *In vivo*, ethopabate reduces lesion score in broiler chickens infected with *E. tenella*, *E. acervulina*, and *E. maxima* field isolates when administered in combination with the antiprotozoal agent amprolium (Item No. 25506). Ethopabate has been found as a contaminant in environmental waters.³ Formulations containing ethopabate have been used to prevent coccidiosis in broiler chickens.

References

1. Mathis, G.F. and McDougald, L.R. Drug responsiveness of field isolates of chicken *Coccidia*. *Poult. Sci.* **61**(1), 38-45 (1982).
2. Oe, O. and Arakawa, A. Effect of feed additive antibiotics on chickens infected with *Eimeria tenella*. *Poult. Sci.* **54**(4) 1008-1018 (1975).
3. Mooney, D., Coxon, C., Richards, K.G., *et al.* A new sensitive method for the simultaneous chromatographic separation and tandem mass spectrometry detection of anticoccidials, including highly polar compounds, in environmental waters. *J. Chromatogr. A* **1618**, 460857 (2020).

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, 12/19/2022

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