

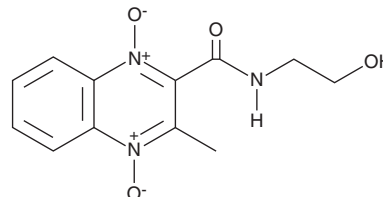
PRODUCT INFORMATION



Olaquinox

Item No. 27725

CAS Registry No.: 23696-28-8
Formal Name: N-(2-hydroxyethyl)-3-methyl-2-quinoxalinecarboxamide 1,4-dioxide
Synonym: NSC 634933
MF: C₁₂H₁₃N₃O₄
FW: 263.3
Purity: ≥98%
UV/Vis.: λ_{max}: 231, 268, 385 nm
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

Olaquinox is supplied as a crystalline solid. A stock solution may be made by dissolving the olaquinox in the solvent of choice.

Description

Olaquinox is an antibiotic and a swine growth regulator.¹ *In vivo*, olaquinox (100 mg/kg) decreases *E. coli*-induced diarrhea and the number of intraepithelial lymphocytes in the ileum and increases body weight in geld piglets. Olaquinox induces bleeding, intramuscular edema, vacuolar degeneration, and deformation of renal tubules in *C. carpio*.² It decreases mean body weight and increases the number of dead fetuses in pregnant rats when administered at a dose of 110 mg/kg and increases the number of external deformities, including bent tail, craniorachischisis, exencephaly, and anophthalmia in rat pups born to exposed pregnant mothers.³ Formulations containing olaquinox have been used as growth promoting agents in bovine livestock.

References

1. Ding, M.X., Yuan, Z.H., Wang, Y.L., *et al.* Olaquinox and cyadox stimulate growth and decrease intestinal mucosal immunity of piglets orally inoculated with *Escherichia coli*. *J. Anim. Physiol. Anim. Nutr. (Berl.)* **90**(5-6), 238-243 (2006).
2. Ye, C., Wang, E., He, S., *et al.* Subchronic toxicity and hepatocyte apoptosis of dietary olaquinox in common carp (*Cyprinus carpio*). *Ecotoxicol. Environ. Saf.* **164**, 131-139 (2018).
3. Kang, H.G., Ku, H.O., Jeong, S.H., *et al.* Evaluation of embryotoxic potential of olaquinox and vitamin A in micromass culture and in rats. *Toxicol. Res.* **26**(3), 209-216 (2010).

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, 11/14/2022

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