

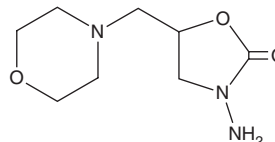
PRODUCT INFORMATION



AMOZ

Item No. 36480

CAS Registry No.: 43056-63-9
Formal Name: 3-amino-5-(4-morpholinylmethyl)-2-oxazolidinone
Synonym: 3-Amino-5-morpholinomethyl-2-oxazolidinone
MF: $C_8H_{15}N_3O_3$
FW: 201.2
Purity: $\geq 95\%$
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

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

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of AMOZ can be prepared by directly dissolving the solid in aqueous buffers. The solubility of AMOZ in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

AMOZ is a metabolite of the nitrofuran antibiotic furaltadone.¹

Reference

1. Verdon, E., Couedor, P., and Sanders, P. Multi-residue monitoring for the simultaneous determination of five nitrofurans (furazolidone, furaltadone, nitrofurazone, nitrofurantoin, nifursol) in poultry muscle tissue through the detection of their five major metabolites (AOZ, AMOZ, SEM, AHD, DNSAH) by liquid chromatography coupled to electrospray tandem mass spectrometry--in-house validation in line with commission decision 657/2002/EC. *Anal. Chim. Acta* **586(1-2)**, 336-347 (2007).

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