

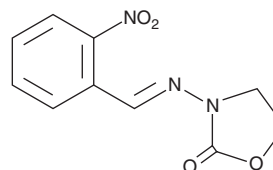
PRODUCT INFORMATION



2-NP-AOZ

Item No. 35575

CAS Registry No.: 19687-73-1
Formal Name: 3-[[[2-nitrophenyl)methylene]amino]-2-oxazolidinone
Synonym: NPAOZ
MF: C₁₀H₉N₃O₄
FW: 235.2
Purity: ≥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

2-NP-AOZ is supplied as a solid. A stock solution may be made by dissolving the 2-NP-AOZ in the solvent of choice, which should be purged with an inert gas. 2-NP-AOZ is slightly soluble in chloroform and methanol.

Description

2-NP-AOZ is a derivatized metabolite of the nitrofurantolone antiprotozoal and antibacterial agent furazolidone (Item No. 34090).¹ It is formed from the tissue-bound furazolidone metabolite 3-amino-2-oxazolidinone (AOZ) after acid hydrolysis and derivatization with *o*-nitrobenzaldehyde (*o*-NBA). 2-NP-AOZ has been used as a marker of furazolidone use in pigs.²

References

1. Cheng, C.-C., Hsieh, K.-H., Lei, Y.-C., *et al.* Development and residue screening of the furazolidone metabolite, 3-amino-2-oxazolidinone (AOZ), in cultured fish by an enzyme-linked immunosorbent assay. *Comparative Study* **57(13)**, 5687-5692 (2009).
2. Hoogenboom, L.A., Berghmans, T.H., Polman, T.H., *et al.* Depletion of protein-bound furazolidone metabolites containing the 3-amino-2-oxazolidinone side-chain from liver, kidney and muscle tissues from pigs. *Food Addit. Contam.* **9(6)**, 623-630 (1992).

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, 12/12/2022

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