

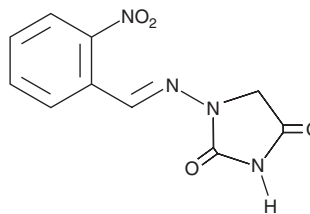
# PRODUCT INFORMATION



## 2-NP-AHD

Item No. 36500

**CAS Registry No.:** 623145-57-3  
**Formal Name:** 1-[[[(2-nitrophenyl)methylene]amino]-2,4-imidazolidinedione  
**MF:** C<sub>10</sub>H<sub>8</sub>N<sub>4</sub>O<sub>4</sub>  
**FW:** 248.2  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 267 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

2-NP-AHD is supplied as a solid. A stock solution may be made by dissolving the 2-NP-AHD in the solvent of choice, which should be purged with an inert gas. 2-NP-AHD is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

### Description

2-NP-AHD is a nitrophenyl-labeled derivative of 1-aminohydantoin (AHD), a metabolite of the nitrofuran antibiotic nitrofurantoin (Item No. 23510).<sup>1</sup> It has been used in immunoassays for the detection of AHD in honey.

### Reference

1. Li, Z., Li, Z., and Xu, D. Simultaneous detection of four nitrofuran metabolites in honey by using a visualized microarray screen assay. *Food Chem.* **221**, 1813-1821 (2017).

#### 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