

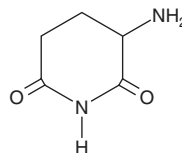
PRODUCT INFORMATION



3-Amino-2,6-Piperidinedione

Item No. 31698

CAS Registry No.: 2353-44-8
Synonyms: α -Aminoglutarimide,
3-Aminoglutarimide, Glutamimide
MF: $C_5H_8N_2O_2$
FW: 128.1
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

3-Amino-2,6-piperidinedione is supplied as a crystalline solid. Aqueous solutions of 3-amino-2,6-piperidinedione can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 3-amino-2,6-piperidinedione in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

3-Amino-2,6-piperidinedione is an active metabolite of (\pm)-thalidomide (Item No. 14610).¹ *In vivo*, 3-amino-2,6-piperidinedione (400 mg/kg) increases hexobarbitone-induced sleeping time in rats.

Reference

1. Fabro, S., Schumacher, H., Smith, R.L., *et al.* The metabolism of thalidomide: Some biological effects of thalidomide and its metabolites. *Br. J. Pharmacol. Chemother.* **25(2)**, 352-362 (1965).

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, 11/02/2022

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