

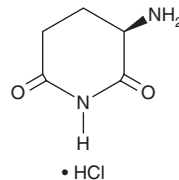
PRODUCT INFORMATION



D-2-Aminoglutarimide (hydrochloride)

Item No. 31697

CAS Registry No.: 1801140-47-5
Formal Name: (3R)-3-amino-2,6-piperidinedione, monohydrochloride
MF: C₅H₈N₂O₂ • HCl
FW: 164.6
Purity: ≥95%
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

D-2-Aminoglutarimide (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the D-2-aminoglutarimide (hydrochloride) in the solvent of choice, which should be purged with an inert gas. D-2-Aminoglutarimide (hydrochloride) is soluble in the organic solvent methanol.

Description

D-2-Aminoglutarimide is a synthetic intermediate.¹ It has been used in the synthesis of hydantoins with anti-angiogenic and cancer cell differentiation-inducing activities.

Reference

1. Yamaguchi, J., Noguchi-Yachide, T., Sakaguchi, Y., *et al.* Synthesis of new hydantoins bearing glutarimide or succinimide moiety and their evaluation for cell differentiation-inducing and anti-angiogenic activities. *Heterocycles* **91(4)**, 764-781 (2015).

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, 11/10/2022

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