

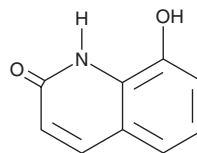
PRODUCT INFORMATION



2,8-Quinolinediol

Item No. 36376

CAS Registry No.: 15450-76-7
Formal Name: 8-hydroxy-2(1H)-quinolinone
Synonyms: 2,8-Dihydroxylquinoline, 2,8-Dihydroxyquinoline, 8-Hydroxycarbostyryl, NSC 108383
MF: C₉H₇NO₂
FW: 161.2
Purity: ≥98%
UV/Vis.: λ_{max}: 234, 258 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

2,8-Quinolinediol is supplied as a solid. A stock solution may be made by dissolving the 2,8-quinolinediol in the solvent of choice, which should be purged with an inert gas. 2,8-Quinolinediol is soluble in the organic solvent DMSO at a concentration of approximately 2 mg/ml.

Description

2,8-Quinolinediol is an endogenous metabolite.¹ Serum levels of 2,8-quinolinediol are increased in hemodialysis patients with mild cognitive decline.

Reference

1. Zhu, B., Shen, J., Jiang, R., *et al.* Abnormalities in gut microbiota and serum metabolites in hemodialysis patients with mild cognitive decline: A single-center observational study. *Psychopharmacology (Berl)* 237(9), 2739-2752 (2020).

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, 02/02/2023

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