

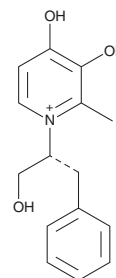
PRODUCT INFORMATION



CN128

Item No. 30833

CAS Registry No.: 2457170-98-6
Formal Name: 3,4-dihydroxy-1-[(1R)-1-(hydroxymethyl)-2-phenylethyl]-2-methyl-pyridinium
Synonym: R-CN128
MF: C₁₅H₁₈NO₃
FW: 260.3
Purity: ≥95%
UV/Vis.: λ_{max}: 287 nm
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

CN128 is supplied as a crystalline solid. A stock solution may be made by dissolving the CN128 in the solvent of choice, which should be purged with an inert gas. CN128 is soluble in organic solvents such as methanol. The solubility of CN128 in methanol is approximately 5 mg/ml.

CN128 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CN128 should first be dissolved in methanol and then diluted with the aqueous buffer of choice. CN128 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of methanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CN128 is an orally bioavailable iron chelator.¹ It contains a side chain hydroxy group that acts as an alternate sacrificial glucuronidation site to alleviate metabolic inactivation at the 3-hydroxy group. CN128 increases iron mobilization by 24.8% in a ⁵⁹Fe-ferritin-loaded rat model of iron overload when administered at doses of 150 and 450 μmol/kg.

Reference

1. Chen, W., Yuan, X., Li, Z., *et al.* CN128: A new orally active hydroxypyridinone iron chelator. *J. Med. Chem.* **63**(8), 4215-4226 (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, 11/07/2025

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