

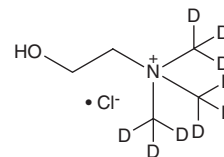
PRODUCT INFORMATION



Choline-d₉ (chloride)

Item No. 36454

CAS Registry No.: 61037-86-3
Formal Name: 2-hydroxy-N,N,N-tri(methyl-d₃)-ethanaminium, monochloride
MF: C₅H₅D₉NO • Cl
FW: 148.7
Chemical Purity: ≥98% (Choline)
Deuterium
Incorporation: ≥99% deuterated forms (d₁-d₉); ≤1% d₀
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

Choline-d₉ (chloride) is intended for use as an internal standard for the quantification of choline (Item No. 31178) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Choline-d₉ (chloride) is supplied as a solid. A stock solution may be made by dissolving the choline-d₉ (chloride) in the solvent of choice, which should be purged with an inert gas. Choline-d₉ (chloride) is soluble in the organic solvent ethanol at a concentration of approximately 2 mg/ml. Choline-d₉ (chloride) is also slightly soluble in DMSO.

Description

Choline is an essential nutrient with roles in liver, neurological, hematological, and skeletal muscle homeostasis.¹⁻⁴ It is a precursor in the biosynthesis of membrane phospholipids, such as phosphatidylcholine, which facilitate cell signaling and transport across the membrane, and a precursor to the neurotransmitter acetylcholine. Choline is required for hepatic lipid transport.¹ Perinatal administration of choline (18.8 mg/kg) improves prenatal alcohol exposure-induced cognitive deficits in rats.² Choline (13 mg/animal per day) improves motor coordination and behavioral deficits in a mouse model of Rett syndrome, as well as improves deficits in recognition memory induced by early-life iron deficiency in rats when administered in the drinking water at a concentration of 5 ppm.³ Deficiencies in choline intake are positively correlated with muscle wasting, and dietary administration of choline (1,000 mg/kg) increases leg and breast muscle protein content in broiler chickens.⁴

References

1. Sanders, L.M. and Zeisel, S.H. Choline: Dietary requirements and role in brain development. *Nutr. Today* **42(4)**, 181-186 (2007).
2. Thomas, J.D., Garrison, M., and O'Neill, T.M. Perinatal choline supplementation attenuates behavioral alterations associated with neonatal alcohol exposure in rats. *Neurotoxicol. Teratol.* **26(1)**, 35-45 (2004).
3. Derbyshire, E. and Obeid, R. Choline, neurological development and brain function: A systematic review focusing on the first 1000 days. *Nutrients* **12(6)**, 1731 (2020).
4. Moretti, A., Paoletta, M., Liguori, S., *et al.* Choline: An essential nutrient for skeletal muscle. *Nutrients* **12(7)**, E2144 (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, 10/26/2022

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