

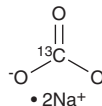
PRODUCT INFORMATION



Sodium carbonate-¹³C

Item No. 44118

CAS Registry No.: 93673-48-4
Formal Name: carbonic-¹³C acid, disodium salt
MF: [¹³C]O₃ • 2Na
FW: 44118
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

Sodium carbonate-¹³C is intended for use as an internal standard for the quantification of sodium carbonate 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).

Description

Sodium carbonate is an alkalinizing agent.¹⁻³ It has been used in the development of alkaline flooding systems for crude oil recovery, to improve the digestibility of lignocellulosic biomass, and in the purification of biogases.

References

1. Sun, J., Sun, L., Liu, W., *et al.* Alkaline consumption mechanisms by crude oil: A comparison of sodium carbonate and sodium hydroxide. *Physiochem. Eng. Aspects* **315**, 38-43 (2008).
2. Mirmohamadsadeghi, S., Chen, Z., and Wan, C. Reducing biomass recalcitrance via mild sodium carbonate pretreatment. *Bioresour. Technol.* **209**, 386-390 (2016).
3. Murialdo, M., H.M., G., Stolaroff, J.K., *et al.* Three-dimensional printable sodium carbonate composite sorbents for efficient biogas upgrading. *Environ. Sci. Technol.* **54(11)**, 6900-6907 (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, 02/06/2026

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