

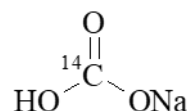
Sodium Bicarbonate, [¹⁴C]-

Product Number:

NEC086H

LOT SPECIFIC INFORMATION

Lot Number:	3250554
Specific Activity:	58 mCi/mmol
	2146 MBq/mmol
Production Date:	07-FEB-2024

M.W. 84.0
CHNaO₃

PACKAGING: 1.0 mCi/ml (37 MBq/ml) steri-packaged in glass ampoules containing aqueous solution, pH 9.5.

STABILITY AND STORAGE: Dilute solutions of sodium bicarbonate, [¹⁴C]- in sterile water have been found to be extremely stable when stored at 5°C (not frozen).

SPECIFIC ACTIVITY RANGE: 40-60 mCi/mmol (1480-2220 MBq/mmol)

IMPORTANT: Since the sodium bicarbonate, [¹⁴C]- solutions are thermally sensitive, the following procedure for using the product is recommended:

- 1) Shake ampoule just prior to opening.
- 2) Gently tap ampoule to remove any solution remaining in the head.
- 3) Open ampoule while still cold.

FAILURE TO FOLLOW THE ABOVE DIRECTIONS MAY RESULT IN THE INCOMPLETE RECOVERY OF SODIUM BICARBONATE, [¹⁴C]-

NOTE: Ensure dilution solvent is basic, and previously degassed with an inert gas. If solution needs to be diluted, the bulk must be stored in an air-tight closed vial under inert gas. Aliquoting the solution into an acidic solvent will result in significant radioactivity loss as gaseous carbon dioxide [¹⁴C]-.

HAZARD INFORMATION: WARNING: This product contains a chemical known to the state of California to cause cancer.

Limited Warranty: The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

NEC086H-R Rev01

revvity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA

(800) 762-4000
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.