

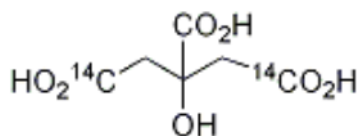
Citric Acid, [1,5-¹⁴C]-

Product Number:

NEC160

LOT SPECIFIC INFORMATION

Lot Number:	3157247
Specific Activity:	116.4 mCi/mmol
	4306.8 MBq/mmol
Production Date:	9 Jun 2023



C₆H₈O₇
M.W. 192.12
CAS 77-92-9

PACKAGING: 0.05 mCi/mL (1.85 MBq/mL) in ethanol : water (2:98), Steri-packaged. Shipped in blue ice.

STABILITY AND STORAGE: When citric acid, [1,5-¹⁴C]- is stored at 5°C in its original solvent and at its original concentration, the rate of decomposition is initially less than 3% per year from the date of purification. The rate of decomposition can accelerate. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.

SPECIFIC ACTIVITY RANGE: 100-124 mCi/mmol (3700-4588 MBq/mmol)

RADIOCHEMICAL PURITY: This product was initially found to be greater than 97% when determined by the following method. It is advisable to check purity prior to use:

High pressure liquid chromatography on an Aminex HPX-87H column using the following mobile phase:

0.01N H₂SO₄

QUALITY CONTROL: The radiochemical purity of citric acid, [1,5-¹⁴C]- is checked at appropriate intervals using the above chromatography method.

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.

NEC160-R Rev01

The Revvity logo consists of the word "revvity" in a lowercase, sans-serif font. The letters are black and have a modern, clean appearance.

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.