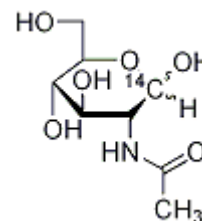


N-Acetyl-D- [1-¹⁴C]- Glucosamine

Product Number: NEC723V

LOT SPECIFIC INFORMATION

Lot Number:	3152268
Specific Activity:	58.1 mCi/mmol
	2149.7 MBq/mmol
Production Date:	4 May 2023



M.W. 221.2
C₈H₁₅NO₆
CAS 7512-17-6

PACKAGING: 0.2 mCi/ml (7.40 MBq/ml) in ethanol : water (2:98), steri-packaged. Shipped in dry ice.

STABILITY AND STORAGE RECOMMENDATIONS: When N-acetyl-D- [1-¹⁴C]- glucosamine is stored at -20°C in its original solvent and at its original concentration, the rate of decomposition is initially 1% per year from 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: 50-62 mCi/mmol (1850-2294 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-87C column at 80°C using the following mobile phase:
Water

QUALITY CONTROL: The radiochemical purity of N-acetyl-D- [1-¹⁴C]- glucosamine is checked at appropriate intervals using the above chromatography method.

SPECIAL INFORMATION:

Removal of solvent: An aliquot or the entire sample can be taken to dryness by directing a gentle stream of inert gas (nitrogen) over the surface of the solution. The temperature of the solution should not be allowed to exceed 30°C during the drying process, and the compound should not be permitted to remain in the solid state any longer than necessary.

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.

NEC723V-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.