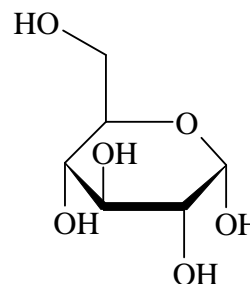


Glucose, D- $^{14}\text{C}(\text{U})$ -

Product Number: NEC042X

LOT SPECIFIC INFORMATION

Lot Number:	2905027
Specific Activity:	265 mCi/mmol
	9805 MBq/mmol
Production Date:	19 Oct 2021

M.W. 180.16  
 $\text{C}_6\text{H}_{12}\text{O}_6$ **PACKAGING:** 0.1 mCi/ml (3.7 MBq/ml) in ethanol : water solution, 9:1.**STABILITY AND STORAGE:** When NEC-042X glucose, D- $^{14}\text{C}(\text{U})$ - is stored at 5°C in its original solvent and at its original concentration, the rate of decomposition is initially 1% for the first 12 months from date of purification. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.**SPECIFIC ACTIVITY RANGE:** 250-360 mCi/mmol (9250-13320 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 Chromotography on an Aminex HPX-87C column using the following mobile phase:  
Water at 85°C.**QUALITY CONTROL:** The radiochemical purity of glucose, D- $^{14}\text{C}(\text{U})$ - is checked at appropriate intervals using the first listed 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.

NEC042X-R Rev01

revvity

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

(800) 762-4000  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.