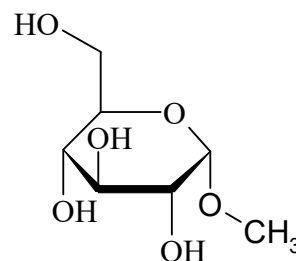


Methyl  $\alpha$ -D-Glucopyranoside, [Glucose-<sup>14</sup>C(U)]-

Product Number: NEC659V

**LOT SPECIFIC INFORMATION**

|                    |               |
|--------------------|---------------|
| Lot Number:        | 3243241       |
| Specific Activity: | 270 mCi/mmol  |
|                    | 9990 MBq/mmol |
| Production Date:   | 10-Jan-2024   |

M.W. 194.2  
C<sub>7</sub>H<sub>14</sub>O<sub>6</sub>**PACKAGING:** 0.2 mCi/ml (7.4 MBq/ml) in ethanol:water (3:97), steri-packaged. Shipped on dry ice.**STABILITY AND STORAGE RECOMMENDATIONS:**

When methyl  $\alpha$ -D-glucopyranoside, [glucose-<sup>14</sup>C(U)]- is stored at -20°C in its original solvent and at its original concentration, the rate of decomposition is initially 2% 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:** 230-370 mCi/mmol (8510-13690 MBq/mmol)**RADIOCHEMICAL PURITY:** This product was initially found to be greater than 97% when determined by the following methods. 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 methyl  $\alpha$ -D-glucopyranoside, [glucose-<sup>14</sup>C(U)]- is checked at appropriate interval using the HPLC 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.

NEC659V-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](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.