

Methyl-D-Glucose, 3-O-[Methyl-¹⁴C]-

Product Number: NEC377

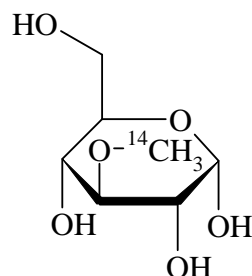
LOT SPECIFIC INFORMATION

Lot Number: 2706926

Specific Activity: 56.4 mCi/mmol

2087 MBq/mmol

Production Date: 23 Mar 2021



M.W. 194.2

C₇H₁₄O₆**PACKAGING:** 0.1 mCi/ml (3.7 MBq/ml) in ethanol : water solution, 9:1.**STABILITY AND STORAGE:** When methyl-D-glucose, 3-O-[methyl-¹⁴C]- is stored at -20°C in its original solvent and at its original concentration, the rate of decomposition is initially 0.5% per year from date of purification. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.**SPECIFIC ACTIVITY RANGE:** 30-60 mCi/mmol (1110-2220 MBq/mmol)**RADIOCHEMICAL PURITY:** This product was initially found to be greater than 97% when determined by the following method. The rate of decomposition can accelerate. It is advisable to check purity prior to use:High performance liquid chromatography on an Aminex HPX-87C column using the following mobile phase:
Water

Paper chromatography on Whatman No. 1 using the following solvent systems:

- n-butanol : ethanol : water, (50:32:18).
- n-butanol : pyridine : water, (6:4:3).
- n-butanol : acetic acid : water, (4:1:5).

QUALITY CONTROL: The radiochemical purity of methyl-D-glucose, 3-O-[methyl-¹⁴C]- is checked at appropriate intervals 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.

NEC377-R Rev01

revvity