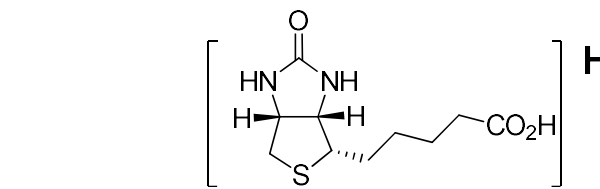


Biotin, d-[³H(G)]-

Product Number: NET1215

LOT SPECIFIC INFORMATION

Lot Number:	3309480
Specific Activity:	14.6 Ci/mmol
	540.2 GBq/mmol
Production Date:	26-Jun-2024



M.W. 244.3
C₁₀H₁₆N₂O₃S

PACKAGING: 1.0 mCi/ml (37 MBq/ml) in aqueous solution containing 50 micromoles of 2-mercaptoethanol per ml, under argon, sterilized by filtration and dispensed under aseptic conditions. Shipped in dry ice.

STABILITY AND STORAGE RECOMMENDATIONS:

- When Biotin, d-[³H(G)]- is stored at -80°C in its original solvent and at its original concentration, the rate of decomposition is initially less than 1% per month from date of purification. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.
- There is a tendency for labeled compounds containing sulfur to undergo oxidation at the sulfur atom. Repeated cycles of thawing and freezing may increase the rate of oxidation. It is recommended that if the entire quantity cannot be used at once the appropriate size aliquots be made and refrozen at -80°C until needed. Storage under argon is recommended.

SPECIFIC ACTIVITY RANGE: >10 Ci/mmol (>370 GBq/mmol)

RADIOCHEMICAL PURITY: This product was initially found to be greater than 95% when determined by the following method. The rate of decomposition can accelerate. It is advisable to check purity prior to use:

High pressure liquid chromatography on a reverse phase column using the following mobile phase:
0.2% trifluoroacetic acid in water : 0.01% trifluoroacetic acid in acetonitrile, (85:15) isocratic for 20 minutes.

QUALITY CONTROL: The radiochemical purity of Biotin, d-[³H(G)]- 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.

NET1215-R Rev01

revvity