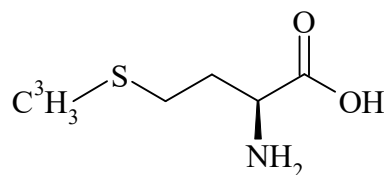


METHIONINE, L-[METHYL- ^3H]-

Product Number: NET061X

LOT SPECIFIC INFORMATION

Lot Number:	3187741
Specific Activity:	82.8 Ci/mmol
	3063.6 GBq/mmol
Production Date:	6-Nov-2023

M.W. 149
 $\text{C}_5\text{H}_{11}\text{NO}_2\text{S}$

PACKAGING: 1.0 mCi/ml (37 MBq/ml) in 50mM Tricine, pH 7.4, containing 10mM 2-mercaptoethanol, under argon. Shipped in dry ice.

STABILITY AND STORAGE RECOMMENDATIONS: When methionine, L-[methyl- ^3H]- is stored at -80°C in its original solvent and at its original concentration, the rate of decomposition is initially 3% per month from date of purification. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.

- Methionine sulfoxide is the major decomposition product.

SPECIFIC ACTIVITY RANGE: 70-85 Ci/mmol (2590-3145 GBq/mmol)

RADIOCHEMICAL PURITY: This product was initially found to be greater than 97% when determined by one or more of the following methods. The rate of decomposition can accelerate. It is advisable to check purity prior to use:

High pressure liquid chromatography on a Zorbax SCX column using the following mobile phase:
potassium phosphate, pH 3.0.

QUALITY CONTROL: The radiochemical purity of methionine, L-[methyl- ^3H]- is checked at appropriate intervals using the listed chromatography method.

SPECIAL INFORMATION: Labeled compounds of high specific activity are often stored in a solvent to inhibit decomposition due to radiolysis. If it is desirable to remove the solvent before use, only an appropriate aliquot should be taken to dryness and the remainder stored in solution as recommended above.

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.

NET061X-R Rev01

revvity