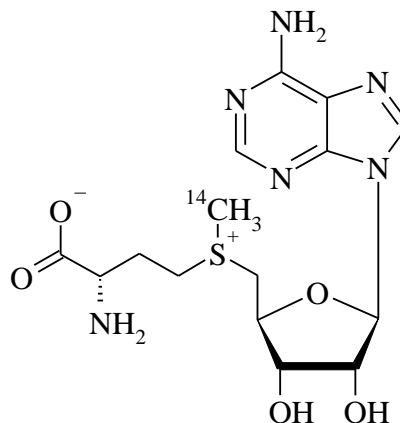


Adenosyl-L-Methionine, S-[Methyl-<sup>14</sup>C]-

Product Number: NEC363

LOT SPECIFIC INFORMATION

Lot Number:	3154705
Specific Activity:	51.4 mCi/mmol
	1901.8 MBq/mmol
Production Date:	6 Jun 2023



M.W. 398.4  
 $C_{15}H_{22}N_6O_5S$

**PACKAGING:** 0.02 mCi/ml (0.740 MBq/ml) in 10mM sulfuric acid : ethanol (9:1). Shipped in dry ice.

**STABILITY AND STORAGE:** When adenosyl-L-methionine, S-[methyl-<sup>14</sup>C]- is stored at -20°C in its original solvent and at its original concentration, the rate of decomposition is initially 1% for 6 months from date of purification. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.

- This product is unstable when stored at neutral or basic pH (1) and is best stored as supplied.
- Repeated freezing and thawing can result in product degradation. If the entire quantity cannot be used at once, it is recommended that the product be slowly thawed at 4°C, aliquoted under N<sub>2</sub> into samples of an appropriate size and then refrozen at -20°C.

**SPECIFIC ACTIVITY RANGE:** 40-60 mCi/mmol (1480-2220 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 pressure liquid chromatography on a strong cation exchange (SCX) column using the following mobile phase:

ammonium formate buffer, pH 2.5.

**QUALITY CONTROL:** The radiochemical purity of adenosyl-L-methionine, S-[methyl-<sup>14</sup>C]- 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.

NEC363-R Rev01

revvity