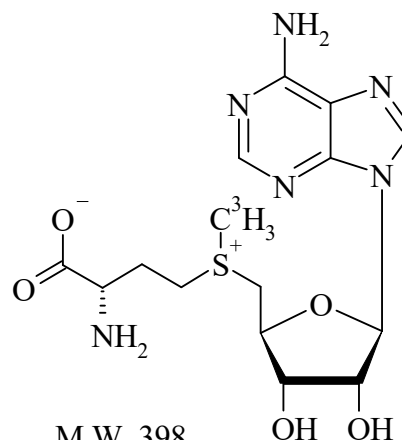


Adenosyl-L-Methionine, S-[Methyl-<sup>3</sup>H]-

Product Number: NET155

LOT SPECIFIC INFORMATION

Lot Number:	3217608
Specific Activity:	13.9 Ci/mmol
	514.3 GBq/mmol
Production Date:	09-Nov-2023



M.W. 398  
C<sub>15</sub>H<sub>22</sub>N<sub>6</sub>O<sub>5</sub>S

**PACKAGING:** 0.55 mCi/ml (20.35 MBq/ml) in 10mM sulfuric acid:ethanol (9:1), under argon. Shipped in dry ice.

**STABILITY AND STORAGE RECOMMENDATIONS:**

- When adenosyl-L-methionine, S-[methyl-<sup>3</sup>H]- is stored at -20°C in its original solvent and at its original concentration, the rate of decomposition is initially 1.5 % per month from date of purification. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.
- Storage at -20°C is recommended. The rate of decomposition increases when adenosyl-L-methionine, S-[methyl-<sup>3</sup>H]- is stored at -80°C.
- To maintain product purity, this sample should be slow thawed at 4°C. Repeated freezing and thawing should be avoided. Store under argon or other inert gas.
- Adenosyl-L-methionine, S-[methyl-<sup>3</sup>H]- is unstable at neutral or basic pH.

**SPECIFIC ACTIVITY RANGE:** 5-15 Ci/mmol (185-555 GBq/mmol)

**ASSAY PROCEDURE:** All samples should be counted in 2.5% Solvable™/Formula 989. Samples for assay are diluted in 0.1 N HCl. in plastic vials. Aquasol® is not recommended.

**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 column using the following mobile phase:  
0.25M ammonium formate, pH 2.5.

**QUALITY CONTROL:** The radiochemical purity of adenosyl-L-methionine, S-[methyl-<sup>3</sup>H]- is checked at appropriate intervals using the 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.

NET155-R Rev01

revvity