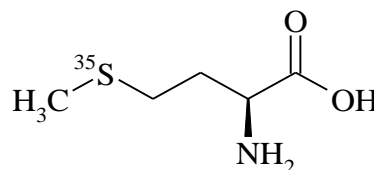


Methionine, L-[³⁵S]-, Cell Labeling Grade

Product Number: NEG009L

LOT SPECIFIC INFORMATION

Lot Number:	04154
Specific Activity:	1165.5 Ci/mmol 43.5 TBq/mmol
Concentration:	31.5 mCi/ml 1165.5 MBq/ml 0.0268 μmol/ml 0.0040 mg/ml
Calibration Date:	10-May-2024



M.W. 149.2
C₅H₁₁NO₂S

PACKAGING: Aqueous solution containing 10mM 2-mercaptoethanol in a NENSURE vial with polypropylene-v-insert, shipped in dry ice.

PRODUCT SPECIFICATIONS: NEG-009L, Methionine, L-[³⁵S]- Cell Labeling Grade is obtained from the protein fraction of bacteria grown in the presence of carrier-free [³⁵S] sulfate. NEG-009L contains only ³⁵S-labeled L-methionine and its oxidation products, L-[³⁵S] methionine sulfoxide and L-[³⁵S] methionine sulfone. All traces of L-[³⁵S] cysteine and other ³⁵S radiolabeled contaminants are removed during processing of hydrolysate.

STABILITY AND STORAGE:

- The rate of decomposition of NEG-009L, Methionine, L-[³⁵S]- Cell Labeling Grade is approximately 8-10% per week when stored at -80°C. Storage at higher temperatures accelerates product decomposition.
- If the entire quantity will not be used at once it is recommended that the product be thawed quickly, aliquoted under N₂ into samples of an appropriate size, refrozen immediately and stored at -20°C or below.
- Methionine, L-[³⁵S] is highly susceptible to oxidation and should always be protected from exposure to air.

HAZARD INFORMATION: WARNING: This product contains a chemical known to the state of California to cause cancer.

QUALITY CONTROL:

Radiochemical Purity: This lot was initially found to be >95% when determined by high performance liquid chromatography on a C18 column.

Chemical Purity: As determined by amino acid analysis, NEG-009A is routinely >99% L-isomer and contains the unlabeled amino acids alanine and valine. Each is present in approximately a 2:1 molar ratio relative to Methionine, L-[³⁵S].

PRODUCT APPLICATION: NEG-009L, Methionine, L-[³⁵S] is intended for use in eukaryotic or bacterial metabolic labeling studies. It is not recommended or warranted for use in cell-free mRNA-directed translations. The use of NEG-009L in the latter application produces irreproducible experimental results, with significant variability in both incorporation and background levels.

SAFE HANDLING: Because this product has been stored at -80°C, it is possible that pressure may develop in the vial during the thawing process. In addition, volatile ³⁵S-labeled decomposition products are generated at a

rate of ~0.01% per week. We recommend that prior to opening, vials are first vented in a fume hood using the following procedure:

- Slide aside the dust cover on the cap to expose the septum.
- Pierce the septum with a cotton-plugged syringe needle or charcoal trap (NENTM NEX-033T), taking care that the tip does not come into contact with the product.
- If the product is frozen, quickly thaw at room temperature or in a 37° C water bath. Any pressure developed will vent through the syringe needle.
- Remove the needle and dispose of as contaminated equipment.

When used for *in vivo* labeling experiments, we recommend that specific steps be taken to minimize incubator and water bath contamination. We suggest using a shallow tray of activated charcoal, charcoal sticks or charcoal filter units to trap ³⁵S volatiles and reduce contamination.

SPECIAL INFORMATION: Visit www.revvy.com to use our online Radioactive Decay Calculator.

Decay of ³⁵S (physical half-life, 87.4 days):

		DECAY FACTORS									
Days BEFORE		0	1	2	3	4	5	6	7	8	9
Assay Date											
30		1.269	1.279	1.289	1.299	1.309	1.320	1.330	1.341	1.352	1.362
20		1.172	1.181	1.191	1.200	1.210	1.219	1.229	1.239	1.249	1.259
10		1.083	1.091	1.100	1.109	1.117	1.135	1.135	1.144	1.153	1.163
0		1.000	1.008	1.016	1.024	1.032	1.049	1.049	1.057	1.066	1.074
Days AFTER		0	1	2	3	4	5	6	7	8	9
Assay Date											
0		1.000	0.992	0.984	0.976	0.969	0.961	0.954	0.946	0.939	0.931
10		0.924	0.916	0.909	0.902	0.895	0.888	0.881	0.874	0.867	0.860
20		0.853	0.847	0.840	0.833	0.827	0.820	0.814	0.807	0.801	0.795
30		0.788	0.782	0.776	0.770	0.764	0.758	0.752	0.746	0.740	0.734

The specific activity at time t (SA_t) may be calculated, using the following equation, from the specific activity at the calibration date (SA_0) and the decay factor (f) given above.

$$SA_t = \frac{f}{1/SA_0 - (1-f)/1494}$$

RELATED PRODUCTS:

- NEG009A Methionine, L-[³⁵S]-
- NEG009C Methionine, L-[³⁵S]-
- NEG022T Cysteine, L-[³⁵S]-

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.

