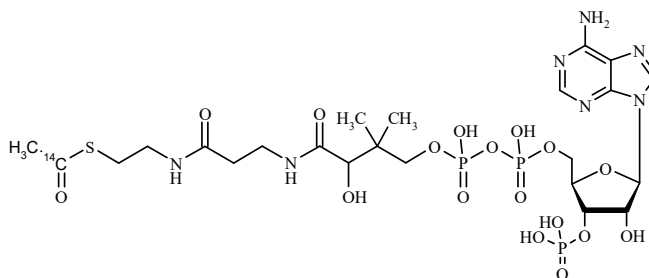


Acetyl Coenzyme A, [Acetyl-1-<sup>14</sup>C]-

Product Number: NEC313

## LOT SPECIFIC INFORMATION

|                    |               |
|--------------------|---------------|
| Lot Number:        | 3384706       |
| Specific Activity: | 58 mCi/mmol   |
|                    | 2146 MBq/mmol |
| Production Date:   | 17-Mar-2025   |



M.W. 809.6  
 $C_{23}H_{38}N_7O_{17}P_3S$

**PACKAGING:** 0.02 mCi/ml (0.74 MBq/ml) in 0.01M sodium acetate (pH 5.0), under nitrogen. Shipped in dry ice.

**STABILITY AND STORAGE:**

- When acetyl coenzyme A, [acetyl-1-<sup>14</sup>C]- is stored at -20°C in its original solvent and at its original concentration, the rate of decomposition is initially 3% over the first six months from date of purification. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.
- Optimal storage of NEC-313 is at pH 5.0, -20°C under nitrogen. This compound is extremely unstable at pH levels greater than 8.5.
- To maintain product purity, this sample should be slowly thawed at 4°C. Repeated freezing and thawing should be avoided.

**SPECIFIC ACTIVITY RANGE:** 40-60 mCi/mmol (1480-2220 MBq/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 C8 or C18 reverse phase column using the following mobile phase: 50mM potassium phosphate buffer (pH 5.3):MeOH, (85:15).

**QUALITY CONTROL:** The radiochemical purity of acetyl coenzyme A, [acetyl-1-<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.

NEC313-R Rev01

revvity