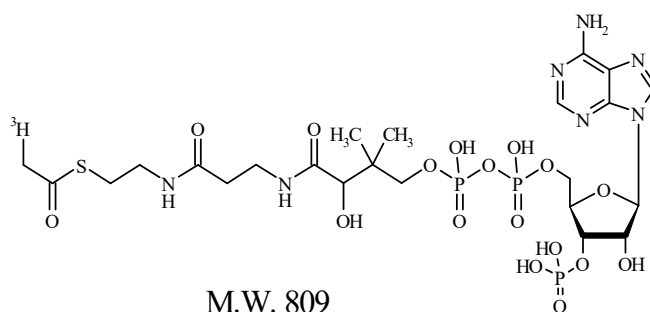


Acetyl Coenzyme A, [Acetyl- $^3\text{H}$ ]-

Product Number: NET290

## LOT SPECIFIC INFORMATION

Lot Number:	3257048
Specific Activity:	3.2 Ci/mmol
	118.4 GBq/mmol
Production Date:	10-Jul-2024

**PACKAGING:** 0.1 mCi/ml (3.7 MBq/ml) in 0.01M sodium acetate (pH 5.0), under nitrogen. Shipped in dry ice.

## STABILITY AND STORAGE RECOMMENDATIONS:

- When acetyl coenzyme A, [acetyl- $^3\text{H}$ ]- is stored at  $-20^\circ\text{C}$  in its original solvent and at its original concentration, the rate of decomposition is initially 1% per month of from date of purification. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.
- Optimal storage of NET-290 is at pH 5.0,  $-20^\circ\text{C}$ , and under  $\text{N}_2$ . This compound is extremely unstable at pH levels greater than 8.5.
- To maintain product purity, this sample should be slow thawed at  $4^\circ\text{C}$ . Repeated freezing and thawing should be avoided. If the amount will not be used at once, it is recommended that the product be thawed at  $4^\circ\text{C}$ , aliquoted into samples of an appropriate size and immediately refrozen at  $-20^\circ\text{C}$ .

**SPECIFIC ACTIVITY RANGE:** 1-10 Ci/mmol (37-370 GBq/mmol)**RADIOCHEMICAL PURITY:** This product was initially found to be greater than 95% when determined by 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 C18 column using the following mobile phase:  
10mM potassium phosphate buffer (pH 5.3) : methanol, (85:15).**QUALITY CONTROL:** The radiochemical purity of acetyl coenzyme A, [acetyl- $^3\text{H}$ ]- 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.

NET290-R Rev01

revvity