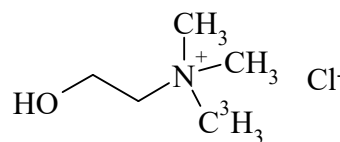


Choline Chloride, [Methyl- ^3H]-

Product Number: NET109

LOT SPECIFIC INFORMATION

Lot Number:	3147572
Specific Activity:	80.8 Ci/mmol
	2989.6 GBq/mmol
Production Date:	12-May-2023

M.W. 139.6
 $\text{C}_5\text{H}_{14}\text{ClNO}$ **PACKAGING:** 1 mCi/ml (37 MBq/ml) in ethanol, in a silanized vial. Shipped on dry ice.**STABILITY AND STORAGE RECOMMENDATIONS:** When choline chloride, [methyl- ^3H]- is stored at -20°C in its original solvent and at its original concentration, the rate of decomposition is initially less than 0.5% per month from date of purification. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.**SPECIFIC ACTIVITY RANGE:** 60-90 Ci/mmol (2220-3330 GBq/mmol)**RADIOCHEMICAL PURITY:** This product was initially found to be greater than 97% when determined by the following methods. The rate of decomposition can accelerate. It is advisable to check purity prior to use:

Thin layer chromatography on Avicel or paper chromatography on Whatman No. 1 using the following solvent system:

n-butanol : ethanol : acetic acid : water, (8:2:1:3).

QUALITY CONTROL: The radiochemical purity of choline chloride, [methyl- ^3H]- is checked at appropriate intervals using the listed chromatography method.**SPECIAL INFORMATION:** Choline chloride, [methyl- ^3H]- has been found to absorb to glass walls. For optimal assay conditions, use silanized glassware and pipets and add cold carrier prior to liquid scintillation counting.**REFERENCE:** L. T. Potter and W. Murphy, *Biochem. Pharmacol.* **16**, 1386-1388 (1967).**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.

NET109-R Rev01

revvity