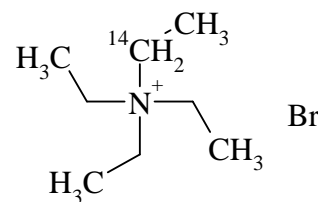


Tetraethylammonium Bromide, [$1\text{-}^{14}\text{C}$]-

Product Number: NEC298

LOT SPECIFIC INFORMATION

Lot Number:	2808844
Specific Activity:	3.5 mCi/mmol
	129.5 MBq/mmol
Production Date:	18 Dec 2020

M.W. 210.2
 $\text{C}_8\text{H}_{20}\text{BrN}$ **PACKAGING:** 0.1 mCi/ml (3.7 MBq/ml) in ethanol, in a silanized vial.**STABILITY AND STORAGE:** When tetraethylammonium bromide, [$1\text{-}^{14}\text{C}$]- is stored at 5°C in its original concentration, the rate of decomposition is initially less than 2% for 12 months from date of purification. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.**SPECIFIC ACTIVITY RANGE:** 1-5 mCi/mmol (37-185 MBq/mmol)**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:

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

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

QUALITY CONTROL: The radiochemical purity of tetraethylammonium bromide, [$1\text{-}^{14}\text{C}$]- is checked at appropriate intervals using the above 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.

NEC298-R Rev01

