

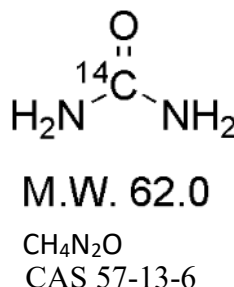
Urea, [^{14}C]-

Product Number:

NEC108V

LOT SPECIFIC INFORMATION

Lot Number:	3371936
Specific Activity:	56.5 mCi/mmol
	2090.5 MBq/mmol
Production Date:	27-Jan-2025



PACKAGING: Crystalline solid under nitrogen. Shipped ambient.

STABILITY AND STORAGE: When urea, [^{14}C]- is stored at -20°C as a dry solid, the rate of decomposition is initially less than 2% per year. The rate of decomposition can accelerate. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.

SPECIFIC ACTIVITY RANGE: 50-62 mCi/mmol (1850-2294 MBq/mmol)

RADIOCHEMICAL PURITY: This product was initially found to be greater than 97% when determined by the following methods. It is advisable to check purity prior to use:

Thin layer chromatography on a Silica Gel plate using the following mobile phase:
Chloroform:Methanol:Water(7:5:1)

Thin layer chromatography on an Avicel plate using the following mobile phase:
n-Butanol:Acetic Acid:Water (4:1:1)

High Pressure liquid chromatography on an Aminex HPX-87C column at 70°C using the following mobile phase:
Water

QUALITY CONTROL: The radiochemical purity of urea, [^{14}C]- is checked at appropriate intervals using the above chromatography methods.

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.

NEC108V-R Rev01

revvity