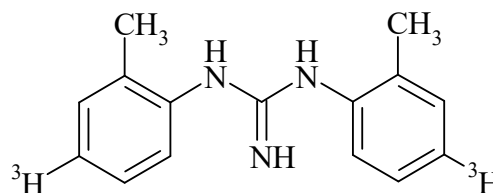


DTG, [5-³H(N)]-(1,3-Di-o-tolylguanidine, [p-ring-³H]-)

Product Number: NET986

LOT SPECIFIC INFORMATION

Lot Number:	3287393
Specific Activity:	37.8 Ci/mmol
	1398.6 GBq/mmol
Production Date:	16-May-2024



M.W. 239
C₁₅H₁₇N₃

PACKAGING: 1.0 mCi/ml (37 MBq/ml) in ethanol. Shipped in dry ice.

STABILITY AND STORAGE RECOMMENDATIONS: The stability of 1,3-di-O-tolylguanidine, di-[p-ring-³H]- is currently being evaluated. Initial studies indicate that when 1,3-di-O-tolylguanidine, di-[p-ring-³H]- is stored at -20°C in its original solvent and at its original concentration, the rate of decomposition is initially 1% for 6 months from date of purification. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.

SPECIFIC ACTIVITY RANGE: 30-60 Ci/mmol (1110-2220 GBq/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:

High pressure liquid chromatography on a Zorbax ODS column using the following mobile phase:
1% triethylamine acetate : acetonitrile, (70:30).

QUALITY CONTROL: The radiochemical purity of 1,3-di-O-tolylguanidine, di-[p-ring-³H]- 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