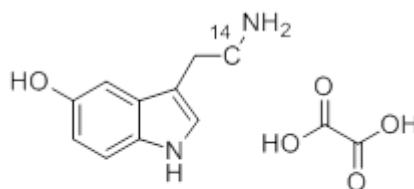


Hydroxytryptamine Binoxalate, 5-[1-¹⁴C]- (Serotonin Binoxalate)

Product Number: NEC855

LOT SPECIFIC INFORMATION

Lot Number:	3365433
Specific Activity:	51.8 mCi/mmol
	1916.6 MBq/mmol
Production Date:	19-Dec-2024



MW: 266.25
C₁₀H₁₂N₂O · C₂H₂O₄

PACKAGING: 0.1 mCi/ml (3.7 MBq/ml) in ethanol : water (2:98), under argon in a vial which protects the contents from UV light.

STABILITY AND STORAGE: When hydroxytryptamine binoxalate, 5-[1-¹⁴C]- is stored at 5°C in its original solvent and at its original concentration, the rate of decomposition is initially less than 1% for 1 month from date of purification. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.

SPECIFIC ACTIVITY RANGE: 40-60 mCi/mmol (1480-2220 MBq/mmol)

RADIOCHEMICAL PURITY: This product was initially found to be greater than 95% 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 Brownlee C18 column using the following mobile phase: methanol : 1% triethylamine acetate, pH 4.0, (5:95), 10 minutes isocratic, then 5 minute gradient to 100% methanol.

Hydroxytryptamine binoxalate, 5-[1-¹⁴C]- is susceptible to oxidation and is sensitive to light.

QUALITY CONTROL: The radiochemical purity of hydroxytryptamine binoxalate, 5-[1-¹⁴C]- 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.

NEC855-R Rev01

