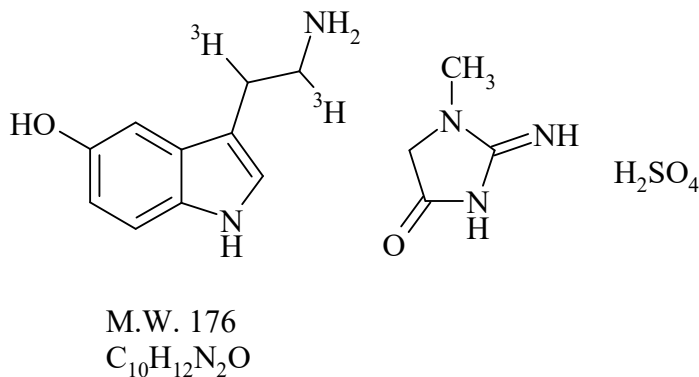


# Hydroxytryptamine Creatinine Sulfate, 5-[1,2- $^3\text{H}$ (N)]- (Serotonin)

Product Number: NET498

## LOT SPECIFIC INFORMATION

Lot Number:	3239860
Specific Activity:	32.8 Ci/mmol
	1214 GBq/mmol
Production Date:	27-Dec-2023



**PACKAGING:** 1 mCi/ml (37 MBq/ml) in ethanol:water (2:98), in a combi-vial, under argon in a vial which protects the contents from UV light. Shipped in blue ice.

**STABILITY AND STORAGE RECOMMENDATIONS:** When hydroxytryptamine creatine sulfate, 5-[1,2- $^3\text{H}$ (N)]- is stored at 5°C in its original solvent and at its original concentration, the rate of decomposition is initially 2% per month for the first three months, after which time the decomposition may accelerate rapidly. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.

**SPECIFIC ACTIVITY RANGE:** >15 Ci/mmol (>555 GBq/mmol)

**RADIOCHEMICAL PURITY:** This product was initially found to be greater than 95% when determined by one of the following method. The rate of decomposition can accelerate. It is advisable to check purity prior to use:

High pressure liquid chromatography on C18 column using the following mobile phase:  
1% triethylamine acetate, pH4 : methanol, (95:5).

**QUALITY CONTROL:** The radiochemical purity of hydroxytryptamine creatine sulfate, 5-[1,2- $^3\text{H}$ (N)]- 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.

NET498-R Rev01

revvity

Revvity, Inc.  
940 Winter Street  
Waltham, MA 02451 USA

(800) 762-4000  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.