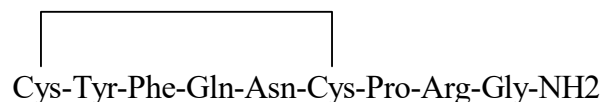


Vasopressin (8-L-Arginine), [Phenylalanyl-3,4,5- $^3\text{H}(\text{N})$ ]-

Product Number: NET800

**LOT SPECIFIC INFORMATION**

Lot Number:	3274567
Specific Activity:	69.2 Ci/mmol
	2560 GBq/mmol
Production Date:	26-Mar-2024



M. W. 1084.4

**PACKAGING:** 1.0 mCi/ml (37 MBq/ml) in ethanol : 0.05N acetic acid (7:3), in a silanized vial, under argon in a vial which protects the contents from UV light. Shipped in dry ice.

**STABILITY AND STORAGE RECOMMENDATIONS:** When vasopressin (8-L-arginine), [phenylalanyl-3,4,5- $^3\text{H}(\text{N})$ ]- is stored at -20°C in its original solvent and at its original concentration, the rate of decomposition is initially 1-2% per month from date of purification. Stability is nonlinear and not correlated to isotope half-life. Lot to lot variation may occur.

- The material is sensitive to light and air.

**SPECIFIC ACTIVITY RANGE:** 35-85 Ci/mmol (1295-3145 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 Vydac C18 column using the following mobile phase:  
0.2% trifluoroacetic acid : acetonitrile, (85:15).

**QUALITY CONTROL:** This radiochemical purity of vasopressin (8-L-arginine), [phenylalanyl-3,4,5- $^3\text{H}(\text{N})$ ]- 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