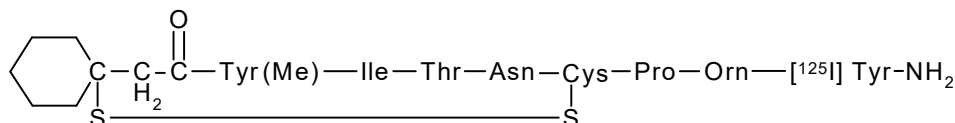


**[<sup>125</sup>I]-ORNITHINE VASOTOCIN ANALOG**

Product Number: NEX254

Vasotocin, d(CH<sub>2</sub>)<sub>5</sub>[Tyr(Me)<sup>2</sup>, Thr<sup>4</sup>, Orn<sup>8</sup>, [<sup>125</sup>I]Tyr<sup>9</sup>-NH<sub>2</sub>]-**[<sup>125</sup>I]-OVTA****LOT SPECIFIC INFORMATION**

**CALCULATED AS OF:** 13-May-2024

**LOT NUMBER:** EP71240

**SPECIFIC ACTIVITY:** 81.4 TBq/mmol  
2200 Ci/mmol  
63.7 MBq/μg  
1723 μCi/μg

**Package Size Information**

Package Size as of 12-Jul-2024
370 kBq 10 μCi
1.85 MBq 50 μCi

**RADIOCHEMICAL PURITY:** ≥ 95%**MOLECULAR WEIGHT:** ~1277

**PACKAGING:** [<sup>125</sup>I]-OVTA is lyophilized from a solution containing 0.04M sodium phosphate, 1M glycine, 0.2M NaCl, 0.25% BSA, 500 KIU/ml Trasylol® at pH 7.2. It is shipped ambient.

**STABILITY AND STORAGE:** The lyophilized [<sup>125</sup>I]-OVTA should be stored at 4°C or lower. Following reconstitution with distilled water to a concentration of approximately 50 μCi/ml on calibration date, aliquot and store at -20°C or lower. Under these conditions the product is stable and usable for at least ten weeks after fresh lot date.

**SPECIFIC ACTIVITY:** The initial specific activity of [<sup>125</sup>I]-OVTA is 2200 Ci/mmol (81 TBq/mmol), 1723 μCi/μg (63.7 MBq/μg). Preparative HPLC is used to separate unlabeled ornithine vasotocin analog from [<sup>125</sup>I]-OVTA. Upon decay, [<sup>125</sup>I]-OVTA undergoes decay catastrophe and the specific activity remains constant with time. However, it is not known what molecular or peptide fragments are generated from the decay event or what functional activity these fragments may have in different assays. References on <sup>125</sup>I decay and decay catastrophe of <sup>125</sup>I labeled compounds are available.<sup>1-5</sup>

**RADIOCHEMICAL PURITY:** Initially greater than 95% radiochemically pure as determined by HPLC.

**PREPARATIVE PROCEDURE:** OVTA is radioiodinated with no carrier added <sup>125</sup>I using a modification of the Hunter and Greenwood method<sup>6</sup> and purified by reversed phase HPLC.

**AVAILABILITY:** [<sup>125</sup>I]-OVTA is routinely available from stock and is prepared fresh and packaged for shipment on the second Mondays of January, March, May, July, September and November. Please inquire for larger package sizes.  
NEX254-R-REV01

**APPLICATIONS:** [<sup>125</sup>I]-OVTA is a selective oxytocin antagonist which exhibits very low cross reactivity with arginine vasopressin receptors<sup>7</sup>. This, together with its high specific activity, makes it the radioligand of choice for the study of oxytocin receptors.

**HAZARD WARNING:** This product contains a chemical (s) known to the state of California to cause cancer. This product also contains a component which is harmful by contact, ingestion and inhalation. It is irritating to the eyes, skin and respiratory tract and is toxic.

**RADIATION UNSHIELDED:** 280mR/hr/mCi at vial surface.

### REFERENCES:

1. Doyle, V.M., Buhler, F.R., Burgisser, E., *Eur. J. Pharm.* 99 353 (1984).
2. Schmidt, J., *J. Biol. Chem.* 259 1660 (1984).
3. Loring, R.H., Jones, S.W., Matthews-Bellinger, J., Salpeter, M.M., *J. Biol. Chem.* 257 1418 (1982).
4. Berridge, M.S., Jiang, V.W., Welch, M.J., *Rad. Res.* 82 467 (1980).
5. Charlton, D.E., *Rad. Res.* 107 163 (1986).
6. Hunter, W.M. and Greenwood, F.C., *Nature* 194 495 (1962).
7. Elands, J., Barberis, C., Jard, S., Tribollet, E., Jean-Jacques, D., Bankowski, K., Manning, M. and Sawyer, W.H., *Eur. J. Pharmacol.* 147 197-207 (1988).

## IODINE-125 DECAY CHART HALF LIFE=60 days

Radiations: Gamma 35.5 keV (7%), X-ray K alpha 27 KeV (112%), K beta 31 keV (24%)

DAYS	0	2	4	6	8	10	12	14	16	18
0	1.000	.977	.955	.933	.912	.891	.871	.851	.831	.812
20	.794	.776	.758	.741	.724	.707	.691	.675	.660	.645
40	.630	.616	.602	.588	.574	.561	.548	.536	.524	.512
60	.500	.489	.477	.467	.456	.445	.435	.425	.416	.406
80	.397	.388	.379	.370	.362	.354	.345	.338	.330	.322
100	.315	.308	.301	.294	.287	.281	.274	.268	.262	.256
120	.250	.244	.239	.233	.228	.223	.218	.213	.208	.203

To obtain the correct radioactive concentration or amount for a date before the calibration date: divide by the decay factor corresponding to the number of days before the calibration date. To obtain the correct radioactive concentration or amount for a date after the calibration date: multiply by the decay factor corresponding to the number of days after the calibration date.

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.

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

(800) 762-4000  
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.