

# [<sup>125</sup>I]Tyr<sup>A14</sup> Insulin (porcine) [<sup>125</sup>I]-Receptor Grade Insulin

Product Number: NEX196

## LOT SPECIFIC INFORMATION:

CALCULATED AS OF: 26-Aug-2024

LOT NUMBER: CZ92740

SPECIFIC ACTIVITY: 81.4 TBq/mmol  
2200.0 Ci/mmol  
13.7 MBq/μg  
371 μCi/μg

CONCENTRATION: 3.0 MBq/ml  
80.8 μCi/ml

RADIOCHEMICAL PURITY: ≥ 95%

## Package Size Information

Package Size as of 27-Sep-2024	Volume
370 kBq 10 μCi	0.20 ml
1.85 MBq 50 μCi	1.00 ml

MOLECULAR WEIGHT: ~5,927

**PACKAGING:** [<sup>125</sup>I]Tyr<sup>A14</sup> Insulin is in a solution containing 0.05M sodium phosphate, 1M glycine, 0.2M NaCl, 0.25% BSA, 500 KIU/ml Trasylol® at pH 7.2. It is shipped on dry ice.

**SPECIAL INFORMATION:** This compound is light sensitive. Exposure to light may hasten decomposition. [<sup>125</sup>I]Tyr<sup>A14</sup> Insulin is supplied in a red NENSURE™ vial which contains a U.V. inhibitor.

**STABILITY AND STORAGE:** [<sup>125</sup>I]Tyr<sup>A14</sup> Insulin should be stored at -20°C or lower in the dark. Under these conditions the product is stable and usable for at least four weeks after fresh lot date.

**SPECIFIC ACTIVITY:** The initial specific activity of [<sup>125</sup>I]Tyr<sup>A14</sup> Insulin is 2200 Ci/mmol (81 TBq/mmol), 371 μCi/μg (13.7 MBq/μg). Preparative HPLC is used to separate unlabeled insulin from [<sup>125</sup>I]-Tyr<sup>A14</sup>Insulin. Upon decay, [<sup>125</sup>I]Tyr<sup>A14</sup>Insulin 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.

**POSITION OF LABELING:** Limited protease digestion of the product followed by HPLC identification of the fragments<sup>6</sup> indicates that it is the Tyr<sup>A14</sup> which is iodinated.

**PREPARATIVE PROCEDURE:** Porcine insulin is radioiodinated with no carrier added <sup>125</sup>I using a modification of the Hunter and Greenwood method<sup>7</sup>, and is purified by reversed phase HPLC. Only the Tyr<sup>A14</sup> iodinated isomer is collected as product.

**APPLICATIONS:** [<sup>125</sup>I]Tyr<sup>A14</sup> Insulin has been proven to be active in a receptor binding assay using cultured human fibroblasts as described.<sup>8</sup> It may also be used in radioimmunoassay.

**AVAILABILITY:** [<sup>125</sup>I]Tyr<sup>A14</sup> Insulin is routinely available from stock and prepared fresh and packaged for shipment on the fourth Monday of each month. Please inquire for larger package sizes

**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 or inhalation. It is irritating to the eyes, skin and respiratory tract. It is toxic and flammable. Target organ is the central nervous system.

**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., *Radiation Research* 82 467 (1980).
5. Charlton, D.E., *Radiation Research*, 107 163 (1986).
6. Frank, B., Beckage, M.J. and Willey, K.J., *Chrom.* 266, 239-248 (1983).
7. Hunter, W.M. and Greenwood, F.C., *Nature* 194, 495 (1962).
8. Podskalny, J. and Kahn, R., *J. Clin. Endocrinol. and Metab.* 54, 261-268 (1982)

**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.