

# [<sup>125</sup>I]-INSULIN-LIKE GROWTH FACTOR-1 (HUMAN, RECOMBINANT)

## [<sup>125</sup>I]-IGF-1

Product Number: NEX241

### LOT SPECIFIC INFORMATION

**CALCULATED AS OF:** 19-Feb-2024

**LOT NUMBER:** FO32240

**SPECIFIC ACTIVITY:** 69.0 TBq/mmol  
1864 Ci/mmol  
8.9 MBq/μg  
240 μCi/μg

**UNBOUND IODIDE:** < 5%

**MOLECULAR WEIGHT** 7773

### Package Size Information

Package Size as of 22-Mar-2024
185 kBq 5 μCi
925 kBq 25 μCi

**PACKAGING:** [<sup>125</sup>I]-IGF-1 is lyophilized from a solution containing 0.05M sodium phosphate, 1M glycine, 0.2M NaCl, and 0.25% BSA at pH 5.2. It is shipped ambient.

**STABILITY AND STORAGE:** The lyophilized [<sup>125</sup>I]-IGF-1 should be stored at 4°C or lower. Following reconstitution with distilled water to a concentration of approximately 25 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 six weeks after fresh lot date.

**SPECIFIC ACTIVITY:** 210-360 μCi/μg (7.8-13.3 MBq/μg), 1630-2800 Ci/mmol (60-103 TBq/mmol) on fresh lot date as determined from <sup>125</sup>I incorporation into insulin-like growth factor-1. Equimolar protein and iodine implies 285 μCi/μg (10.5 MBq/μg). Specific activity decays with time.

**RADIOCHEMICAL PURITY:** Initially less than 5% unbound iodide as determined by thin layer chromatography

**PREPARATIVE PROCEDURE:** Insulin-like growth factor-1 is radioiodinated with no carrier added <sup>125</sup>I by a lactoperoxidase procedure<sup>1</sup> and purified by reversed phase HPLC. This method predominantly labels tyrosine residues.

**AVAILABILITY:** [<sup>125</sup>I]-IGF-1 is routinely available from stock and is prepared fresh and packaged for shipment on the third Monday of each month. Please inquire for larger package sizes.

**APPLICATIONS:** Insulin-like growth factor-1 (an anabolic peptide hormone) stimulates growth, metabolism, and repair of damaged tissues. [<sup>125</sup>I]-IGF-1 finds use in receptor binding, radioimmunoassay, and autoradiography.

**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 ingestion or inhalation. It is irritating to the

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.

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

**REFERENCES:**

1. Morrison, M., Bayse, G. S., Webster, R. G., *Immunochemistry* 8, 289-297 (1971).

**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%)

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