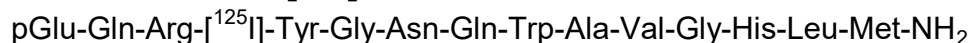


# [<sup>125</sup>I]-Tyr<sup>4</sup>-BOMBESIN

Product Number: NEX258

## [<sup>125</sup>I]-Bombesin



### LOT SPECIFIC INFORMATION

**CALCULATED AS OF:** 25-Nov-2024

**LOT NUMBER:** EUB2740

**SPECIFIC ACTIVITY:**

81.4	TBq/mmol
2200	Ci/mmol
45.4	MBq/μg
1226	μCi/μg

**CONCENTRATION:**

3.20	MBq/ml
86.4	μCi/ml

**RADIOCHEMICAL PURITY:** ≥ 95%

**MOLECULAR WEIGHT:** ~1,794

**PACKAGING:** [<sup>125</sup>I]-Bombesin is in a solution containing 0.05M sodium phosphate, 0.25% BSA, and a stabilizer (pH 3.4): n-propanol, 1:1. It is shipped on dry ice.

**STABILITY AND STORAGE:** [<sup>125</sup>I]-Bombesin should be stored 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:** The initial specific activity of [<sup>125</sup>I]-Bombesin is 2200 Ci/mmol, (81 TBq/mmol), 1230 μCi/μg (46 MBq/μg). Preparative HPLC is used to separate unlabeled bombesin from [<sup>125</sup>I]-Bombesin. Upon decay, [<sup>125</sup>I]-Bombesin undergoes decay catastrophe and the specific activity remains constant with time. However, it is not known what molecular 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:** [Tyr<sup>4</sup>]-bombesin is radioiodinated using no carrier added <sup>125</sup>I by a lactoperoxidase procedure and purified by HPLC. Amino acid analysis indicates that this product is monoiodinated exclusively on Tyr<sup>4</sup>.

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

**APPLICATIONS:** [<sup>125</sup>I]-Bombesin will be useful as a tracer in RIA's for bombesin like peptides, and as a high specific activity ligand in receptor studies.

**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 flammable and explosive may be fatal if swallowed or inhaled

### Package Size Information

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

(4000 ppm dangerous to life or health). This component causes central nervous depression and may be poisonous upon decomposition.

**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 1160 (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).

**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	0.977	0.955	0.933	0.912	0.891	0.871	0.851	0.831	0.812
20	0.794	0.776	0.758	0.741	0.724	0.707	0.691	0.675	0.66	0.645
40	0.63	0.616	0.602	0.588	0.574	0.561	0.548	0.536	0.524	0.512
60	0.5	0.489	0.477	0.467	0.456	0.445	0.435	0.425	0.416	0.406
80	0.397	0.388	0.379	0.37	0.362	0.354	0.345	0.338	0.33	0.322
100	0.315	0.308	0.301	0.294	0.287	0.281	0.274	0.268	0.262	0.256
120	0.25	0.244	0.239	0.233	0.228	0.223	0.218	0.213	0.208	0.203

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.