

[¹²⁵I]-Goat Anti-Rabbit IgG

Product Number: NEX155

LOT SPECIFIC INFORMATION:

CALCULATED AS OF: 5-Feb-2024

LOT NUMBER: BZ30840

SPECIFIC ACTIVITY: 51.9 TBq/mmol
 1402 Ci/mmol
 0.3 MBq/μg
 9.3 μCi/μg

CONCENTRATION: 5.00 MBq/ml
 135.3 μCi/ml

UNBOUND IODIDE: <5% unbound iodine

MOLECULAR WEIGHT: ~150,000

PACKAGING: [¹²⁵I]-Goat anti-rabbit IgG is in a solution containing 0.05M sodium phosphate, 0.15M NaCl, 0.1% BSA, and 0.2% Tween 80® at pH 7.4. It is shipped on dry ice.

STABILITY AND STORAGE: [¹²⁵I]-Goat anti-rabbit IgG should be stored at -20°C or lower. It should be aliquoted in appropriate volumes to avoid repeated freeze-thaw cycles. Under these conditions, the product is stable and usable for at least four weeks after fresh lot date.

SPECIFIC ACTIVITY: 2-10 μCi/ μg (74-370 kBq/ μg) on fresh lot date as determined from ¹²⁵I incorporation into goat anti-rabbit IgG. Specific activity decays with time.

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

PREPARATIVE PROCEDURE: Affinity purified goat anti-rabbit IgG is radioiodinated with no carrier added ¹²⁵I using a modification of the Hunter and Greenwood method¹ and is purified by gel filtration chromatography. This method predominantly labels tyrosine residues.

AVAILABILITY: [¹²⁵I]-Goat anti-rabbit IgG is routinely available from stock and is prepared fresh and packaged for shipment on the first 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.

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

Package Size Information

Package Size as of 8-Mar-2024	Volume
3.70 MBq 100 μCi	1.25 ml
9.25 MBq 250 μCi	3.125 ml

REFERENCES:

1. Hunter, W.M. and Greenwood, F.C., *Nature* 194 495 (1962).

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

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
Copyright ©2023, Revvity, Inc. All rights reserved.