

[¹²⁵I]-Goat Anti-Rabbit IgG

Product Number: NEX155

LOT SPECIFIC INFORMATION:

CALCULATED AS OF: 1-Apr-2024

LOT NUMBER: BZ51040

SPECIFIC ACTIVITY: 0.0 TBq/mmol
0 Ci/mmol
0.0 MBq/μg
0.0 μCi/μg

CONCENTRATION: 5.45 MBq/ml
147.2 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% 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 of 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 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	
10-May-2024	
3.70 mBq	
100 μCi	
9.25 MBq	
250 μCi	