

[¹²⁵I]-[NLe⁴, D-Phe⁷]-α-MELANOCYTE STIMULATING HORMONE

Product Number: NEX352

[¹²⁵I]-[NLe⁴, D-Phe⁷]-α-MSH**LOT SPECIFIC INFORMATION**

CALCULATED AS OF: 19-Feb-2024

LOT NUMBER: IM32240

SPECIFIC ACTIVITY: 81.4 TBq/mmol
 2200 Ci/mmol
 46 MBq/μg
 1243 μCi/μg

CONCENTRATION: 3.0 MBq/ml
 82.4 μCi/ml

RADIOCHEMICAL PURITY: ≥ 95%

MOLECULAR WEIGHT: ~1770

Package Size Information

Package Size as of 22-Mar-2024	Volume
370 kBq 10 μCi	0.20 mL
1.85 MBq 50 μCi	1.00 mL

PACKAGING: [¹²⁵I]-[NLe⁴, D-Phe⁷]-α-MSH is in a solution containing 0.08M Tris HCl, 0.08M NaCl, 0.05M β-mercaptoethanol, 0.25% BSA, 50 KIU/ml Trasylo[®], (pH 8.5); ethanol, 5:1. It is shipped on dry ice.

STABILITY AND STORAGE: [¹²⁵I]-[NLe⁴, D-Phe⁷]-α-MSH should be stored at -20°C. 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 [¹²⁵I]-[NLe⁴, D-Phe⁷]-α-MSH is 2200 Ci/mmol, (81 TBq/mmol), 1243 μCi/μg (46 MBq/μg). Preparative HPLC is used to separate unlabeled [NLe⁴, D-Phe⁷]-α-MSH from [¹²⁵I]-[NLe⁴, D-Phe⁷]-α-MSH. Upon decay, [¹²⁵I]-[NLe⁴, D-Phe⁷]-α-MSH 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 ¹²⁵I decay and decay catastrophe of ¹²⁵I labeled compounds are available.¹⁻⁵

RADIOCHEMICAL PURITY: Initially greater than 95% radiochemically pure as determined by HPLC.

