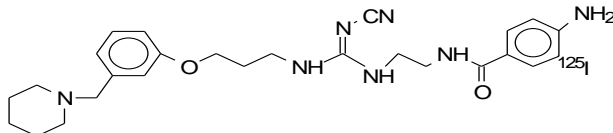


[¹²⁵I] AMINOPOTENTIDINE

Product Number: NEX431

**LOT SPECIFIC INFORMATION**

CALCULATED AS OF: 16-Sep-2024

LOT NUMBER: LN02540

SPECIFIC ACTIVITY: 81.4 TBq/mmol
 2200 Ci/mmol
 135 MBq/μg
 3657 μCi/μg

CONCENTRATION: 6.4 MBq/ml
 172.4 μCi/ml

RADIOCHEMICAL PURITY >95%

MOLECULAR WEIGHT: 601.6

Package Size Information

Package Size as of 25-Oct-2024	Volume
370 kBq 10 μCi	0.10 mL
925 kBq 25 μCi	0.25 mL
92.5 MBq 250 μCi	2.5 mL

PACKAGING: [¹²⁵I]-Aminopotentidine is supplied in an ethanol and water solution and dispensed in TPX vials. It is shipped ambient.

STABILITY AND STORAGE: [¹²⁵I]-Aminopotentidine should be stored at 2-8°C. Under these conditions the product has been shown to stable for at least four weeks after the fresh lot date.

SPECIFIC ACTIVITY: The initial specific activity is 2200 Ci/mmol, (81.4 TBq/mmol), 3657 μCi/μg (135 MBq/μg). Upon decay, the product 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.

PREPARATIVE PROCEDURE: [¹²⁵I]-Aminopotentidine is radioiodinated with no carrier added ¹²⁵I using a modification of the Hunter and Greenwood method⁶ (chloramine-T) and purified by reversed phase HPLC.

AVAILABILITY: [¹²⁵I]-Aminopotentidine 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.
 NEX431-R-REV01

packaged for shipment on the third Monday of each month. Please inquire for larger package sizes.

APPLICATIONS: [¹²⁵I]-Aminopotentidine has been shown to work in receptor binding and autoradiography and is an antagonist for histamine (H₂) receptors.⁷

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 contact, ingestion and inhalation. It is irritating to the eyes, skin and respiratory tract and is toxic.

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).
6. Hunter, W.M. and Greenwood, F.C., *Nature* 194 495 (1962).
7. Ruat, M., Traiffort, E., et al, *Proc. Natl. Acad. Sci. USA, Vol 87, p.1658 March 1990*

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
0	1	0.977	0.955	0.933	0.912	0.891	0.871	0.851
20	0.794	0.776	0.758	0.741	0.724	0.707	0.691	0.675
40	0.63	0.616	0.602	0.588	0.574	0.561	0.548	0.536
60	0.5	0.489	0.477	0.467	0.456	0.445	0.435	0.425
80	0.397	0.388	0.379	0.37	0.362	0.354	0.345	0.338
100	0.315	0.308	0.301	0.294	0.287	0.281	0.274	0.268
120	0.25	0.244	0.239	0.233	0.228	0.223	0.218	0.213

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 constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity" 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 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
Copyright ©2023, Revvity, Inc. All rights reserved.



16
0.831
0.66
0.524
0.416
0.33
0.262
0.208

d does not
/") do not
AIMS ALL
ORAL OR
ON WITH

