

GET TO KNOW OUR ULTRAPURE YTTRIUM-90

90
Y
yttrium

What makes our NEN® Yttrium-90 (⁹⁰Y)* radionuclide unique? It's extracted from Strontium-90 using a proprietary process, developed by PerkinElmer, in compliance with GMP**.

In labs like yours, ⁹⁰Y radionuclide is combined with monoclonal antibodies and peptides to create potential "smart drugs" with specific targets. A promising therapeutic application, this process is being studied for treating cancers such as ovarian, lung, breast, colon, prostate, brain, and non-Hodgkin's lymphoma.

Every lot of ⁹⁰Y is quality tested to ensure consistency. Documentation of test results at release, including ICP analysis for metals contamination, is provided with each customer shipment. Our ⁹⁰Y is fresh every week, and each vial produced is custom dispensed to meet the requirements of each unique application.

Expect quality products and a global distribution network to ensure timely delivery.

Features and Benefits of ⁹⁰Y

- High purity – low levels of contaminants
- High radiochemical concentration – can be diluted to meet specific lab requirements
- Excellent labeling efficiency – high incorporation
- Available weekly – reliable production schedule
- Global distribution capability – assures worldwide delivery soon after production
- Custom dispensed – flexibility to meet your needs
- GMP** Compliant Manufacturing – assurance of consistent product quality
- Unique NEN® product packaging - safe, secure, and maximizes recovery of material while ensuring minimal personal exposure

*Yttrium-90 (⁹⁰Y), is for investigational use only and not for use in humans. PerkinElmer makes no clinical use representations.

**Manufactured under an ISO 9001:2015 Quality Management System registration and consistent with ICH Q7 Good Manufacturing Practice Guidance for Active Pharmaceutical Ingredients, August 2001.

Yttrium-90 (⁹⁰Y) Specifications

Product Name	Yttrium-90* (Y-90) GMP** Compliant
Catalog Number	NEZ306, NEZ306A, NEZ306B, NEZ306C, NEZ306D
Fresh Lot	Each week; calibrated for Monday (NEZ306), Tuesday (NEZ306A), Wednesday (NEZ306B), Thursday (NEZ306C), or Friday (NEZ306D).
Half-life	64 hours
Decay Mode	Beta, 2.28MeV Maximum
Chemical Form	As Yttrium chloride in ~0.05 N HCl
Specific Activity	~500 Ci/mg (carrier free)
Radiochemical Concentration	~5,000 mCi/mL at production
Radionuclide Purity	< 1.8E-7Ci/Ci at production
Chemical Purity	µg of analyte/Ci of Y-90:
	< 100 µg/Ci Ca < 20 µg/Ci Fe
	< 20 µg/Ci Zn < 20 µg/Ci Cr
	< 20 µg/Ci Al < 20 µg/Ci Pb
	< 20 µg/Ci Cu < 100 µg/Ci P
	< 20 µg/Ci Zn+Fe+Cu+Pb

*Yttrium-90 (⁹⁰Y), is for investigational use only and not for use in humans. PerkinElmer makes no clinical use representations.

**Manufactured under an ISO 9001:2015 Quality Management System registration and consistent with ICH Q7 Good Manufacturing Practice Guidance for Active Pharmaceutical Ingredients, August 2001.

**For more information on ⁹⁰Y or to access product decay tables please visit us at
www.perkinelmer.com/radiotherapeutics**

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

014610_01

PKI