

HUMAN HEALTH

ENVIRONMENTAL HEALTH




HOW WELL
DO YOU KNOW
RAD
DETECTION?

Radiochemical Custom Synthesis

Over fifty years of experience in labeling technologies allows us to label virtually any biomolecule. Our expertise in assay development and custom synthesis help streamline your projects and save valuable time. Rely on PerkinElmer to have what you need, when you need it.

	ISO 9001:2000	Technical Data Sheets	Total Activity Measurement	Specific Activity Measurement	Radiochemical Purity or Synthesis	Chiral Separation	GMP (upon Request)
¹⁴ C, ³ H, Custom Synthesis	■	■	■	■	■	■	■
³³ P, ³⁵ S, Custom Synthesis	■	■	■	■	■	■	
¹²⁵ I, ¹³¹ I, Radioiodination	■	■	■	■	■	■	
Tritium Labeling Plus	■	■	■	■	■	■	
Tritium Labeling Services							
Catalytic Reduction with Tritium Gas	■	■	■				
Reduction with Sodium Borohydride, ^{[3} H]	■	■	■				
Catalytic Exchange with Tritium Gas	■	■	■				
Tritium Gas Exposure	■	■	■				
¹⁴ C-Labeled Precursors	■	■	■	■	■	■	

NEN® GMP* Custom Synthesis

Customers developing pharmaceutical products are faced with the challenge to provide clinical study materials which meet escalating GMP requirements. As proof of our commitment to you, PerkinElmer is able to help you meet this goal by offering GMP custom services.

Why choose GMP Custom Synthesis from PerkinElmer?

- Dedicated GMP Custom Synthesis Lab
- Complete documentation
- Project team environment
- Experience and expertise
- Manufacturing excellence
- Guaranteed confidentiality

* Compliance with Q7A Good Manufacturing Practice Guidance for Active Pharmaceutical Ingredients (issued 8/2001), Section XIX, APIs for use in clinical trials.

¹⁴C, ³H, ³⁵S and ³³P Custom Synthesis Services

Our custom synthesis scientists are experts in dealing with the technically challenging synthesis of radiolabeled biochemicals, such as peptides, steroids, lipids, sulfonamides and others. They may collaborate with chemists from our Analytical Laboratory to ensure the successful purification and analysis of final products.

Tritium Labeling Service

Our Tritium Labeling Service is a cost-effective way to prepare a ³H-labeled radiochemical. The process includes:

- Reaction of the precursor you submit, according to the labeling method you select.
- Removal of the catalyst and labile tritium from the resulting crude product.
- Assay of the resulting crude product for total radioactivity.

To request a custom quote or for further information, please visit our Web site at www.perkinelmer.com/customsynthesis.

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 ©2009-2010, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

008241D_01

Printed in USA

¹²⁵I and ¹³¹I Radioiodination Services

PerkinElmer will label your compound with ¹²⁵I or ¹³¹I using the most appropriate radioiodination method and reagent quantities to meet your specifications. We can perform large scale iodinations (>10 mCi).

Available methods:

- Bolton-Hunter Reagent
- Chloramine-T
- IODOGEN® Reagent
- Lactoperoxidase
- Exchange Labeling with Sodium Iodide
- Custom Labeling Methods

³³P and ³⁵S Nucleotide and Oligonucleotide Synthesis

PerkinElmer will synthesize and package radiolabeled mono-, di- and triphosphate nucleotides, oligonucleotides and other related biochemicals to your specifications on a make-to-order basis. We have extensive experience in labeling nucleotides in different positions (e.g., beta position) and synthesizing labeled oligos for hybridization.

¹⁴C-Labeled Precursor Materials

To assist in your in-house synthesis of labeled compounds, PerkinElmer offers a number of ¹⁴C-labeled precursors that can be purchased in large quantities.

These precursors are routinely produced and are readily available with the same high purity specifications as comparable catalog compounds. We will provide custom packaging in special glassware to meet your unique protocol requirements.

