



HOW WELL DO YOU KNOW  
RAD DETECTION?

## Radiochemical Custom Synthesis



Over sixty 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 9001:2015	Technical Data Sheets	Total Activity Measurement	Specific Activity Measurement	Radiochemical Purity or Synthesis	Chiral Separation	GMP (upon Request)
<sup>14</sup> C, <sup>3</sup> H, Custom Synthesis	■	■	■	■	■	■	■
<sup>32</sup> P, <sup>35</sup> S, Custom Synthesis	■	■	■	■	■	■	
<sup>125</sup> I, <sup>131</sup> I, Radioiodination							
Bolton-Hunter Reagent	■	■	■	■	■	■	
Chloramine-T	■	■	■	■	■	■	
IODOGEN® Reagent	■	■	■	■	■	■	
Lactoperoxidase	■	■	■	■	■	■	
Custom Labeling Methods	■	■	■	■	■	■	
<b>Tritium Labeling Services</b>							
Catalytic Reduction with Tritium Gas	■	■	■				
Reduction with Sodium Borohydride, [ <sup>3</sup> H]	■	■	■				
Catalytic Exchange with Tritium Gas	■	■	■				
Tritium Gas Exposure	■	■	■				
<sup>14</sup> C-Labeled Precursors	■	■	■	■	■	■	

## 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, we are 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 ICH Q7 Good Manufacturing Practice Guidance for Active Pharmaceutical Ingredients (Section 19), August 2001, APIs for use in clinical trials.

## <sup>14</sup>C, <sup>3</sup>H, and <sup>35</sup>S 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 <sup>3</sup>H-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

## <sup>125</sup>I and <sup>131</sup>I Radioiodination Services

PerkinElmer will label your compound with <sup>125</sup>I or <sup>131</sup>I 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
- Custom Labeling Methods

## <sup>32</sup>P and <sup>35</sup>S Nucleotide Synthesis

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

## <sup>14</sup>C-Labeled Precursor Materials

To assist in your in-house synthesis of labeled compounds, PerkinElmer offers a number of <sup>14</sup>C-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.

To request a custom quote or for further information, please visit our Web site at <http://www.perkinelmer.com/category/custom-radiosynthesis-radiolabeling-services>

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



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2009-2018, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

008241E\_01

PKI