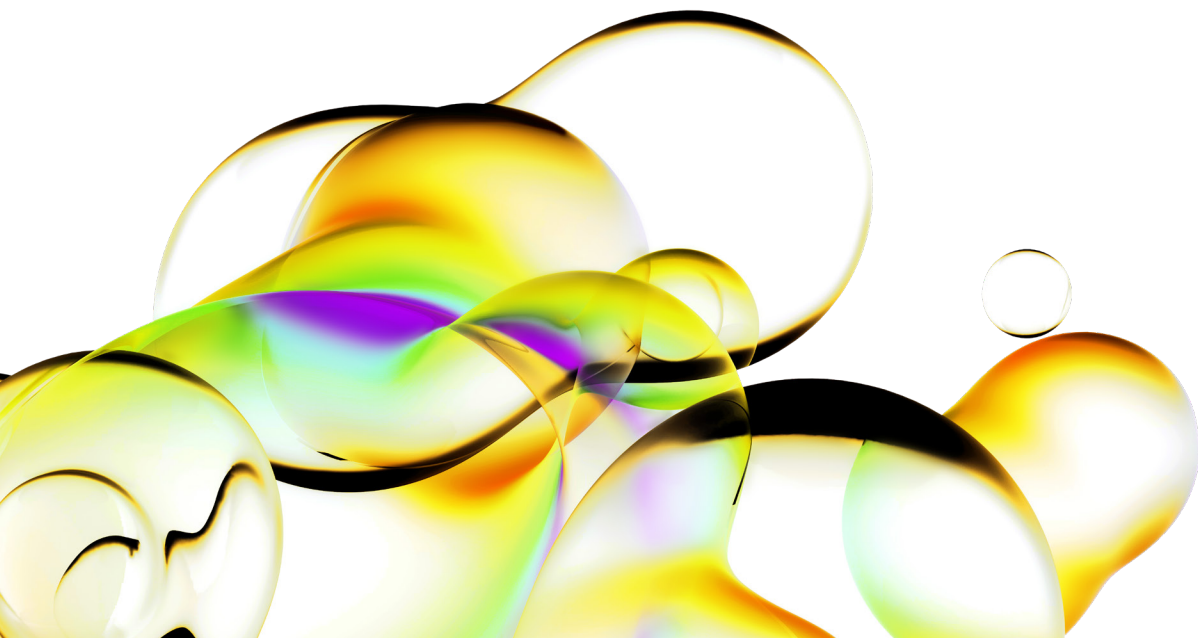




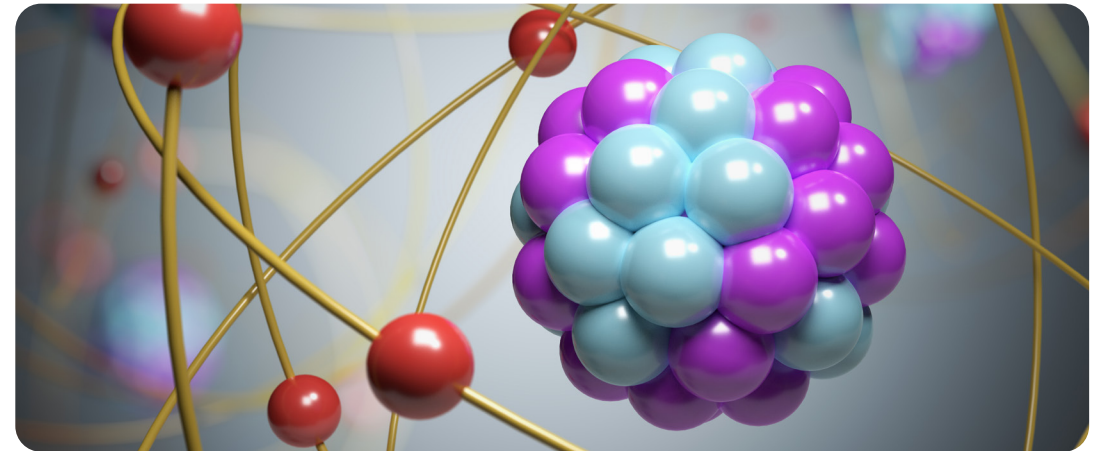
You discover,
we customize.



Radiochemical custom synthesis

Over sixty years of experience in labeling technologies allows us to label many biomolecules. Our expertise in custom radiosynthesis and labeling help streamline your projects and save valuable time. Rely on Revvity to have what you need, when you need it.

To request a custom quote or for further information, please visit our website at <http://www.revvity.com>



	ISO 9001:2015	Technical data sheets	Total activity measurement	Specific activity measurement	Radiochemical purity	Chiral separation	GMP (Upon Request)
¹⁴C Custom synthesis	▪	▪	▪	▪	▪	▪	▪
¹²⁵I Radioiodination							
Bolton-hunter reagent	▪	▪	▪	▪	▪	▪	
Chloramine-T	▪	▪	▪	▪	▪	▪	
IODOGEN® Reagent	▪	▪	▪	▪	▪	▪	
Lactoperoxidase	▪	▪	▪	▪	▪	▪	
Custom labeling methods	▪	▪	▪	▪	▪	▪	
Tritium labeling services							
Catalytic reduction with tritium gas	▪	▪	▪				
Reduction with sodium Borohydride, ^[3H]	▪	▪	▪				
Catalytic exchange with tritium gas	▪	▪	▪				
Tritium gas exposure	▪	▪	▪				



¹⁴C 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.



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.

- 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), September 2016, APIs for use in clinical trials.



¹²⁵I radioiodination services

Revvity will label your compound with ¹²⁵I 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



³⁵S services

Revvity will synthesize and package ³⁵S nucleotides labeled on the alpha or gamma phosphate and provide other related reagents to your specifications on a made-to-order basis.



³H custom labeling services

Our tritium labeling service is a cost-effective way to prepare a ³H-labeled radiochemical. The process includes:

- Reaction of the material 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

