

Revvity: Empowering research
in common and rare diseases.



revvity



Reliable and early diagnosis is the cornerstone of disease management. With cutting-edge technologies and a commitment to innovation, we provide comprehensive solutions to support non-profit organizations and biopharmaceuticals dedicated to unraveling the mysteries of common and rare diseases.

Assay development

- Range of assay modalities available in-house.
- Technical expertise across life sciences and diagnostics.
- Manufacturing of custom reference materials to support assay QC.

Preclinical services

- CRISPR based screening.
- Viral Vector design & manufacturing.

- Global reagents and instruments footprint.
- Accredited sites.
- Secure raw materials supply.

Clinical services

- Global lab network.
- Population screening.
- Preclinical testing and clinical trials.
- Sponsored testing programs.
- Clinical expertise.

Regulatory expertise

- Experienced IVD, CE IVDR & CDx filing support for multiple regulatory agencies.
- Global regulatory expertise with footprint in more than 40 accredited sites.

Manufacturing

- Global commercial footprint for deep reach.
- Diagnostics path to market.
- Established global sales team.

Commercial channel

Explore our solutions

Innovative, scalable, end-to-end solutions that are uniquely positioned to meet your needs.

Goals

Revvity solutions

STEP 1 Discovery & development

1. Target identification
2. Assay development

- Assay development for biomarker identification.
- Cell line engineering services for generation of cell models.
- CRISPR, RNAi, base editing, and gene expression platforms for determining gene function and target ID.
- Analysis solutions to reliably detect, analyze and characterize cells, protein or nucleic acids.

STEP 2 Preclinical research

1. Authentication of identified target
2. Lead development and optimization
3. *In silico* assay screening

- Combine functional genomic and cell panel screening services to identify and confirm targets.
- Test and optimize different formulations or oligos using high-throughput platforms.
- Functional genomic screening services to identify drug mechanisms.
- Immune response and toxicology insights with immune cell assay services.

STEP 3 Clinical development

1. Sample collection, storage & logistics
2. Individual stratification
3. Testing for clinical trials

- Providing end-to-end support required for sample collection to clinical testing solutions using next generation sequencing technologies.
- State-of-the-art bioinformatics pipelines designed to deliver the most reliable data processing and interpretation.
- Fluorospot and Flow cytometry to measure the efficacy of your vaccine or immunotherapy.
- Individual stratification for clinical trials through our global laboratory network.

STEP 4 FDA review

1. Regulatory compliance
2. FDA approval application

- Guidance for regulatory and compliance approval.
- Through a dynamic process, we consult with regulatory bodies to facilitate the introduction of innovative products to market.

STEP 5 Companion Diagnostics

1. Manufacturing and commercialization of CDx kits

- Ability to develop and commercialize the kits for companion diagnostic assays for targeted therapies.

STEP 6 Study expansion

1. Program management
2. Global laboratory network

- Dedicate program management support through out the lifespan of the study.
- Global commercial footprint for deeper reach.
- Support to scale up the study through the global lab network.

Our commitment to common and rare diseases research

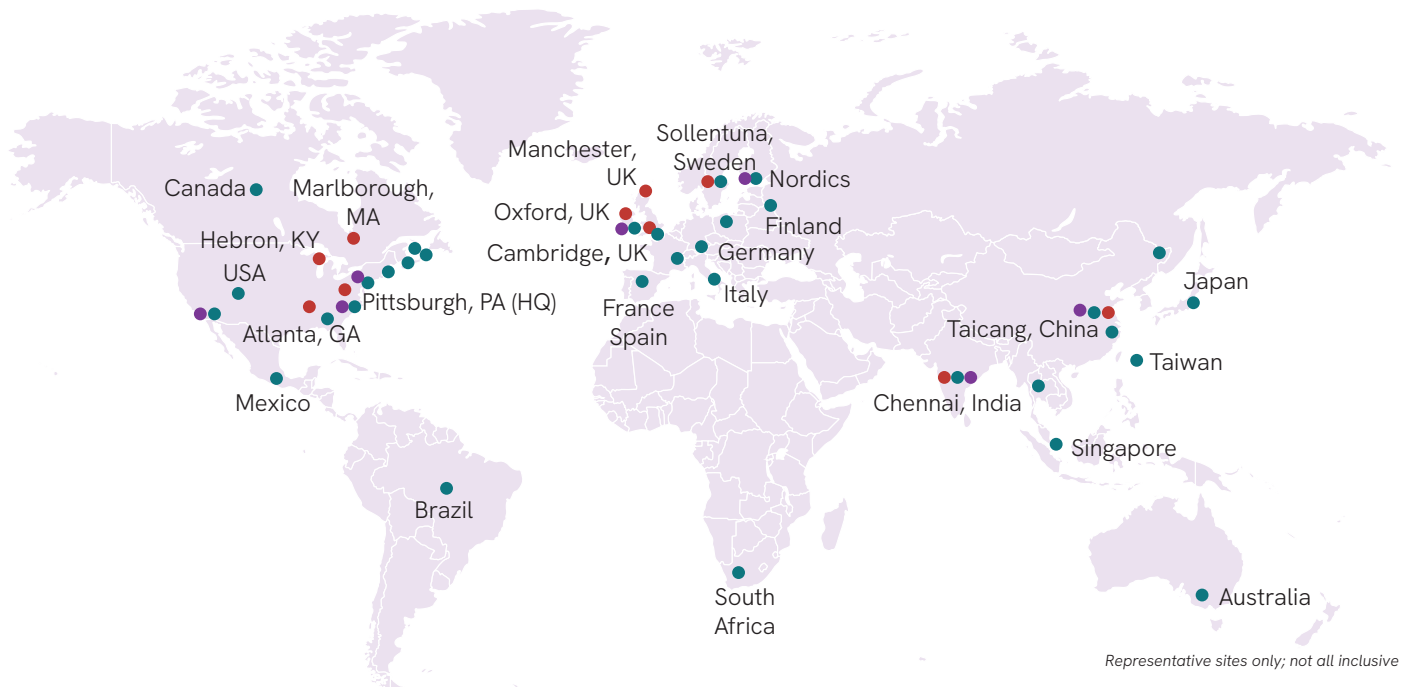
Manufacturing

- Global footprint of **40+ sites**, with **multiple dedicated manufacturing facilities**.
- **Research use only, CE-IVD and FDA cleared** reagents, instruments and consumables, including sample preparation, molecular and NGS assays, and lab automation.
- Deep **in-territory-for-territory** expertise.

Quality & Regulatory

- Global footprint of **40+ accredited sites** offering experience and deep territory expertise across all regulatory authorities.
- QRC (Quality, Regulatory, and Clinical Affairs) plays an **integral role in the product lifecycle** from concept to commercial launch, plus post market surveillance to drive continuous improvement.
- QRC, as a **cross-functional shared service**, uses dynamic resource allocation, providing the right people and skill sets at the right place and right time.
- Through a dynamic process, we **consult with regulatory bodies** to facilitate the introduction of innovative products to market.

A global network



Laboratory sites



QRC sites Manufacturing sites



Disclaimer: This testing service has not been cleared or approved by the U.S. Food and Drug Administration. Testing services may not be licensed in accordance with the laws in all countries. Laboratory licensing and accreditation, as well as the availability of specific test offerings, is dependent upon laboratory location. The content of this pamphlet is provided for informational purposes only, not as medical advice. It is not intended to substitute the consultation, diagnosis, and/or treatment provided by a qualified licensed physician or other medical professional.

