

Start up, Speed up, and Scale up Your Discovery

Enzo is an integrated life sciences company, diagnostic manufacturer, and clinical service provider delivering faster, better, and cost-effective clinical trial management services. Our services rely on 45 years of innovation, technology development, and manufacturing know-how to support your discovery, no matter how small or complex they may be.

Put our expertise in labeling and detection, genomic analysis, protein analysis, cell analysis, tissue analysis, and small molecule chemistry to work for you.

Life Sciences Contract Services



Custom Manufacturing Services

- Volumes
- Packaging
- Formulations
- Kit configuration



Assay & Validation Services

- Validation services
- Stability studies
- Western blot services
- ELISA services



GMP Services

- Small Molecules
- Proteins, cytokines, & growth factors
- Antibodies



ISH Custom Probes

- AMPVIEW™ RNA probes design and development
- Tissue & cell optimization
- Target validation



TECHNICAL EXPERTISE

FLEXIBLE

COST-EFFECTIVE

COLLABORATION

QUALITY



Start up, Speed up, and Scale up Your Discovery



Accelerate Your Drug Discovery with a Trusted Partner

Technical Expertise	Flexibility	Cost-Effective	Quality	Collaboration
Broad expertise in genomics, protein analysis, cell analysis, tissue analysis and small molecules chemistry	Experienced with small and simple projects to multi-million-dollar complex projects	Supply chain security from IP to commercialization	Committed to GxP, ISO 9001, and ISO 13485	Our SCIENTISTS ENABLING HEALTHCARE™ are ready to plan and support your goals and deliverables
Bringing to our partners over four decades of innovation and manufacturing know-how	Assisting with small scale to large scale	A unique integrated structure as a life science company, diagnostic manufacturer, and clinical service provider	Providing quality and reliable products to the drug discovery market	Our dedicated project managers will communicate with our partners throughout the process to ensure we meet our shared goals
High quality insight provided by our subject matter experts		Reducing cost and timeframe for our partners by doing more in-house	We provide clinical services in a New York State Department of Health regulated, CAP accredited, and CLIA certified lab	

Learn More at: enzolifesciences.com/life-sciences-contract-services



Global Headquarters
ENZO LIFE SCIENCES, INC.
 Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES (ELS) AG
 Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg
 Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
 Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
 Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
 Phone (UK customers):
 0845 601 1488
 Phone: +44 1392 825900
info-uk@enzolifesciences.com

For local distributors and detailed product information visit us online: enzolifesciences.com
 All product names, logos, and brands are the property of their respective owners.