

Point of Care Device Testing Services

From Benchtop to Bedside, We'll Help You Get There

With the addition of four donor centers, we now offer point of care (POC) device testing services. Let us operationalize your protocol and bring it to life. We will optimize our donor cohorts to get your POC device testing done quickly and efficiently.



Your Time is Valuable, Let Us Do the Work

- US and UK donor centers with on-site collection and testing services
- Access to specimen analysis instrumentation
- Recruitment expertise for clinical indications



Versatility is Key to Your Success

- We offer two ways to get your testing done:
 - Train our staff on your products and let us do the device testing for you
 - Rent our space and have your own operators perform the testing
- Generate a range of data sets for precision, accuracy, specificity and more!
- Diverse donor base with thousands of healthy donors of biofluid samples



Licensed and Certified Facilities

- Qualified support staff operating under:
 - FDA (US) and HTA (UK) registration and license
 - IRB-approval
 - CLIA Certificate of Compliance



BioIVT's medical director and scientific advisor have 40+ years of experience overseeing clinical trials and POC test validations under IRB protocols.

Collection Types

Blood (finger stick or venous draw), oral swab, nasal swab, throat swab, skin swab, saliva, urine, feces and more.

Donor Cohorts

We are able to recruit specific donor cohorts for your individual POC testing needs including cardiovascular diseases, endocrine/metabolic disorders, autoimmune and allergy conditions, infectious disease and more.

BioIVT's POC Testing Services Cover Technology and Non-technology Components

Considerable progress in the POC arena has come through new and expanded technologies, especially in today's digital age. At our donor centers, we can ensure that key components such as assay chemistry and microfluidics function as designed. In addition, we can integrate other components to guarantee a seamless POC ecosystem.

Technology Components				
Assay Chemistry Does your POC isolate, detect and quantify the target molecule as expected?	Microfluidics Does your POC seamlessly integrate and function as expected?	Connected Instrumentation Does your POC interact and function in real world connectivity?	Data Analytics Does your POC transfer information properly and with proper security?	Systems Integration Does your POC function in different POC settings?

Non-Technology	
Clinical Workflow Does your POC fit into diverse settings and have ease of use?	Regulatory Guidance Does your POC ensure privacy and security of protected health-related information?

"We have relied on quality testing by BioIVT to assess our prototypes and validate our POC devices for a number of years."

Mary Donadi
 Clinical Trials Manager, OraSure

