



Understanding your vaccine and immunotherapy efficacy.

Certimmune's global immunology services are brought to you by Revvity, Inc.

Revvity is a pioneer in the development of more reliable, high throughput, standardized tests for measuring the adaptive immune system. Certimmune is built on the foundations of Revvity nearly 20 years' experience in this area.

Our team of experts can support your **T and B cell immunoassay** needs from start to finish of your **clinical trials**, including:

- Assay design, development and validation
- High-volume testing implementation and operations
- Data interpretation and reporting
- Regulatory responses and quality management
- Global partner lab onboarding and monitoring
- Sample cryopreservation and long-term storage



Partner with the experts

Gain valuable efficacy data, faster, for regulatory approval of your vaccine or immunotherapy. Scan QR code to learn more.

