



Unlock immune research possibilities: T-SPOT.Flex.

T-SPOT™.Flex delivers a trusted interferon-gamma (IFN- γ) ELISPOT kit combining WHO-acknowledged methodology¹ with ready-to-use components, providing a customizable immune monitoring platform that reduces hands-on time while enhancing result quality.

Why choose T-SPOT.Flex?

- **Proven foundation:** a WHO-acknowledged technology¹ supported by Revvity's 20+ years of T cell expertise and 500+ peer-reviewed publications
- **Flexible application:** adaptable platform for diverse research applications providing complete freedom to explore any research area with your own preferred targets
- **Reliable results:** ready-to-use reagents save a full day* of prep work while minimizing background noise for enhanced accuracy of results

T-SPOT.Flex enables you to:

- Design customized assays with your own antigens
- Generate reliable immune response data across diverse research areas
- Benefit from established T cell processing expertise with ready to use reagents
- Maintain workflow efficiency with potential automation compatibility and extended sample stability

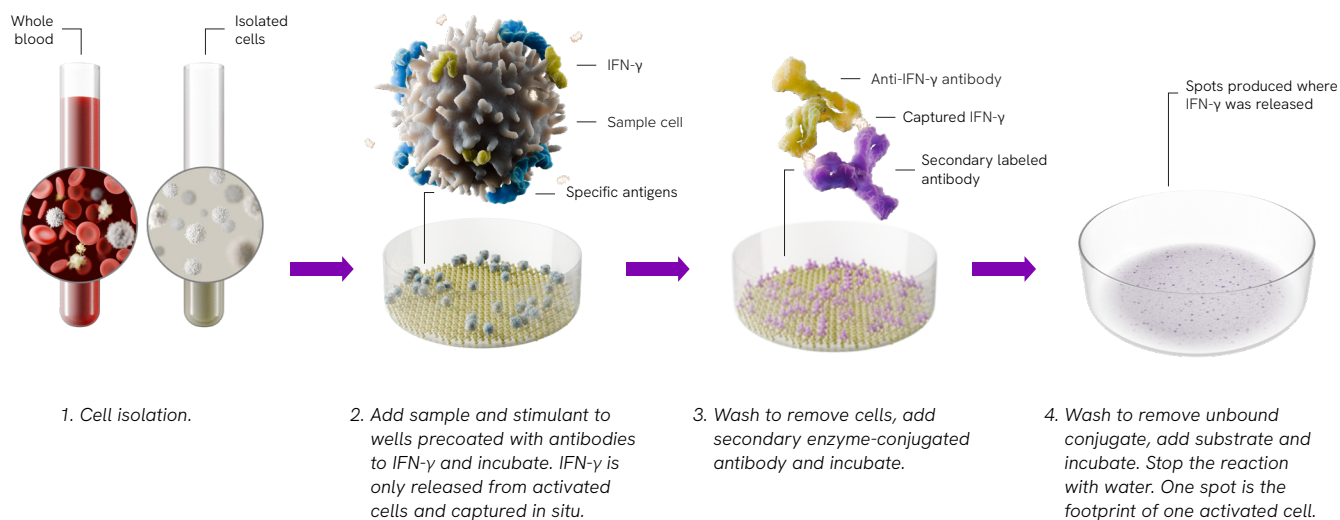


The T-SPOT technology

T cells are crucial components of the adaptive immune system, playing a central role in orchestrating almost all adaptive immune responses. By utilizing the T-SPOT technology (based on the ELISPOT methodology) to

perform an interferon-gamma release assay (IGRA), T-SPOT.Flex can help you accurately identify and analyze T cell responses to specific pathogens, aiding in the detection and monitoring of various immune responses.

T-SPOT technology workflow



* No plate coating or control panel reconstitution necessary

¹⁾ WHO operational handbook on tuberculosis. Module 3: diagnosis. Geneva: World Health Organization; 2025. Licence: CC BY-NC-SA 3.0 IGO.

Please visit our product webpage to access instructions for use and ordering information



Learn more

revvity