



Tag-lite services to evaluate your biosimilar or biobetter samples.

Development of biosimilar or biobetter antibodies requires extensive *in vitro* characterization of the product. This process must include relevant readouts, in particular binding to the representative isoforms of the 3 Fcγ receptors involved in T cell mediated ADCC (FcγRI, FcγRII and FcγRIII) and to FcRn.

Revvity's binding profile services will characterize your samples on all Fcγ and FcRn receptors (CD32A, CD32B, CD64, and CD16) using our cellbased assays.

Benefit from the expertise of Revvity binding service to profile and study your Fcγ receptor

- Direct relationship with a Project Manager for discussion and data review
- Reliable data and fast delivery
- Flexibility: your samples, in any isoform
- Full report: complete dose response curves and data



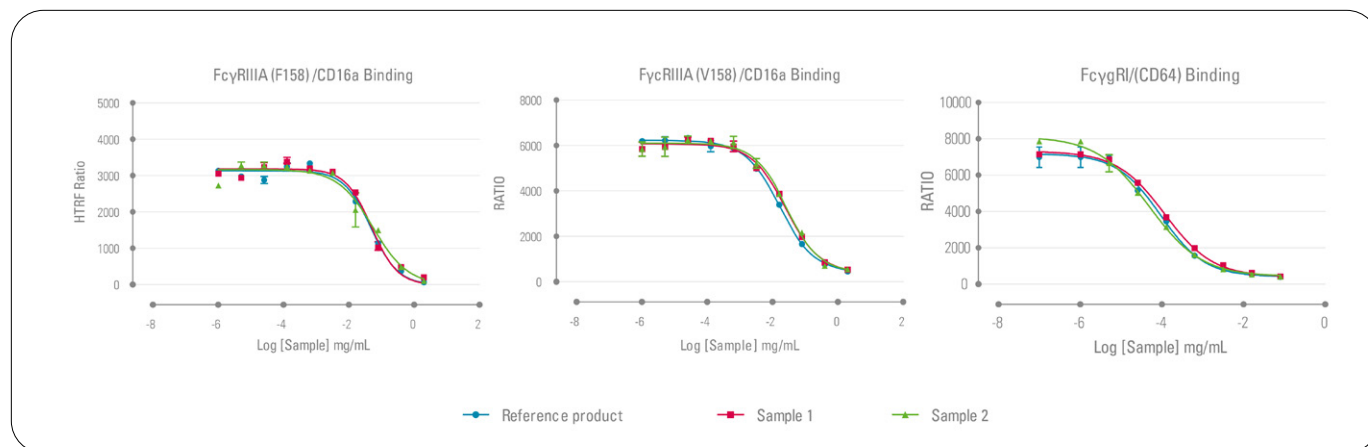
Revvity's team of experienced experts works hand-in-hand with you to characterize your products.

How it works

Profiling

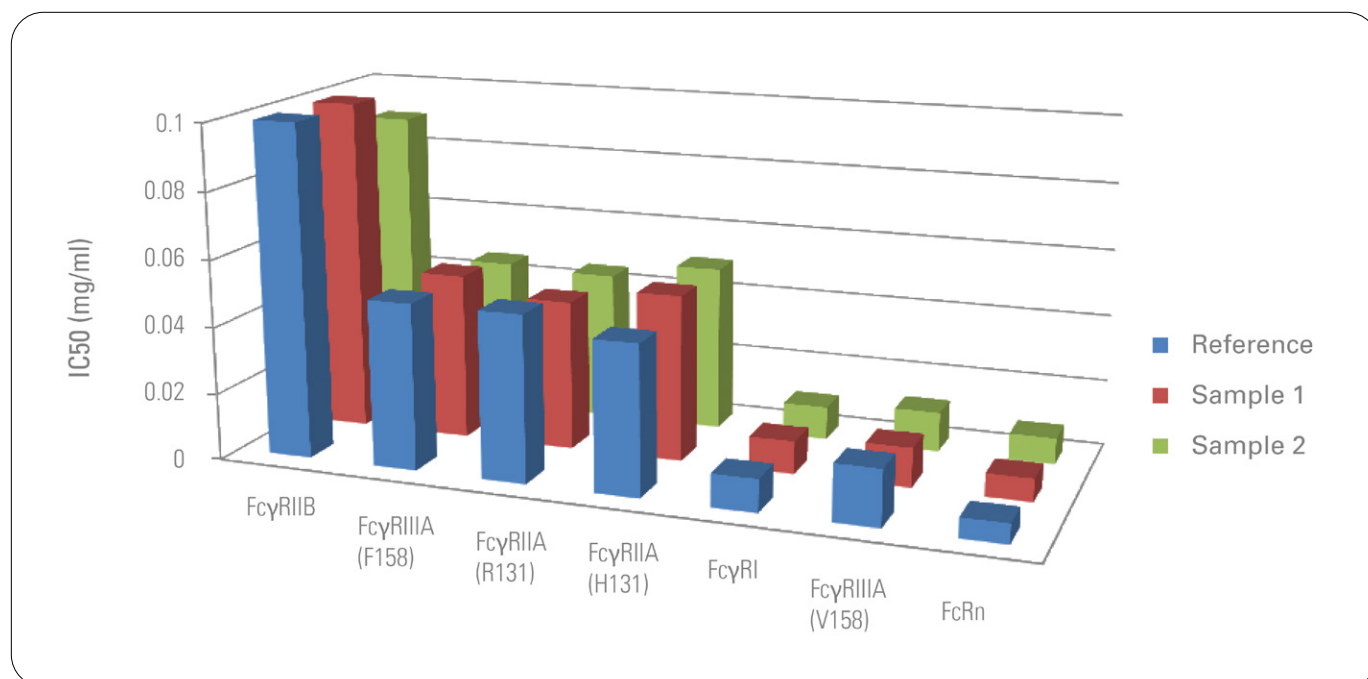
In this case study, 3 samples (2 samples and 1 reference product) were assessed for binding affinity to FcγRI, FcγRIIA, FcγRIIB, FcγRIIA and FcRn using the Revvity cell-based platform.

All samples were tested in dose response (8 concentrations) on each receptor. Binding potency was then determined and compared to establish comparability.



Analysis

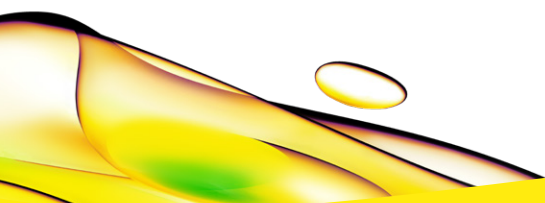
IC50 values were determined for each sample and each receptor tested. Sample 1, 2 and reference show comparable binding profiles for all receptors and all variants tested.



Expert services to satisfy your exact needs

In addition to providing ready-to-use solutions, Revvity's team of experienced experts also works hand-in-hand with clients to create custom solutions. From compound characterization, to small- and large-scale reagent preparation and full assay development, Revvity is the ideal trusted partner to support all your needs.

Keep your therapeutic research moving forward with faster, more robust, physiologically relevant results. Call us or visit www.revvity.com.



revvity