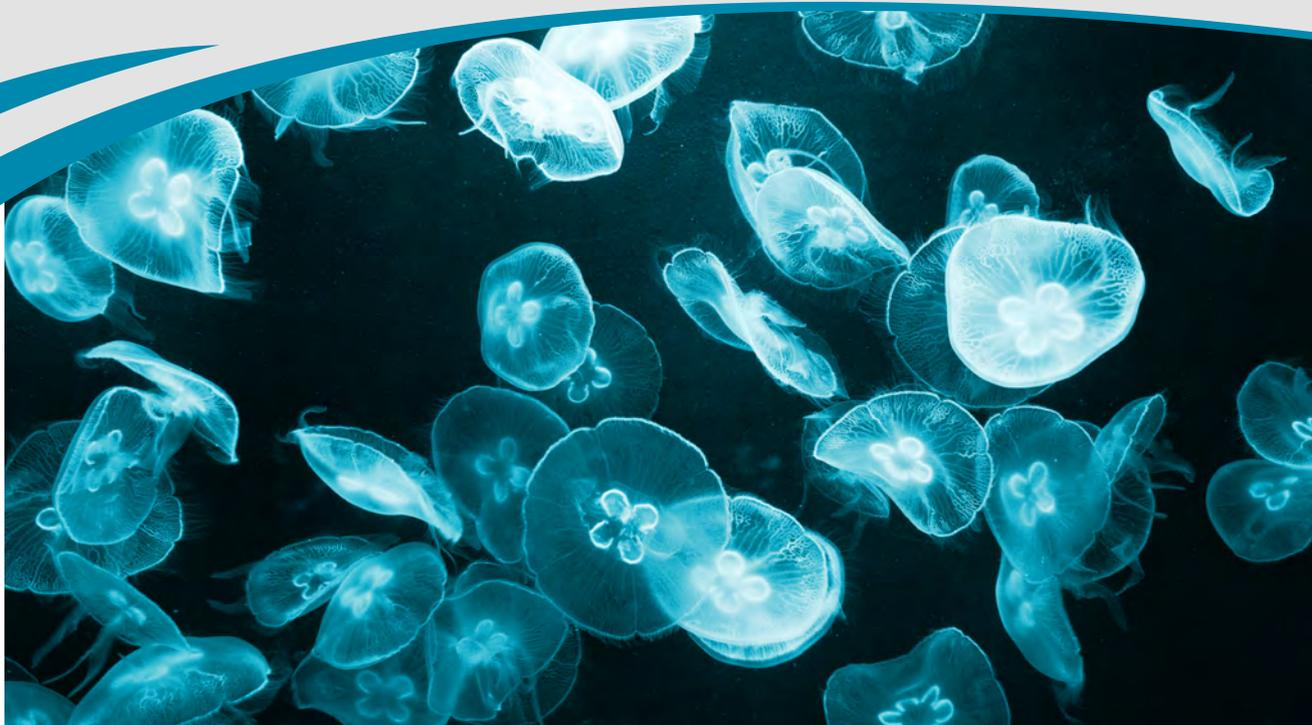


# Tag-lite<sup>®</sup> services to evaluate your Biosimilar or Biobetter samples



BIOTHERAPEUTICS

## THE MOST EFFECTIVE WAY TO FIND THE ONE YOU'RE LOOKING FOR

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 Fc $\gamma$  receptors involved in T cell mediated ADCC (Fc $\gamma$ RI, Fc $\gamma$ RII and Fc $\gamma$ RIII) and to FcRn.

Cisbio's binding profile services will characterize your samples on all Fc $\gamma$  and FcRn receptors (CD32A, CD32B, CD64, and CD16) using our cell-based assays.

### Benefit from the expertise of Cisbio binding service to profile and study your Fc $\gamma$ 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

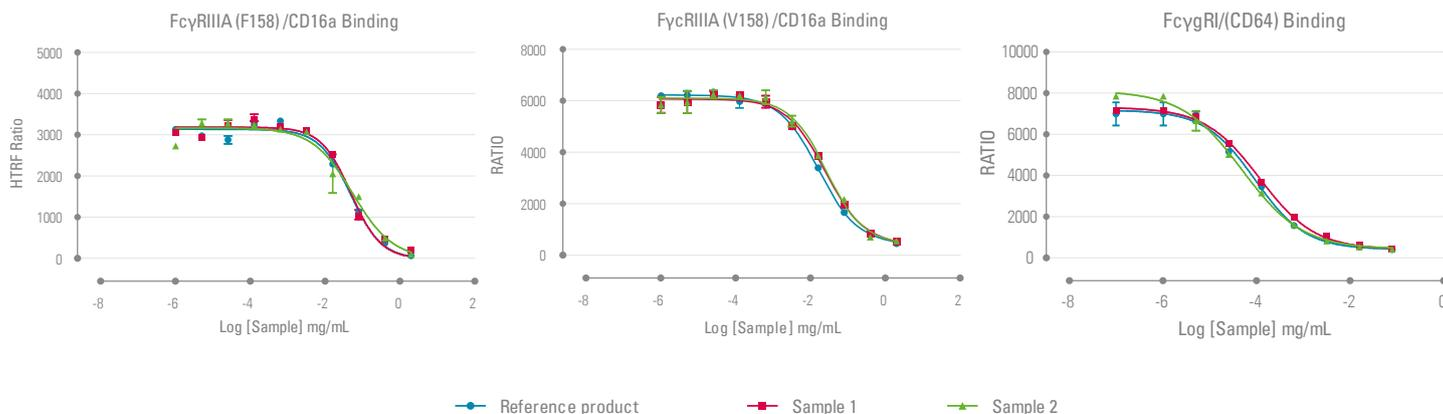


Cisbio's team of experienced experts works hand-in-hand with you to characterize your products

## HOW IT WORKS PROFILING

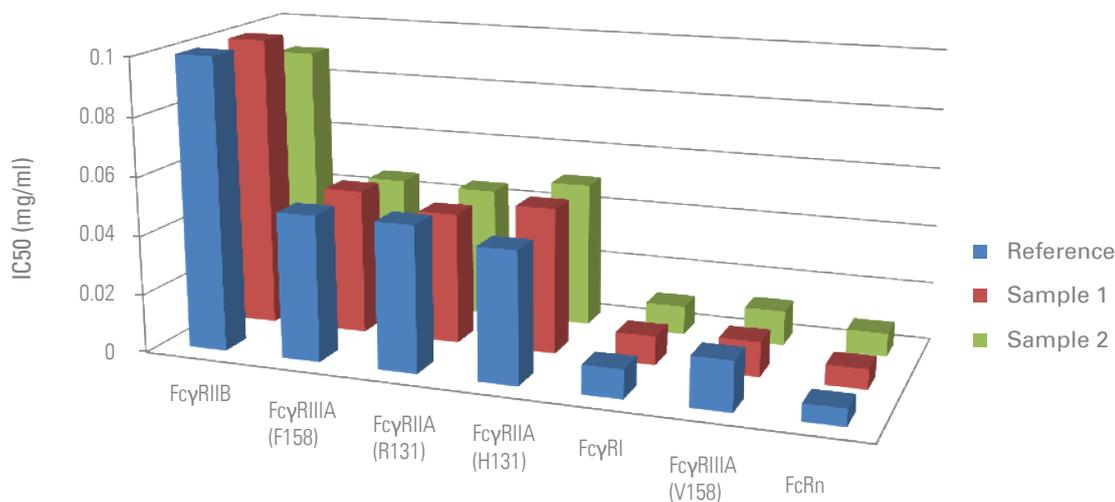
In this case study, 3 samples (2 samples and 1 reference product) were assessed for binding affinity to FcγRI, FcγRIIA, FcγRIIB, FcγRIIIA and FcRn using the Cisbio cell-based platform.

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, Cisbio'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, Cisbio 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.cisbio.com/tag-lite-binding-assays](http://www.cisbio.com/tag-lite-binding-assays)

### FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567

China +86 10 8080 9288 Japan +81-(0)43-306-8712

Visit [www.cisbio.com](http://www.cisbio.com) to find a list of our regional distributors



cisbio