

# Alpha technologies for antibody detection and characterization.

## Alpha technology: Biotherapeutics applications

### Introduction

Therapeutic proteins, namely antibodies, are growing in importance in many drug discovery pipelines. Since the last decade, many therapeutic antibodies were approved by the FDA (Table 1)<sup>1</sup>.

Many laboratories developing and producing antibodies still rely on traditional enzyme-linked immunosorbent assay (ELISA) to perform clonal selection and characterization despite the fact that this proven technology often suffers

### Key benefits

- Consistent, reproducible results
- Simplified workflow
- Fast antibody detection and characterization

from lack of sensitivity and reproducibility due to its heterogenous nature. The numerous wash and blocking steps required to perform ELISA assays also makes them particularly difficult to automate. This process suffers from low throughput and potential high affinity and selective antibodies can be missed.

Table 1: Example FDA approved therapeutic monoclonal antibodies.

Antibody	Brand name	Type	Indication
Abciximab	ReoPro®	chimeric	Cardiovascular disease
Adalimumab	Humira®	Human	Auto-immune disorders
Alemtuzumab	Campath®	humanized	Chronic lymphocytic leukemia
Basiliximab	Simulect®	chimeric	Transplant rejection
Bevacizumab	Avastin®	humanized	Colorectal cancer, Age related macular degeneration
Cetuximab	Erbix®	chimeric	Colorectal cancer, Head and neck cancer
Certolizumab pegol	Cimzia®	humanized	Crohn's disease
Daclizumab	Zenapax®	humanized	Transplant rejection
Eculizumab	Soliris®	humanized	Paroxysmal nocturnal hemoglobi nuria
Efalizumab	Raptiva®	humanized	Psoriasis
Gemtuzumab	Mylotarg®	humanized	Acute myelogenous leukemia
Ibritumomab tiuxetan	Zevalin®	Murine	Non-Hodgkin lymphoma
Infliximab	Remicade®	chimeric	Several autoimmune disorders

Table 1: Example FDA approved therapeutic monoclonal antibodies. (continued)

Antibody	Brand name	Type	Indication
Muromonab-CD3	Orthoclone® OKT3	Murine	Transplant rejection
Natalizumab	Tysabri®	humanized	Multiple sclerosis and Crohn's disease
Omalizumab	Xolair®	humanized	Allergy-related asthma
Palivizumab	Synagis®	humanized	Respiratory Syncytial Virus
Panitumumab	Vectibix®	Human	Colorectal cancer
Ranibizumab	Lucentis®	humanized	Macular degeneration
Rituximab	Rituxan®, Mabthera®	chimeric	Non-Hodgkin lymphoma
Tositumomab	Bexxar®	Murine	Non-Hodgkin lymphoma
Trastuzumab	Herceptin®	humanized	Breast cancer

Over the years, Alpha technology (i.e. AlphaScreen™ and AlphaLISA™) became an established detection technology in many academic and industrial laboratories. Alpha technology is homogeneous and non-radiometric with distinct features that makes it enabling in comparison to other proximity assays. These features include: high sensitivity, reproducibility, robustness and large distance for proximity. The latter is allowed by a unique mode of energy transfer prevailing between Donor and Acceptor beads based on singlet oxygen. Singlet oxygen lifetime allows this molecule to travel up to 200 nm in solution before decaying to its ground state, hence offering an unsurpassed equivalent distance for proximity. Such a large distance for proximity allows one to capture very large molecules and study intractable interactions while using either simple or complex assay configurations, a task which is often difficult to accomplish using other popular homogeneous proximity technologies such as TR-FRET.

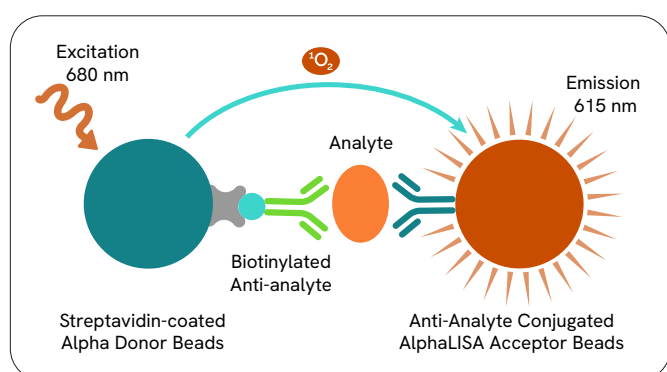


Figure 1: Principles of AlphaLISA bridging assays. Streptavidin-coated Donor beads are used to capture a biotinylated antibody specific to a precise epitope present on the analyte. Another antibody, also specific to an epitope found on the analyte, is directly conjugated to the Acceptor beads. Both the Donor and Acceptor beads are brought into proximity in the presence of the analyte. An AlphaLISA signal, directly proportional to the concentration of the analyte in solution, is then generated after laser excitation at 680 nm.

Based on these benefits, Alpha technologies represent powerful means of detecting and characterizing a wide range of proteins including antibodies. An extensive toolbox was thus developed for this purpose. This toolbox include various beads coated with a series of anti-tag antibodies, anti-species antibodies and other generic capture proteins such as protein A and G. Using these tools, researchers are now capable of converting their traditional ELISAs into more sensitive, reproducible and higher throughput assays; therefore facilitating therapeutic protein discovery and characterization.

The following review will describe how Alpha technologies can be used to enable antibody selection and characterization.

### Antibody clonal selection

Vainshtein et al<sup>2</sup> used AlphaScreen to improve the automation of the antibody clonal selection process and then speed up the discovery of the best IgG clones (Figure 2). Using AlphaScreen toolbox reagents, the authors developed a broad variety of applications for hybridoma fusion screens, serum titers and antibody quantifications. Competitive immunoassays were also performed to determine the affinity of the IgG clones identified during screens. The authors state that AlphaScreen proved to be more sensitive than ELISA and allowed for identification of a larger number of hits. Miniaturization and the homogeneous nature of the assay simplified the work flow, significantly shortened assay run time and made the screening operations more cost-effective.

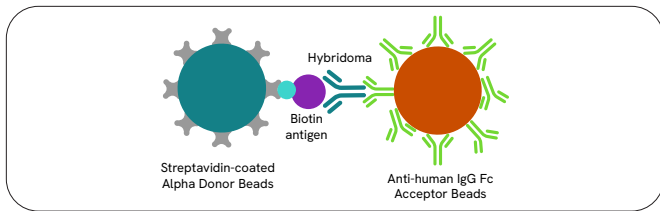


Figure 2: To perform hybridoma screening, Vainshtein et al used anti-human IgG Fc-coated Acceptor beads and Streptavidin-coated Donor beads to capture a biotinylated derivative of the target antigen of interest. In the presence of antibodies specific to that antigen, both beads are brought into proximity and a robust Alpha signal is emitted.

The authors also compared the performance of AlphaScreen to that of ELISA during hybridoma screening (Figure 3). A total of 148 samples were screened: 23 and 55 positives were identified with ELISA and AlphaScreen respectively. Interestingly, all ELISA hits were a subset of the Alpha hits. AlphaScreen identified more hits due to higher sensitivity due to the detection of low affinity antibodies.

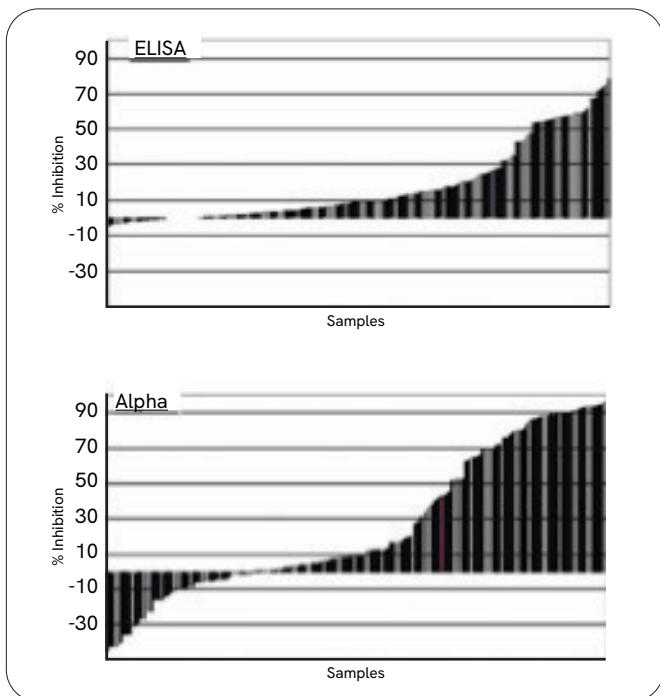


Figure 3: Hit comparison between ELISA and AlphaScreen. Due to its higher sensitivity, AlphaScreen generated more hits.

### Measuring antibody affinity

Performing molecular biology re-engineering allows one to modify antibody Fc or Fab sequences to enhance their specificity and/or affinity to their targeted antigen. Using a combination of computational structure-based protein design methods coupled with high-throughput protein screening, Lazar et al<sup>3</sup> modified Fc sequences of human IgGs to obtain higher affinity antibodies binding to Fc $\gamma$  receptors. Affinity of the resulting Fc variants were estimated with AlphaScreen in competition assays where a biotinylated derivative of the parent antibody (trastuzumab) was used as the tracer.

It is worth mentioning that competition assays are the best means to measure affinities with AlphaScreen. As predicted by the Cheng and Prusoff<sup>4</sup> equation, when both proteins used as tracers and targets are used in negligible concentrations compared to their putative Kd values, IC<sub>50</sub> values obtained from competition assays will closely match the Kd. Affinities obtained with AlphaScreen closely matched those estimated with surface Plasmon resonance (SPR).

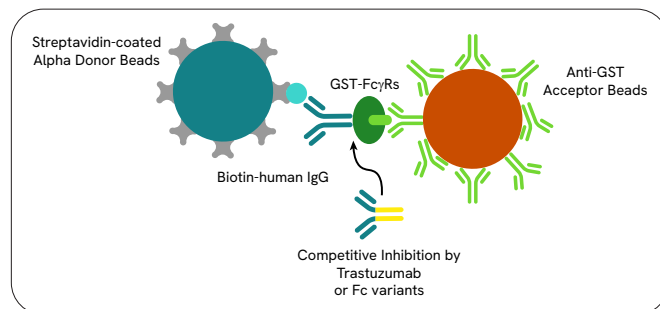


Figure 4: To measure the affinity of different IgG clones bearing Fc variants, Lazar et al. developed a competition assay using biotinylated trastuzumab as a tracer. Anti-GST coated Acceptor beads are used to capture recombinant Fc $\gamma$  receptors expressed as GST-fusion proteins. In absence of Fc variants, biotinylated trastuzumab binds to its receptor so both the Donor and Acceptor beads are brought into proximity to produce a signal. In the presence of increasing Fc variant concentrations, a progressive signal decrease is observed.

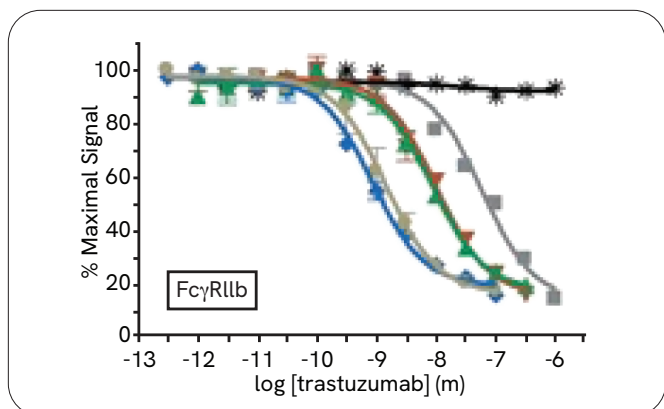


Figure 5: Examples of Fc variant competition isotherms. Black asterisk, buffer; gray squares, WT; black diamonds, S298A/E333A/K334A; green triangles, S239D; red inverted triangles, I332E; blue diamonds, S239D/I332E; and tan circles, S239D/I332E/A3301.

Using AlphaScreen, the authors identified and characterized a series of Fc variants with optimized Fc $\gamma$  receptor affinity and specificity. These variants showed more than 2 orders of magnitude enhancement of *in vitro* effector function. They were efficacious against cells expressing low levels of target antigen, producing increased response in an *in vivo* preclinical model. Re-engineering Fc regions offer a means to improve therapeutic antibodies and have the potential to broaden the diversity of antigens that can be targeted for antibody-based tumor therapy.

### Selecting & matching antibody pairs for immuno-sandwich assays

Careful antibody selection is mandatory to select the best pair to produce highly sensitive and selective sandwich immunoassays. The identification of pairing antibodies is a tedious and laborintensive process. Using protein A coated Donor and Acceptor beads, Bembenek et al<sup>5</sup> developed a robust and high-throughput method for identifying pairing complementary antibodies derived either from commercial sources or identified during a rabbit hybridoma monoclonal screening. This group demonstrated the value and effectiveness of their assay with different protein targets, including: Akt2, ATF3, and NAE $\beta$  (the  $\beta$ -subunit of the neddylation activation enzyme).

Using a matrix approach, the authors systematically tested a wide range of antibody samples against each other by pre-coating them on both Donor and Acceptor beads (Figure 6). To make sure that the beads were optimally loaded with antibodies, saturating amounts of commercial IgGs or hybridoma supernatants were mixed with beads. Beads were then centrifuged and washed multiple times to remove unbound antibodies.

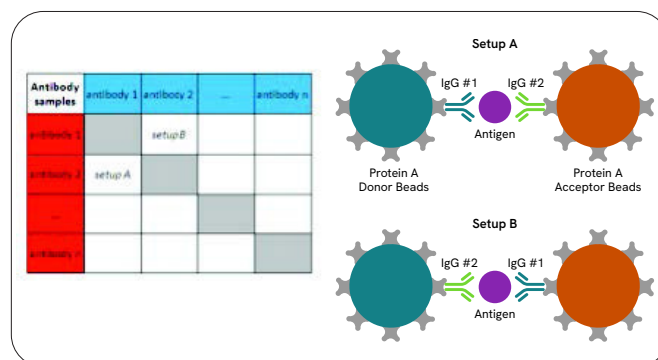


Figure 6: Left: Antibody samples (hybridomas) were tested in pairs using a matrix approach. Right: Protein A coated Donor and Acceptor beads were first pre-loaded with distinct antibody samples. Each antibody combination was tested on both Donor and Acceptor bead to find the optimal assay configuration (i.e. setup A or B). Gray shaded boxes represent assay setups where the same antibody is found on both Donor and Acceptor beads. When the antibody sample is of monoclonal nature, no significant antigen capture is possible and therefore only background signal is measured.

Representative matrix design for 18 positive anti-NAE $\beta$  rabbit monoclonal antibody-producing hybridoma clones is shown on figure 7. Results represent averages of duplicate samples with complementary antibody pairs. Positives are highlighted in yellow boxes while same antibodies tested against themselves are highlighted in gray along the diagonal. The average background values obtained across the entire matrix in the absence of antigen was approximately 1000 counts. In that example, clones 18 and 81 were those allowing for the most efficient antibody pairing. Based on total counts (155,000 cps), best setup was obtained when clones 87 and 81 were immobilized on Donor and Acceptor beads respectively. Interestingly, only 60,000 counts were generated when the same antibodies were switch from Donor to Acceptor beads and vice-versa. Difference in levels of protein A conjugation or antibody capture capacity between the Donor and the Acceptor beads may account for the difference.

Hybridoma number	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30	31	32	33	34	35				
1	4825	4548	4877	3845	3939	4189	468	4336	2764	4377	3906	4539	396	3948	2367	4674	4387	4335																					
2	4748	4772	2309	3886	2252	2521	4957	4872	4836	4814	2254	4784	4276	4848	2893	4845	4635	2454																					
3	486	436	4387	4817	486	4575	4845	461	2904	4847	436	222	484	4742	4797	422	437	4287																					
4	4383	4247	4867	432	4818	473	4855	2389	2141	436	3173	2888	401	438	4848	424	441	488																					
5	481	4855	484	4825	4826	4825	458	4864	2518	4813	4883	475	458	4842	4812	4382	478	438																					
6	2267	4307	4767	461	4475	488	427	4813	441	4814	4888	2123	468	471	4811	2714	2798	2188																					
7	4358	4387	4813	475	4745	4792	484	484	4331	468	4882	4874	474	481	4838	437	465	4887																					
8	4314	4216	4848	487	4845	4458	478	4338	4857	4216	4473	4358	469	425	4388	437	488	4236																					
9	474	4279	4798	486	4818	4814	422	4814	469	2714	4888	486	448	486	4311	4318	4473	48818																					
10	4342	4382	4888	484	4818	478	478	4373	4869	468	4672	4768	486	438	4876	4883	4338	4574																					
11	4811	4853	2384	3838	4425	4647	478	4381	2484	42178	4395	4454	4647	4311	2718	438	422	4328																					
12	4392	4387	4888	482	4545	2287	488	4387	4385	2818	4425	4634	484	438	4811	4845	481	4843																					
13	4887	444	4587	46	4413	4813	427	488	481	4867	4883	4378	462	431	4318	4826	4318	4338																					
14	4799	4384	4788	2878	4718	4411	4855	24318	4634	4814	4748	4878	431	486	4311	4311	4311	4311																					
15	479	4345	4813	4818	4881	4841	4528	465	4888	2784	438	4817	4378	4811	4811	4811	4314	4844																					
16	4415	4463	4377	4818	4811	2738	427	437	4789	438	4388	4388	431	431	4818	438	438	4843																					
17	4313	4178	4844	474	4848	2422	4836	433	4413	4718	4378	4312	437	4313	4818	4318	4318	4844																					
18	4338	4415	4387	4848	444	4314	481	4528	4818	4818	4781	4888	4312	4677	4384	438	4811	4842																					

Figure 7: Representative matrix design for 18 positive anti-NAEβ rabbit monoclonal antibody-producing hybridoma clones.

Utilizing and characterizing antibodies present in complex matrices

During the course of antibody development, the detection and characterization of specific clones is complicated by the presence of non-specific antibodies. This is the case for polyclonal antibodies generated in animals where only a very small proportion of the antibodies present in serum samples are specific to an antigen of interest. Sometimes, monoclonal antibodies also represent a very minor proportion of specific antibodies in ascites fluids.

ELISAs allow one to isolate and characterize clones of interest by immobilizing the antigen of interest onto the surface of a solid matrix (i.e. gel or microplate). After a series of blocking and wash steps, non-specific antibodies are washed out leaving specific antibodies bound to the solid support and available for subsequent detection. Once again, this process suffers from low throughput and potential high affinity and selective antibodies can be missed if a sufficient amount of samples can't be screened.

To speed up the characterization of low abundance antibodies in serum samples, Poulsen et al<sup>6</sup>. developed a semi-homogeneous approach using AlphaLISA. As shown in Figure 8, a 5-step protocol was developed and performed as follows:

- Acceptor beads conjugated with antibodies specific to the analyte of interest are first immobilized onto the wells of "high binding" capacity microplates
- Antigen samples are then added followed by a single wash step

- The antibody to characterize and present in a complex matrix (e.g. serum, ascites) is added followed by another single wash step
- Donor beads conjugated with an antibody (or any other binding protein) specific to the antibody to characterize are finally added
- Microplates are incubated and read

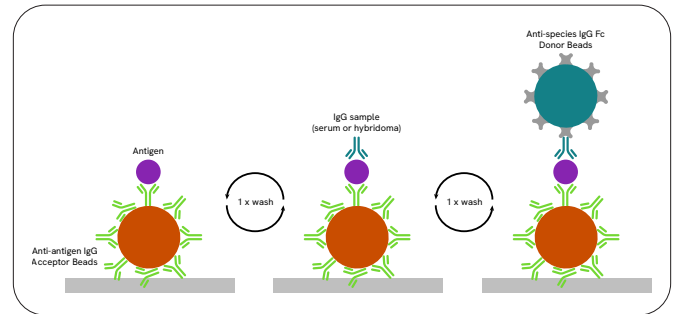


Figure 8: AlphaLISA semi-wash assay configuration used by Poulsen et al to characterize low abundance antibodies present in serum samples.

As shown in Figure 9, the semi-wash AlphaLISA protocol allows one to detect low levels of specific antigens using unpurified primary IgG preparations from serum or hybridoma. In this specific case, insulin was detectable with a LOD of 0.3 pM while the standard homogeneous protocol was less sensitive; allowing one to detect insulin with a LOD of 50 pM.

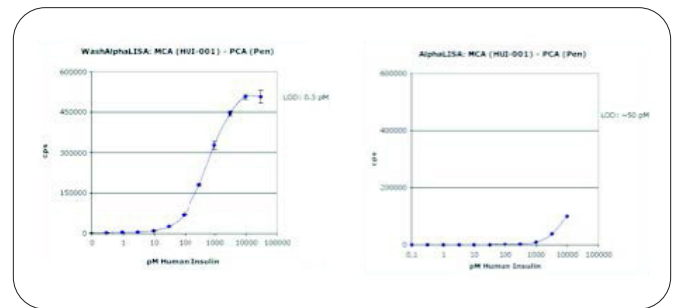


Figure 9: Characterization of an unpurified polyclonal antibody against insulin (PCA-Pen) using AlphaLISA. The performance of the AlphaLISA semi-wash protocol (wash AlphaLISA, left panel) was compared to that of the standard homogeneous AlphaLISA protocol (right panel). The semi-wash protocol allowed for a better performance compared to the standard protocol: key parameters such as the lowest limit of detection (LOD) and dynamic range were improved by at least 2 orders of magnitude.

Performing semi-wash assays with Alpha technologies is one way to expand their use and versatility. Polyclonal antibody preparations with low content of specific IgG may be characterized without the needs of performing affinity chromatography. Alpha assays developed with this alternative still show broad dynamic range and excellent sensitivity while no hook effect due to antigen oversaturation is observed. It is worth mentioning that Alpha homogeneous and semi-wash assays can be run with exactly the same reagents and equipment (liquid handler and reader).

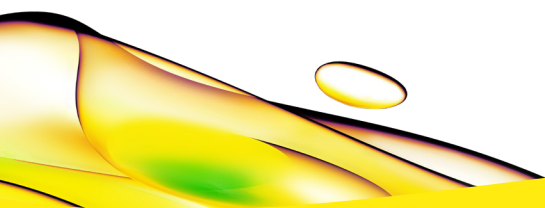
## Summary

Antibodies represent an emerging class of therapeutic drugs and clinical tools. Higher throughput and more sensitive technologies are needed to speed up antibody selection and characterization. The versatility of AlphaScreen and AlphaLISA makes these technologies very appropriate to detect, measure and characterize antibodies present in different matrices.

Using AlphaScreen or AlphaLISA reagents, scientists showed that the antibody clonal selection process could be significantly improved. Other groups used competitive immunoassay setups and demonstrated that it was possible to determine antibody affinities with values closely matching those obtained using SPR. Using Alpha toolbox reagents ([www.revvy.com](http://www.revvy.com)), straightforward protocols were developed to select and match antibody pairs for immuno-sandwich assays. Alternate semi-wash protocols were also designed to utilize and characterize antibodies present in complex matrices such as serum or ascites fluid, expanding the usability of Alpha technologies further.

## References

1. Waldmann, TA. Immunotherapy: past, present and future. *Nat. Med.* 9 (3) 2, 269-277 (2003).
2. Vainshtein I, Kurose S, Vickroy J, Russell Grove J, Liang M. Fast and efficient soluble antigen-antibody screening using homogeneous AlphaScreen assay. Poster Presentation. SBS 11th Annual Conference. September 2004, Orlando (USA)
3. Lazar GA, Dang W, Karki S, Vafa O, Peng JS, Hyun L, Chan C, Chung HS, Eivazi A, Yoder SC, Vielmetter J, Carmichael DF, Hayes RJ, Dahiyat BI. Engineered antibody Fc variants with enhanced effector function. *Proc Natl Acad Sci.* 103 (11):4005-1(2006)
4. Cheng Y, Prusoff WH. Relationship between the inhibition constant (K<sub>1</sub>) and the concentration of inhibitor which causes 50 percent inhibition (I<sub>50</sub>) of an enzymatic reaction. *Biochem Pharmacol.* 22(23):3099-108 (1973).
5. Bembenek ME, Burkhardt A, Ma J, Li Z, Loke HK, Wu D, Xu Q, Tayber O, Xie L, Li P, Li L. *Anal Biochem.* Determination of complementary antibody pairs using protein A capture with the AlphaScreen assay format. 408(2):321-7 (2011).
6. Poulsen F. Wash AlphaLISA-an extension of the AlphaLISA/ LOCI technology. Revvity Technology Forum - University of Copenhagen, Denmark October 6th 2010.



revvity