

Key Benefits

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

Alpha Technologies for Antibody Detection and Characterization

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)¹.

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 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.

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 hemoglobinuria
Efalizumab	Raptiva®	humanized	Psoriasis
Gemtuzumab	Mylotarg®	humanized	Acute myelogenous leukemia
Ibritumomab tiuxetan	Zevalin®	Murine	Non-Hodgkin lymphoma
Infliximab	Remicade®	chimeric	Several autoimmune disorders
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

Table 1. Example FDA approved therapeutic monoclonal antibodies.

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.

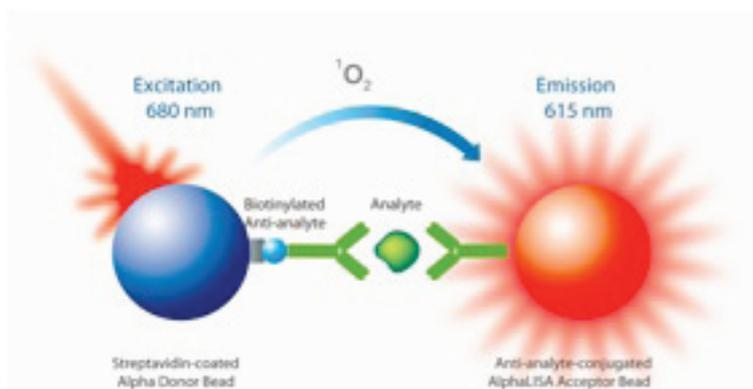


Figure 1. Principles of AlphaLISA bridging assays. Streptavidin-coated Donor beads are used to capture a biotinylated antibody specific to a precise epitope present on the analyte. Another antibody, also specific to an epitope found on the analyte, is directly conjugated to the Acceptor beads. Both the 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*² 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.

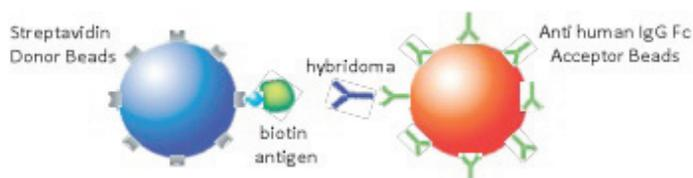


Figure 2. To perform hybridoma screening, Vainshtein *et al* used anti-human IgG Fc-coated Acceptor beads and Streptavidin-coated Donor beads to capture a biotinylated derivative of the target antigen of interest. In the presence of antibodies specific to that antigen, both beads are brought into proximity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

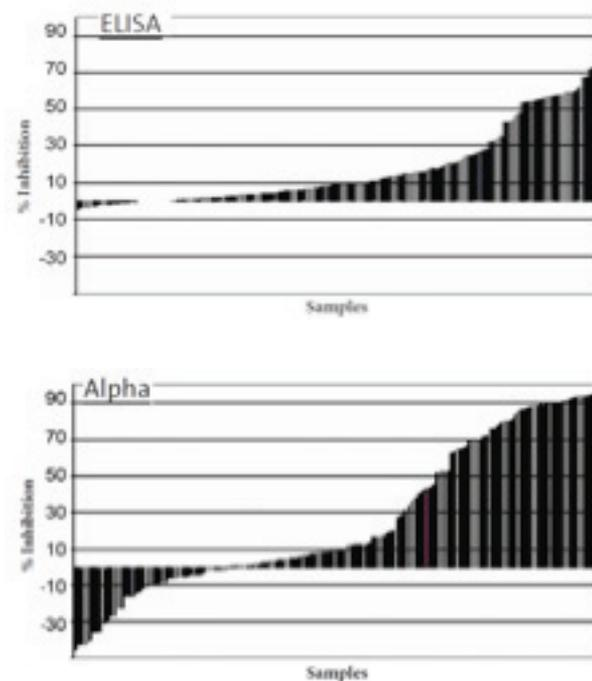


Figure 3. Hit comparison between ELISA and AlphaScreen. Due to its higher sensitivity, AlphaScreen generated more hits.

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*³ modified Fc sequences of human IgGs to obtain higher affinity antibodies binding to Fc γ 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 Prusoff⁴ equation, when both proteins used as tracers and targets are used in negligible concentrations compared to their putative K_d values, IC₅₀ values obtained from competition assays will closely match the K_d. Affinities obtained with AlphaScreen closely matched those estimated with surface Plasmon resonance (SPR).

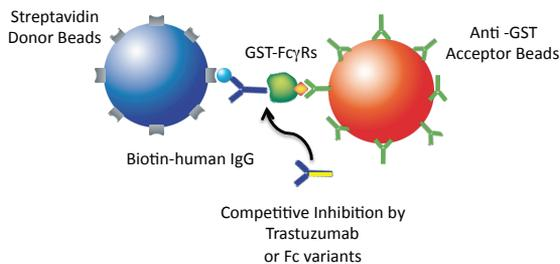


Figure 4. To measure the affinity of different IgG clones bearing Fc variants, Lazar *et al.* developed a competition assay using biotinylated trastuzumab as a tracer. Anti-GST coated Acceptor beads are used to capture recombinant Fc γ receptors expressed as GST-fusion proteins. In absence of Fc variants, 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.

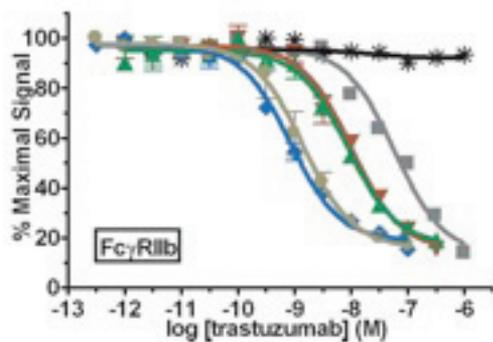


Figure 5. Examples of Fc variant competition isotherms. Black asterisk, buffer; gray squares, WT; black diamonds, S298A/E333A/K334A; green triangles, S239D; red inverted triangles, I332E; blue diamonds, S239D/I332E; and tan circles, S239D/I332E/A3301.

Using AlphaScreen, the authors identified and characterized a series of Fc variants with optimized Fc γ 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 labor-intensive process. Using protein A coated Donor and Acceptor beads, Bembenek *et al*⁵ 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 β (the β -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.

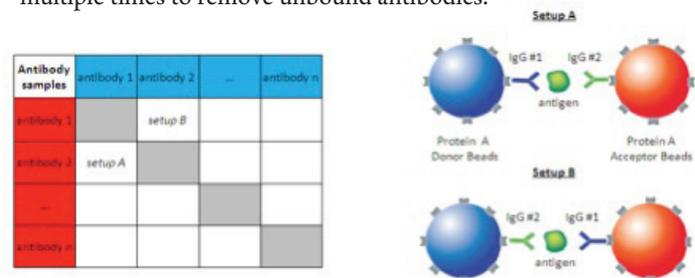


Figure 6. Left: Antibody samples (hybridomas) were tested in pairs using a matrix approach. Right: Protein A coated Donor and Acceptor beads were first pre-loaded with distinct antibody samples. Each antibody combination was tested on both Donor and Acceptor bead to find the optimal assay configuration (ie. 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 β 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	4	9	14	18	25	28	31	32	39	57	79	81	87	97	99	99
1	1835	858	1577	5966	1919	5180	968	8330	27864	13177	8986	9539	950	10045	2567	9878	1087	1425
2	1748	8777	2309	3386	2252	25913	8957	8072	14336	4014	2214	6791	1236	13636	2893	8845	1638	2451
4	884	836	1397	4817	846	16775	1045	969	29004	96897	958	222	884	96142	1997	122	817	1287
9	2083	4247	3867	732	3838	713	8055	2280	2147	836	3173	2888	931	1138	5828	614	141	4038
14	903	8955	9864	4826	1626	46520	958	1084	25916	80013	8883	175	950	32853	5492	1582	158	1188
18	22871	83070	37867	903	14476	108	627	10103	5443	10314	20940	23237	968	7953	148713	2764	2096	17369
25	1958	8287	6893	625	1245	1270	894	684	1131	560	8883	6874	874	1369	9538	117	945	1887
28	1831	8216	1869	1197	9845	3438	778	8330	8957	1216	1473	1159	969	2252	1368	127	188	1236
31	7634	12797	12938	1486	10878	5824	102	4094	365	2214	8898	7186	1148	1036	1391	1038	1473	10858
32	4347	7382	7688	881	5938	1878	778	3785	3869	960	5673	4719	846	2518	18748	5883	1938	4378
39	1891	8853	2386	5138	1625	16875	798	1391	20849	12178	8395	1434	8845	13511	2256	198	122	1929
57	1492	1587	1588	8812	1549	2387	988	4387	4385	2638	1435	1634	884	4788	8991	8845	183	1843
79	8987	1444	1587	968	1463	4893	127	988	8817	8882	8883	1378	1622	1011	1038	1626	1438	1536
81	25293	71464	78952	2479	67578	14387	8095	24319	16264	4878	57546	14879	1311	7884	175178	13172	2679	10143
87	979	8145	1482	8710	1083	168372	4528	905	10985	27184	1319	8817	1378	14872	1886	1574	1644	1739
97	1615	1463	1577	1159	1615	2796	107	817	1719	1568	1638	1318	1031	2471	1039	1348	1636	1443
99	1311	1178	1444	1214	1348	2122	1036	1131	1158	1178	1112	1127	1013	1119	1119	1119	1119	1149
99	1338	1425	1387	7186	1436	47544	783	1529	48388	18736	1791	1036	1132	14717	1264	158	1893	1482

Figure 7. Representative matrix design for 18 positive anti-NAE β 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.*⁶ 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

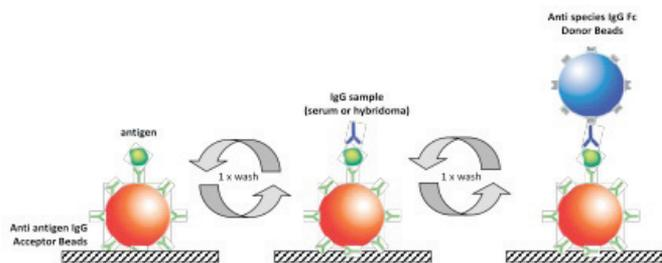


Figure 8. AlphaLISA semi-wash assay configuration used by Poulsen *et al* to characterize low abundance antibodies present in serum samples.

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.

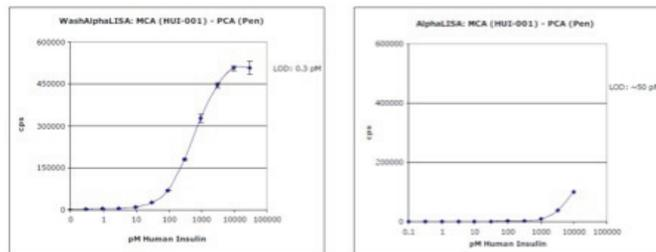


Figure 9. Characterization of an unpurified polyclonal antibody against insulin (PCA-Pen) using AlphaLISA. The performance of the AlphaLISA semi-wash protocol (wash AlphaLISA, left panel) was compared to that of the standard homogeneous AlphaLISA protocol (right panel). The semi-wash protocol allowed 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.perkinelmer.com/AlphaToolbox), 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_1) and the concentration of inhibitor which causes 50 percent inhibition (IS_0) 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. PerkinElmer Technology Forum – University of Copenhagen, Denmark October 6th 2010.