



## References to Fortebio's Products in Published Literature.

### **2009**

A Global Benchmark Study Using Affinity-based Biosensors.

Rich, R.L. and others,

*Analytical Biochemistry*, 2009, 386 (2), 194-216.

Exploring Blocking Assays Using Octet, ProteOn, and Biacore Biosensors.

Abdiche, Y. N.; Malashock, D. S.; Pinkerton, A; Pons, J.

*Analytical Biochemistry*, 2009, 386(2), 172-180.

High Affinity Human Antibodies to Human Nerve Growth Factor.

MacDonald, L.; Torres, R.; Morra, M. R.; Martin, J. H.; Reinhardt, J. C.

Regeneron Pharmaceuticals Inc., United States Patent Application

20090041717.

Antibodies to IL-6 and Use Thereof.

Garcia-Martinez, L.; Jensen, A. E. C.; Olson, K.; Dutzar, B.; Ojala, E.;

Kovacevich, B.; Latham, J.; Smith, J. T. L.

Alder Biopharmaceuticals Inc., United States Patent Application 20090028784.

Antibodies to TNF Alpha and Use Thereof.

Olson, K.; Dutzar, B.; Jensen, A. E. C.; Garcia-Martinez, L.; Ojala, E.; Latham, J.

Alder Biopharmaceuticals Inc., United States Patent Application 20090022659.

Human Antibodies to Human Delta-like Ligand 4.

Papadopoulos, N. J.; Martin, J. H.; Smith, E.; Noguera-Troise, I.; Thurston, G.

Regeneron Pharmaceuticals, Inc., United States Patent Application

20090017035.

Human Antibodies to Human Delta-like Ligand 4.

Papadopoulos, N. J.; Martin, J. H.; Smith, E.; Noguera-Troise, I.; Thurston, G.

Regeneron Pharmaceuticals, Inc., United States Patent 7,488,806.

### **2008**

Probing the Binding Mechanism and Affinity of Tanezumab, a Recombinant Humanized Anti-NGF Monoclonal Antibody, Using a Repertoire of Biosensors.

Abdiche, Y. N.; Malashock, D. S.; Pons, J.

*Protein Science*, 2008, 17(8), 1326-1335.



Determining Kinetics and Affinities of Protein Interactions Using a Parallel Real-time Label-free Biosensor, the Octet.

Abdiche, Y.N.; Malashock, D. S.; Pinkerton, A; Pons, J.  
*Analytical Biochemistry*, 2008, 377(2), 209-217.

A Rapid Method for Determining Dynamic Binding Capacity of Resins for the Purification of Proteins.

Do, T.; Ho, F.; Heidecker, B.; Witte, K.; Chang, L.; Lerner, L.  
*Protein Expression and Purification*, 2008, 60, 147–150.

The Surface Protein TIGIT Suppresses T Cell Activation by Promoting the Generation of Mature Immunoregulatory Dendritic Cells.

Yu, X.; Harden, K.; Gonzalez, L. C.; Francesco, M.; Chiang, E.; Irving, B.; Tom, I.; Ivelja, S.; Refino, C. J.; Clark, H.; Eaton, D.; Grogan, J. L.  
*Nature Immunology*, 2008, 10, 48-57.

Advantages and Application of Label-free Detection Assays in Drug Screening.

Cunningham, B. T.; Laing, L. G.  
*Expert Opinion on Drug Discovery*, 2008, 3(8), 891-901.

Humaneered Anti-factor B Antibody.

Emlen, W.; Holers, V. M.; Flynn, P.  
Taligen Therapeutics, Inc., United States Patent Application 20080299114.

Human Antibodies to Human Delta-like Ligand 4.

Papadopoulos, N. J.; Martin, J. H.; Smith, E.; Noguera-Troise, I.; Thurston, G.  
Regeneron Pharmaceuticals, Inc., United States Patent Application 20080181899.

High Affinity Human Antibodies to Human IL-4 Receptor.

Stevens, S.; Huang, T. T.; Martin, J. H.; Farihurst, J. L.; Rafique, A.; Torres, M.; Pobursky, K. J.; Leidich, R. W.; Windsor, J. A.; Mikulka, W. R.; Ahrens, D. M.; Shi, E.; Papadopoulos, N. J.  
Regeneron Pharmaceuticals, Inc., United States Patent Application 20080160035.

WNT Antagonists and Their Use in the Diagnosis and Treatment of WNT-mediated Disorders.

Ernest, J. A.; Polakis, P.; Rubinfeld, B.; Dealmeida, V. I.  
Genentech, Inc., United States Patent Application 20080299136.



## **2007**

Detecting DNA: Getting and Begetting Cancer.

Farkas, A. M.; Kilgore, T. M.; Lotze, M. T.

*Current Opinion in Investigational Drugs*, 2007, 8, 981-986.

Culture Method for Obtaining a Clonal Population of Antigen-specific B Cells.

Jensen, A. E. C.; Garcia, L.; Ojala, E.; Latham, J.

Alder Biopharmaceuticals, Inc., United States Patent Application 20070269868.

Label-free Optical Biosensors in Drug Discovery.

Fang, Y.

*Trends in Biopharmaceutical Industry*, 2007, 3, 34-38.

Higher-throughput, label-free, real-time molecular interaction analysis.

Rich, R. L.; Myszka, D. G.

*Analytical Biochemistry*, 2007, 361, 1-6.

## **2006**

Optical Biosensors: Where Next and How Soon?

Cooper, M. A.

*Drug Discovery Today*, 2006, 11 (23-24), 1061-1067.

Current Biosensor Technologies in Drug Discovery.

Cooper, M. A.

*Drug Discovery World*, 2006, 68-82.