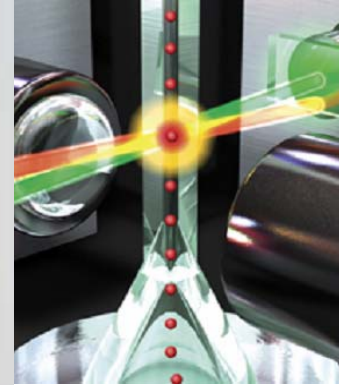


Sherry A. Dunbar, Ph.D.
sdunbar@luminexcorp.com

Luminex

The **x**MAP[®] Technology Platform

FLEXMAP^{3D}
THE NEXT DIMENSION IN MULTIPLEXING

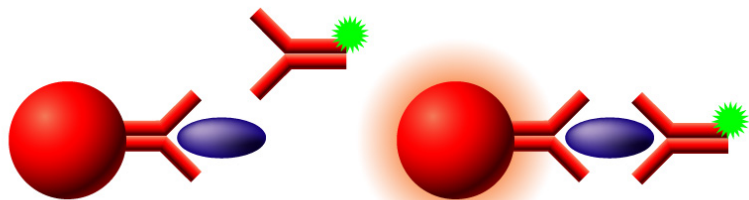


Luminex

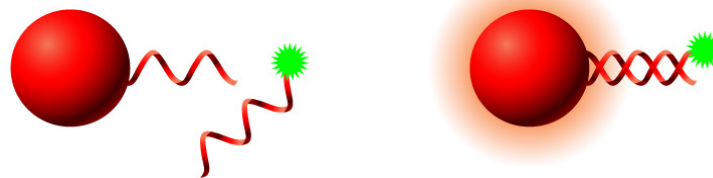
- Applications and Menu



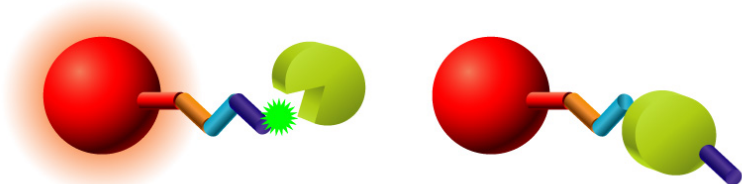
Immunoassay



Nucleic Acid Assay



Enzyme Assay



Receptor-Ligand



● Protein Expression Profiling

● Cancer Markers

- AFP, CA 15-3, CA 19-9, CA 125, CEA, PSA ...

● Cardiac Markers

- APO A1, APO E, Fibrinogen, Adiponectin ...

● Cell Signaling

- Akt, c-Jun, NF-kappaB ...

● Cytokines, Chemokines, Growth Factors

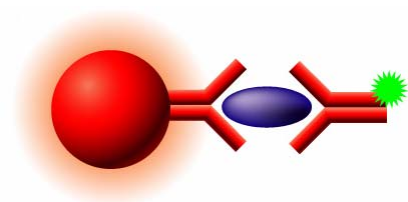
- ILs, IFN, TNF, MIP, RANTES ...

● Endocrine

- Amylin, Glucagon, Insulin, Leptin ...

Novagen®

RULES • BASED MEDICINE, INC.



- Protein Expression Profiling
 - Isotyping
 - IgG1,2,3,4; IgA; IgM
 - Matrix Metalloproteinases
 - MMP-1, MMP-2, MMP-3 ...
 - Metabolic Markers
 - Collagen Type 1, Collagen Type 4 ...
 - Neurobiology
 - Abeta 40, BDNF, Tau ...
 - Transcription Factors, Nuclear Receptors
 - ATF, EGR, Myc, STAT1, STAT3



RULES • BASED MEDICINE, INC.

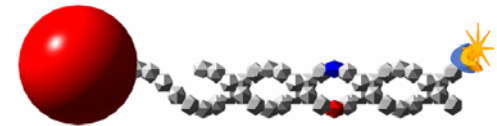
Novagen®



invitrogen™



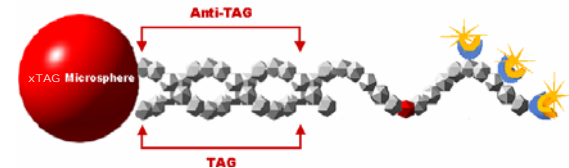
- Gene Expression Profiling
 - GAPDH, PGK1, IL8, VEGF ...
- miRNA Profiling
 - FlexmiR >300 miRNA targets, FlexmiR Select
- Genotyping
 - IL6, ACE, GST, VDR, VKOR ...
 - Y SNPs



Genetic Diseases

Luminex

- Cystic Fibrosis
 - 39 mutations (IVD)
- Cytochrome P450
 - CYP2C19, CYP2D6, CYP2C9
- Ashkenazi Jewish Panel
 - Tay-Sachs
 - Canavan
 - Familial Dysautonomia
 - Gaucher Disease
 - Bloom's Syndrome
 - Fanconi Anemia
 - Niemann-Pick
 - Mucopolipidosis Type IV
- Coagulation
 - Factor V
 - Factor II
 - MTHFR
- Leukemia Translocation Mutations

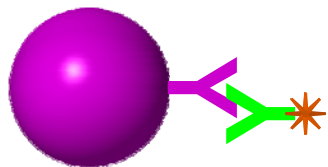


Luminex
MOLECULAR DIAGNOSTICS

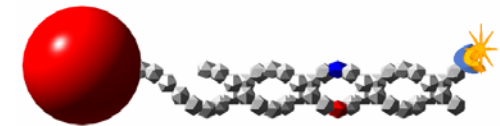
Asuragen



- Allergy Testing
 - ImmuneTech – service
- Autoimmune Disease
 - SSA, SSB, Scl-70, Jo-1 ...
- HLA
 - Class I & Class II antigens
 - HLA-A, HLA-B, HLA-C, DQA, DQB, DRB ...
- Infectious Diseases
 - Pneumococcus
 - EVB
 - Syphilis



- Respiratory Pathogens
 - Respiratory Virus Panel – IVD
 - ResPlex
 - MultiCode PLx
- Staphylococci



- Several Partners provide testing & custom assay services

RULES • BASED MEDICINE, INC.



- Certified Assay Developers



- See Partner Listings for complete menus

<http://www.luminexcorp.com/partners/index.html>

Thank you

Luminex

