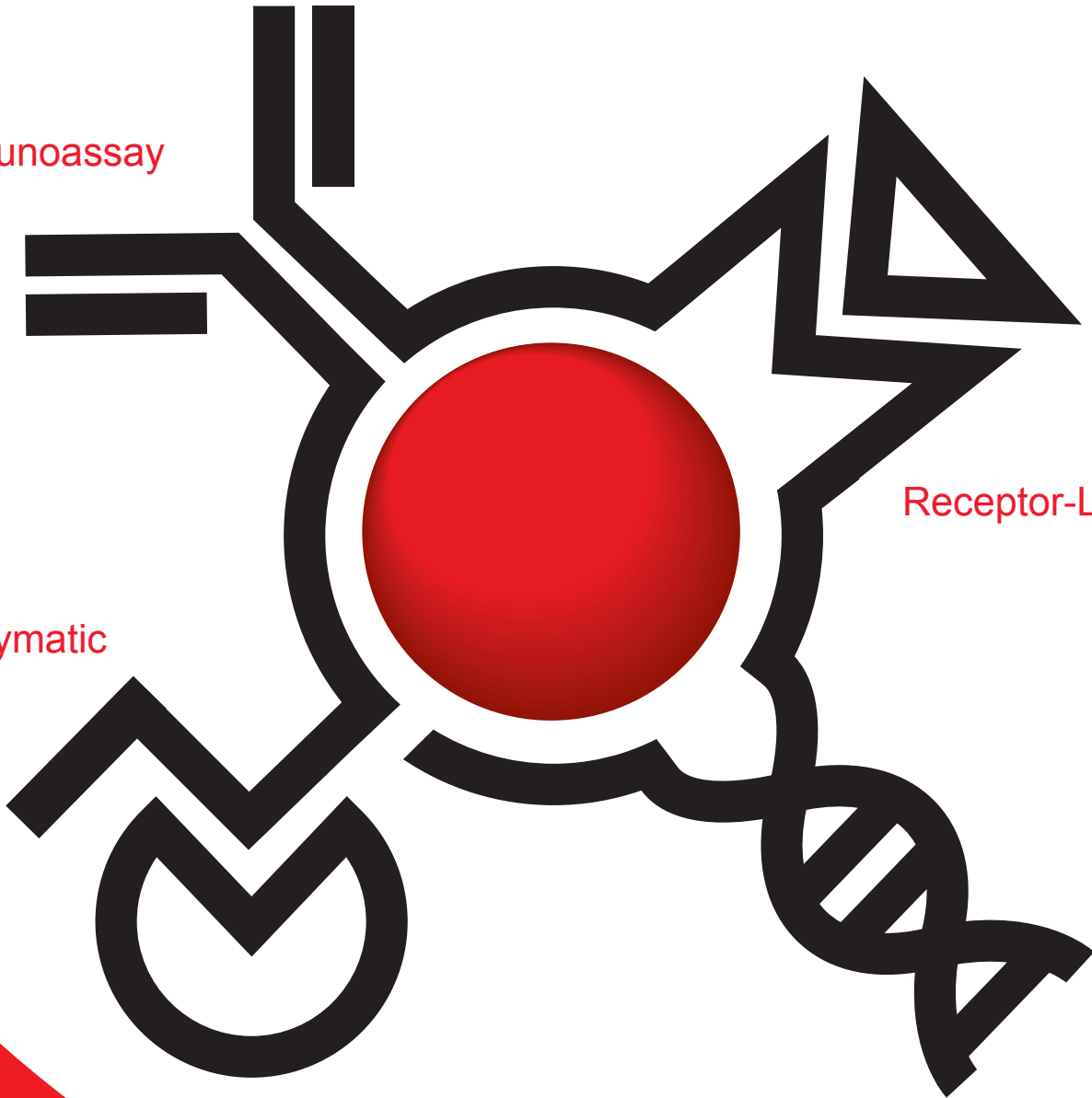


Immunoassay



Receptor-Ligand

Enzymatic

Molecular-Biology

xMAP[®] Technology Analyte Menu



Research Menu

C=Canine, G=General, GP=Guinea Pig, H=Human, M=Mouse, Po=Porcine, Pr=Primate, R=Rat

PROTEIN EXPRESSION PROFILING

Acute Phase Inflammation

beta2 Microglobulin (H)
CRP (H)
Gc-globulin (H)
Haptoglobin (H)
Serum Amyloid A (SAA) (H)

Apoptosis

Active Caspase 3 (H, M)
Akt-PKB (ser473) (H, M)
BAD (H)
c-Jun (ser7) (H, M)
JNK (Thr183/Tyr185) (H)
MIF (H)
PAI-1 (total) (H)
PARP (active, Cleaved) (H)
s Fas (H)
s Fas Ligand (H)
s ICAM-1 (H)
s VCAM-1 (H)

Autoimmune Disease

HSP-27 pS82 (G)
HSP-27 Total (G)

Cancer Markers

alpha Fetoprotein (H)
CA-125 (H)
Cancer Antigen 125 (H)
Carcinoembryonic Antigen (H)
IGF-II (H)
Leptin (H)
MIF (H)
Osteopontin (H)
Prolactin (H)
PSA, Free (H)

Cardiac Markers

Adiponectin (H, M, R)
Apo A1 (M)
Apo E (M)
BNP (R)
Creatine Kinase-MB (H)
CRP (H, M, R)
Endothelin-1 (M)
Fibrinogen (H, M, R)
Haptoglobin (H)
IFN-gamma (H)
IL-10 (H)
IL-1beta (H)
IL-6 (H, R)
IL-8 (H)
MCP-1 (H, R)
MMP-9 (H, M)
MPO (H, R)
NT-proBNP (H)
PAI-1 (total) (H, M, R)
PAP (H)
sE-Selectin (H, M, R)
Serum Amyloid A (H)
Serum Amyloid P (H)
SGOT (H, M)
sICAM-1 (H, M, R)
sIL-2Ralpha (H)
sIL-6Ralpha (H)
sVCAM-1 (H, M)
TIMP-1 (H, M, R)

TNF-alpha (H)
Tnl (R)
TnT (R)
VEGF (H, R)
vWF (R)

Cellular Signaling

3BP2 (G)
ABL2 (G)
Active beta-Catenin (G)
Active Caspase 3 (H, M)
Akt (G)
Akt (Ser473) (G)
Akt (total) (G)
Akt-PKB (ser473) (H, M)
Akt/PKBalpha (Thr308) (H)
ATF2 (Thr71) (G)
ATF2 (total) (G)
B-Raf (pS446) (H, M)
BAD (H)
BTK (G)
c-Jun (Ser63) (G)
c-Jun (ser7) (H, M)
c-Jun (total) (G)
c-Kit (G)
Caspase3 (active) (H)
Caspase8 (active) (H)
CD3epsilon (G)
CRCB (Ser133) (G)
CRCB (total) (G)
CREB (Ser 133) (G)
CRK (G)
CRKL (G)
CSK (G)
DAPP1 (G)
EGF-Receptor (G)
EGFR (Tyr) (G)
Erk 1/2 (Total) (G)
Erk-2 (G)
Erk/MAPK 1/2 (Thr185/Tyr187) (G)
ERK1 (pT202/pY204) (H, M)
Erk1 (Thr202/Tyr204) (G)
ERK1 Total (H, M)
Erk1/2 (Thr202/Tyr204, Thr185/Tyr187) (G)
ERK2 (pT185/pY187) (H, M)
Erk2 (Thr185/Tyr187) (G)
Erk2 (total) (G)
ERK2 Total (H, M)
FYN (G)
GAPDH (total) (H)
GRAP (G)
GRB10 (G)
GRB14 (G)
GRB2 (G)
GSK-3a/b (Ser21/Ser9) (G)
GSK-3beta (G)
GSK3 (H)
Histone H3 (Ser10) (G)
HSP-27 (Ser78) (G)
HSP-27 Total (G)
IkappaB-alpha (phospho) (H)
IkappaB-alpha (Ser32/Ser36) (G)
IkappaB-alpha Total (G)
IkappaBalpha (serS32) (G)
IRS-1 (Ser636/Ser639) (G)
IRS1 (G)
JNK (pTpy183/185) (G)
JNK (Thr183/Tyr185) (H)
JNK Total (G)
JNKP (Thr183/Y182) (G)
LAT (G)
LCK (G)
lck (G)

LCP2 (G)
MAPKAP K2 (G)
MATK (G)
MEK1 (Ser217/Ser221) (G)
MEK1 (total) (G)
MEK1 Total (H, M)
MEK1/2 (pS222/pS226) (H, M)
MEK2 Total (H, M)
NF-kappaB p65 (Ser536) (G)
NFkB (H)
NSP1 (G)
p38 (Thr180/Tyr182) (G)
p38 MAPK (Thr150/Tyr182) (G)
p38 MAPK (total) (G)
p38 MAPK pT180/pY182 (G)
p53 pS15 (G)
P55G-D1 (G)
p70 S6 Kinase (Thr421) (G)
p70 S6 kinase (Thr421/Ser424) (G)
P85A-D1 (G)
P85A-D2 (G)
P85B-D1 (G)
P85B-D2 (G)
p90RSK (Thr359/Ser363) (G)
p90RSK (total) (G)
PARP (active, Cleaved) (H)
PARP (cleaved) (H)
PARP (total) (H)
PDGF Receptor-alpha (G)
PDGF Receptor-beta (G)
PDGF receptor-beta (Tyr751) (G)
Phospho PKC alpha (G)
PKB-alpha (G)
PKC (G)
PLCG1-D1 (G)
PTPN11-D2 (G)
PTPN6-D2 (G)
pTYR EGFR (H)
pTYR HER-2 (H)
pTYR HGFR (H)
pTYR IGF-1R (H)
pTYR PDGFRbeta (H)
pTYR Tie-2 (H)
pTYR VEGFR2 (H)
Raf1 (pS338) (H, M)
Raf1 Total (H, M)
Rsk1 (Ser380) (G)
S6 Ribosomal Protein (Ser235/236) (G)
SAPK1 (G)
SAPK1a/JNK2 (G)
SAPK4 (G)
SHC1 (G)
SOCS2 (G)
STAP2 (G)
STAT1 Total (H, M)
STAT2 (Tyr689) (G)
STAT3 (Ser727) (G)
STAT3 (Tyr705) (G)
STAT5A/B (Tyr694/699) (G)
STAT6 (Tyr641) (G)
Syk (G)
SYK-D2 (G)
TNS (G)
Total EGFR (H)
Total HER-2 (H)
Total HGFR (H)
Total IGF-1R (H)
Total PDGFRbeta (H)
Total VEGFR2 (H)
TrkA (Tyr490) (G)
VAV3 (G)
ZAP-70 (G)

Cytokines, Chemokines & Growth Factors

Abeta 40 (H)
Abeta 42 (H)
Adiponectin (H)
BDNF (H)
beta-2 Microglobulin (H)
beta-NGF (H)
C-Reactive Protein (H)
CCL11/Eotaxin (H)
CCL2/MCP-1 (H)
CCL3/MIP-1 alpha (H)
CCL4/MIP-1 beta (H)
CCL5/RANTES (H)
CD40 Ligand/TNFSF5 (H)
CINC-2 alpha/beta (R)
CINC-3 (R)
CRP (H)
CTACK (H)
CXCL10/IP-10/CRG-2 (H)
CXCL11/I-TAC (H)
CXCL5/ENA (H)
CXCL8/IL-8 (H)
DR5 (H)
EGF (H, M)
ENA-78 (H)
Eotaxin (H, M, R)
Eotaxin/CCL11 (H)
Fatty Acid Binding Protein (H)
FGF-2 (H, M)
FGF-basic (H, M)
FLT-3 ligand (H)
Fractalkine (H)
G-CSF (H, M, Pr, R)
Gc-globulin (H)
GCP-2 (M)
GDNF (H)
GM-CSF (C, H, M, Pr, R)
GM-CSF Ultrasensitive (H)
GRO (H)
GRO alpha (H)
GRO-KC (R)
Haptoglobin (H)
Hepatocyte GF (H)
HGF (H, M)
I-CAM (R)
I-TAC (H)
ICAM-1 (H)
IFN-alpha (H)
IFN-alpha2 (H)
IFN-gamma (C, H, M, Pr, R)
IFN-gamma Ultrasensitive (H)
IL-10 (C, H, M, Pr, R)
IL-10 Ultrasensitive (H)
IL-11 (M)
IL-12 (P, R)
IL-12 p40 (H, M)
IL-12 p40/p70 (H, M, R)
IL-12 p70 (H, M, R)
IL-12/23 p40 (Pr)
IL-13 (H, M, Pr, R)
IL-15 (C, H, M, Pr)
IL-16 (H)
IL-17 (H, M, Pr, R)
IL-18 (C, H, M, Pr, R)
IL-18/IL-1F4 (R)
IL-1a (H, Pr)
IL-1alpha (H, M, R)
IL-1alpha/IL-1F1 (H, R)
IL-1alpha/IL-1F2 (H, M, R)
IL-1beta (H, M, Pr, R)
IL-1beta Ultrasensitive (H)
IL-1ra (H)
IL-1ra/IL-1F3 (H)

IL-2 (C, H, M, Pr, R)
IL-2 Ultrasensitive (H)
IL-2R (H)
IL-2Ralpha (H)
IL-3 (H, M)
IL-4 (C, H, M, Pr, R)
IL-4 Ultrasensitive (H)
IL-5 (H, M, Pr, R)
IL-5 Ultrasensitive (H)
IL-6 (C, H, M, Pr, R)
IL-6 Ultrasensitive (H)
IL-6R (H)
IL-7 (C, H, M)
IL-8 (C, H, Pr)
IL-8 Ultrasensitive (H)
IL-9 (H, M, R)
Insulin (H)
IP-10 (C, H, M, R)
JE/MCP-1 (M)
KC (C, M, R)
KC/GR0a (M)
L-Selectin/CD62L (R)
Leptin (H, R)
LIF (H, M)
LIX (M)
Lymphotacin (H, M)
M-CSF (H, M)
MCP-1 (C, H, M, Pr, R)
MCP-1 (MCAF) (M)
MCP-1(MCAF) (H)
MCP-1/CCL2 (H, M, Pr)
MCP-2 (H)
MCP-2/CCL8 (H)
MCP-3 (H, M)
MCP-3/CCL7 (H)
MCP-5 (M)
MDC (H, M)
MDC/CCL22 (H)
MIF (H)
MIG (H, M)
MIG/CXCL9 (H, M)
MIP-1 alpha (H, M, Pr)
MIP-1 beta (H, M, Pr)
MIP-1 gamma (M)
MIP-1-alpha (R)
MIP-1alpha/CCL3 (H, M, Pr)
MIP-1beta/CCL4 (H, M, Pr)
MIP-2 (M)
MIP-3 beta (M)
MIP-3beta/CCL19 (M)
MMP-13 (H)
MMP-3 (H)
MMP-9 (H)
Neurobiology buffer kit (H)
NGF (H)
OSM (M)
PDGF-AA (H)
PDGF-AB/BB (H)
PDGF-BB (H, M)
RANTES (H, M, R)
RANTES/CCL5 (H, M, Pr)
Rb (pT821) (H)
Resistin (H)
sCD30 (H, M)
sCD40L (H)
SCF (H)
SCGF-beta (H)
SDF-1alpha (H)
sE-Selectin/CD62E (H)
sEGFR (H, M)
Serum Amyloid A (SAA) (H)
sFAS-ligand (H)
sgp130 (H, M)
sICAM-1/CD54 (H, R)
sIL-1RI (H, M)

sIL-1RII (H, M)
 sIL-2Ralpha (H, M)
 sIL-4R (H, M)
 sIL-6R (H, M)
 sP-Selectin/CD62P (H)
 sRAGE (H, M)
 sTNFRI (H, M)
 sTNFRII (H, M)
 sVCAM-1/CD106 (H)
 sVEGFR1 (H, M)
 sVEGFR2 (H, M)
 sVEGFR3 (H, M)
 Tau (pS214) (H)
 Tau (pS396) (H)
 TGF alpha (H, Pr)
 TGF-beta1 (activated/treated) (H, M, Pr, R)
 TGFbeta1 (H, M, Po, R)
 Thrombopoietin (Tpo) (H)
 TIMP-1 (R)
 Tissue Factor (H, M)
 TNF-alpha (C, GP, H, M, Pr, R)
 TNF-alpha Ultrasensitive (H)
 TNF-alpha/TNFSF1A (H, M, R)
 TNF-beta (H)
 TNF-RI (H)
 TNF-RII (H)
 TNFSF5 (H)
 total Tau (H)
 TRAIL (H)
 V-CAM (R)
 VCAM-1 (H, M)
 VEGF (H, M, Pr, R)

Endocrine

ACTH (H, R)
 Adiponectin (H, M)
 Adiponectin/Acrp30 (H)
 Amylin (active) (H, M, R)
 Amylin (total) (H, M, R)
 BDNF (R)
 C-Peptide (H)
 C-Reactive Protein (H)
 Calcitonin (H)
 Complement Factor D (H)
 Corticosterone (R)
 CRF (H)
 FGF-9 (M)
 FSH (H, R)
 GH (H, R)
 Ghrelin (active) (H)
 Ghrelin (total) (M, R)
 GIP (total) (H, M, R)
 GLP-1 (C, H, M, R)
 GLP-1 (active) (H, M, R)
 Glucagon (C, H, M, R)
 Growth Hormone (H, M)
 IGF-1 (M, R)
 Insulin (C, H, M, R)
 Leptin (C, H, M, R)
 LH (H, R)
 Lipoprotein (a) (H)
 Melatonin (R)

PAI-1 (active) (H)
 PAI-1 (total) (H, M)
 Pancreatic Polypeptide (H, M, R)
 Prolactin (H, R)
 PYY (total) (H, M, R)
 Resistin (H, M, R)
 Serpin E1/PAI-1 (H)
 T3 (H, R)
 T4 (H, R)
 TBG (H)
 Thyroglobulin (H)
 TSH (H, R)

Isotyping

IgA (H, M)
 IgA1 (H)
 IgA2 (H)
 IgE (H, M)
 IgG (H, M)
 IgG1 (H, M)
 IgG2 (H)
 IgG2alpha (M)
 IgG2beta (M)
 IgG3 (H, M)
 IgG4 (H)
 IgM (H, M)
 light chain (kappa or gamma) (M)

Matrix Metalloproteinases

MMP (H)
 MMP-1 (H)
 MMP-12 (H)
 MMP-13 (H)
 MMP-2 (H)
 MMP-3 (H)
 MMP-7 (H)
 MMP-8 (H)
 MMP-9 (H)

Metabolic Markers

Apolipoprotein A-1 (M)
 Apolipoprotein A-I (H)
 Apolipoprotein A-II (H)
 Apolipoprotein B (H)
 Apolipoprotein C-II (H)
 Apolipoprotein C-III (H)
 Apolipoprotein E (H)
 beta-2 Glycoprotein (H, M)
 Collagen Type 1 (H)
 Collagen Type 2 (H)
 Collagen Type 4 (H)
 Collagen Type 6 (H)
 Apolipoprotein E (H)
 beta-2 Glycoprotein (H, M)
 Collagen Type 1 (H)
 Collagen Type 2 (H)
 Collagen Type 4 (H)
 Collagen Type 6 (H)
 Glutathione S-Transferase (H, M)
 Pancreatic Islet Cells (H)
 tTG (Celiac Disease) (H)

Neurobiology

Aggregated A β (H)
 Aggregated α -Synuclein (H)
 A α 40 (H)
 A α 42 (H)

BDNF (H)
 GDNF (H)
 PDGF-BB (H)
 PDGF-BB (M)
 Tau (total) (H)
 Tau [pS199] (H)
 Tau [pT181] (H)
 IL-8 (H)
 IL-6 (H)
 IL-1 α (H)
 TNF- α (H)

Other

ACTH (M, R)
 Adiponectin (H)
 Cortisol (H)
 Fibronectin (H)
 IL-1beta (M)
 IL-6 (M)
 Insulin (M, R)
 Involucrin (H)
 Keratin-1, 10, 11 (H)
 Keratin-6 (H)
 Leptin (M, R)
 OPG (M, R)
 OPN (H)
 Osteocalcin (M, R)
 PTH (H)
 RANKL (H, M, R)
 TNF-alpha (M)

Signal Transduction

Akt (pS473) (G)
 Akt (total) (G)
 c-Kit (pY823) (G)
 c-Kit (total) (G)
 c-Met (pY1349) (G)
 c-Met (total) (G)
 Caspase-3 (175/176) (G)
 CREB (pS133) (G)
 CREB (Total) (G)
 Cytochrome-c (G)
 EGFR (pY1068) (G)
 EGFR (total) (G)
 Erk 1/2 (pTpY185/187) (G)
 Erk 1/2 (Total) (G)
 GS-1 (pSpS641/645) (G)
 GS-1 (Total) (G)
 GSK-3beta (pS9) (G)
 GSK-3beta (total) (G)
 HSP27 (pS82) (G)
 HSP27 (total) (G)
 IGF-1R (pYpY1135/1136) (G)
 IGF-1R (total) (G)
 I κ Balpha (pS32) (G)
 I κ Balpha (total) (G)
 IR (pYpY1162/1163) (G)
 IR (total) (G)
 IRS-1 (pS312) (G)
 JNK 1/2 (pTpY183/185) (G)
 JNK 1/2 (total) (G)
 p38 MAPK (pTpY180/182) (G)
 p38 MAPK (total) (G)

p53 (pS15) (G)
 p53 (total) (G)
 p70S6K (pTpS421/424) (G)
 p70S6K (total) (G)
 PARP Cleaved (G)
 PIM-1 (pS8) (G)
 PIM-1 (total) (G)
 PRAS40 (pT246) (G)
 PRAS40 (total) (G)
 Retinoblastoma (Rb) (pSpT249/252) (G)
 Retinoblastoma (Rb) (pT821) (G)
 Retinoblastoma (Rb) (total) (G)
 STAT1 (pY701) (G)
 STAT1 (total) (G)
 STAT3 (pY705) (G)
 STAT5a/b (pY694) (G)
 Tau (pS199) (G)
 Tau (pT181) (G)
 Tau (total) (G)
 VEGFR2 (pY1059) (G)

Transcription Factors/Nuclear Receptors

AP-1 (G)
 AP-2 (G)
 AR (G)
 ATF2 (G)
 BRN3 (G)
 C-MYB (G)
 CEBP (G)
 CRE-ATF1 (G)
 CREB (G)
 E2F1 (G)
 E2F1-5 (G)
 E2F6 (G)
 EGR (G, H)
 ELK1 (G)
 ER (G)
 ETS/PEA (G)
 FAST1 (G)
 FKHR-1 (G)
 GATA (G)
 GR/PR (G)
 HIF-1 (H)
 HNF1 (G)
 HNF4 (G)
 IRF1 (G)
 ISRE (G)
 MEF-2 (G)
 Myc-Max (G)
 MYOD (G)
 NF-1 (G, H)
 NF-E1(Y1) (G)
 NF-E2 (G)
 NF-Y (G)
 NFAT (H)
 NF κ B (G)
 NF κ B Gene Family (H)
 NKX-2.5 (G)
 OCT (G)
 p53 (G)
 PAX3 (G)

PAX5 (G)
 PPAR (G, H)
 RUNX/AML (G)
 SMAD (G)
 SMAD 2/3 (G)
 SP1 (G)
 SRE (H)
 STAT1 (G)
 STAT3 (G)
 STAT4 (G)
 STAT5 (G)
 YY1 (G, H)

GENOMIC RESEARCH

Gene Expression Profiling

1L6R (H)
 ACTB (H)
 BAD (H)
 BAK1 (BAK) (H)
 BCL2 (H)
 BCL2L1 (BCL-XL) (H)
 CDKN1A (CDKN1) (H)
 CFLAR (CFLIP) (H)
 CSF2 (H)
 Customer Defined Assays (C, G, GP, H, M, Po, Pr, R)
 IFN-gamma (H)
 IL-10 (H)
 IL-2 (H)
 NFKB2 (H)
 NFKBIA (NFKIA) (H)
 NKFB1 (H)
 PPIB (H)
 Plk2B (RAFTK) (H)
 RELA (H)
 RELB (H)
 TNF (H)
 TNFAIP3 (A20) (H)
 TNFRSF6 (FAS) (H)
 TNFSF6 (FASL) (H)
 VEGF (H)

Genotyping

Customer Defined Assays (C, G, GP, H, M, Po, Pr, R)
 Mitochondrial DNA Screening (H)
 Y-SNP Identification (H)

MicroRNA

Customer Defined Assays (C, G, GP, H, M, Po, Pr, R)
 Equivalent to Sanger miRBase (H, M, R)

Diagnostic Menu

AUTOIMMUNE TESTING

ANA Panel® (Bio-Rad)

- dsDNA, quantitative
- Chromatin, semi-quantitative
- Ribosomal Protein
- SSA
- SSB
- Sm
- Sm/RNP
- RNP
- Scl-70
- Jo-1
- Centromere b

ANA Panel* (Zeus/Inverness)

- SSA
- SSB
- Sm
- RNP
- Scl-70
- Jo-1
- Centromere b
- Histone
- IgG class antibody to dsDNA, quantitative
- IgG class antibody to ANA, qualitative

ANCA Panel* (Zeus/Inverness, INOVA)

- Myeloperoxidase, IgG class antibody, qualitative and semi-quantitative (Zeus/Inverness)
- Proteinase 3, IgG class antibody, qualitative and semi-quantitative (Zeus/Inverness)
- Myeloperoxidase (INOVA)
- Proteinase 3 (INOVA)

Celiac IgA Panel* (INOVA)

- IgA (htTG), semi-quantitative
- Anti-DGP

Celiac (htTG) Connective Panel* (BMD)

- DS dna
- SSA 60 kDa
- SSA 52 kDa
- SSB
- Sm
- Sm/RNP
- Scl70
- Jo-1
- Ribosome
- Centromere

ENA Panel* (INOVA)

- Sm
- RNP
- SSA
- SSB
- Scl-70
- Jo-1

Rheumatoid Factor Panel* (Zeus/Inverness)

- RF IgM, qualitative or quantitative

SLE Profile* (INOVA)

- Sm
- RNP
- SSA
- SSB
- Scl-70
- Jo-1
- Ribosome-P
- Chromatin

Thyroid Panel* (BMD, Zeus/Inverness)

- TPO IgG, qualitative and semi-quantitative (Zeus/Inverness)
- Tg, IgG, qualitative and semi-quantitative (Zeus/Inverness)
- TPO, semi-quantitative (BMD)
- Tg, semi-quantitative (BMD)

Vasculitis Panel* (BMD)

- ANCA
- Anti-MPO
- Anti-PR3
- GBM

HUMAN GENETICS TESTING

Ashkenazi Jewish Panel (Asuragen, Fisher HealthCare)

- Tay-Sachs Syndrome
- Canavan Syndrome
- Gaucher Syndrome
- Bloom Syndrome
- Fanconi Anemia (Group C)
- Familial Disautonomia Syndrome
- Niemann-Pick Syndrome
- Mucopolidosis Type IV

Coagulation Panel (Fisher HealthCare)

- Factor II
- Factor V
- MTHFR

Cystic Fibrosis 39 IVD Panel* (Fisher HealthCare)

Cystic Fibrosis 70 Panel (Fisher HealthCare)

Leukemia Translocations (Asuragen)

Pharmacogenetics (Fisher HealthCare)

- CYP450 2C9
- CYP450 2C9 + VKORC1
- CYP450 2C19
- CYP450 2D6

HLA TESTING* (One Lambda, Tepnel Lifecodes)

- HLA Class I and II Human
- HLA Class I Single Antigen Antibody, Group 1 Human
- HLA Class I Single Antigen Antibody, Group 2 Human
- PRA Class I Human
- PRA Class I and II Human
- PRA Class II Human
- SSO Class I HLA-A Human
- SSO Class I HLA-B Human
- SSO Class I HLA-C Human
- SSO Class II DP Human
- SSO Class II DQB1 Human
- SSO Class II DRB1 Human

INFECTIOUS DISEASE TESTING

Epstein-Barr Virus

EBV IgM Panel* (Bio-Rad)

- EBV VCA antibodies
- Heterophile antibodies in human serum

EBV IgG Panel* (Bio-Rad)

- EBV-NA-1
- EBV VCA
- EBV EA-D

EBV Panel* (Zeus/Inverness)

- EBV VCA IgM, qualitative
- EBV VCA IgG, qualitative
- EBV-VCA gp-125
- Total EBV early antigens
- Recombinant EBV nuc anti-1

Herpes Panel* (Focus Diagnostics)

- HSV 1 IgG, qualitative
- HSV 2 IgG, qualitative

MMRV Panel* (Zeus/Inverness)

- Measles, IgM, qualitative
- Mumps, IgM, qualitative
- Rubella, IgG, qualitative
- Varicella IgG, qualitative

Syphilis Panel* (Bio-Rad)

- IgG, qualitative

• HSV Panel (Zeus/Inverness)

MOLECULAR INFECTIOUS DISEASE

Respiratory Viral Panel* (Fisher Healthcare/Abbott Molecular)

ResPlex I (QIAGEN)

ResPlex II (QIAGEN)

StaphPlex (QIAGEN)

NEURODEGENERATIVE DISEASE

Alzheimer's Panel (Innogenetics)

- Aβ1-42
- Total Tau
- P-Tau181P

*510(k) cleared

**Pending FDA clearance

UNITED STATES

Luminex Corporation
Austin, Texas
Tel: +1.512.219.8020
Fax: +1.512.219.5195

CANADA

Luminex Molecular Diagnostics
Toronto, Ontario
Tel: +1.416.593.4323
Fax: +1.416.593.1066

EUROPE

Luminex B.V.
Oosterhout, The Netherlands
Tel: +31.162.408333
Fax: +31.162.408337

CHINA

Luminex Shanghai Trading Co.
Shanghai, China
Tel: +21.616.50809
Fax: +21.616.50811

JAPAN

Luminex Japan Corporation, Ltd.
Tokyo, Japan
Tel: +81.3.5545.7440
Fax: +81.3.5545.0451

AUSTRALIA

BSD Robotics, A Luminex Company
Brisbane, Queensland
Tel: +61.7.3273.0273
Fax: +61.7.3273.0274

www.luminexcorp.com

info@luminexcorp.com