

Simple Plex Assay Menu



Impurity Testing

CHO-HCP 3G-1
Endonuclease
Green Fluorescent Protein (GFP)
HEK 293 HCP 3G
Protein A

Viral Titer Quantification

AAV1
AAV2
AAV6
AAV8

Human Analytes

Adiponectin/Acrp30, total
AFP
Alpha 2-Macroglobulin
alpha-Synuclein
Amphiregulin
Amyloid Beta aa1-40
Amyloid Beta aa1-42
Angiogenin
Angiopoietin-1
Angiopoietin-2
Angiopoietin-like 3 (ANGPTL3)
Angiopoietin-like 4 (ANGPTL4)
ASC
Axl
BAFF
BCMA/TNFRSF17
BDNF, free

beta-Synuclein
BLC/BCA-1/CXCL13
BMP-2
BMP-9
C-Reactive Protein (CRP)
CA125
CA9
Caspase-1
CCL16/HCC-4
CCL17/TARC
CCL18/PARC
CCL19/MIP-3 beta
CCL2/MCP-1
CCL3/MIP-1 alpha
CCL5/RANTES
CD14
CD21



Automated ELISA For Single and Multiplex Detection

Scan the QR Code or Visit:
bio-techne.com/simple-plex



Human Analytes

CD23/Fc epsilon RII *Discovery*

CD25/IL-2 R alpha

CD27

CD28 *Discovery*

CD40 Ligand

CD40/TNFRSF5

CD155/PVR *Discovery*

CD163

CG beta *Discovery*

Chitinase 3-like 1/YKL-40

Clusterin

COMP/Thrombospondin-5

Coagulation Factor III / Tissue Factor *Discovery*

CTGF/CCN2

CTLA-4

CXCL1

CX3CL1/Fractalkine

CXCL9/MIG

CXCL10/IP-10

CXCL12/SDF-1 alpha

Cyr61 (CCN1)

Cystatin C

Cytokeratin 18

D-Dimer

DcR3/TNFSF6B

DKK-1

DLL1 *Discovery*

DPPIV/CD26 *Discovery*

E-Cadherin

E-Selectin

EGF

ENA-78/CXCL5

Endoglin/CD105

Endothelin-1

ENPP-2

Eotaxin-3/CCL26

EPO

ErbB2

FABP2/I-FABP

FAP

Fas/TNFRSF6

Fas Ligand/TNFSF6

Ferritin

FGF-19

FGF-21

FGF-23

FGF-7/KGF

Fibronectin

Flt-3 Ligand

Follistatin

Fractalkine/CX3CL1

G-CSF

Galectin-1 *Discovery*

Galectin-3

Galectin-3BP/MAC-2BP *Discovery*

Galectin-9 *Discovery*

GAS6

GDF-8

GDF-15

GDNF *Discovery*

GFAP

GM-CSF

gp130

Granzyme A

Granzyme B

Growth Hormone (hGH)

HE4/WFDC2

Hepcidin

HGF

HGF R/c-MET

HIV-1 Gag p24

HIV-1 p24

HVEM

ICAM-1

IFN-alpha (multi-subtype)

IFN-alpha 2

IFN-beta

IFN-gamma

IGFBP-1

IGFBP-2

IGFBP-rp1/IGFBP-7

IL-1-alpha

Human Analytes

IL-1-beta
IL-1ra/IL-1F3
IL-2
IL-3 *Discovery*
IL-4
IL-5
IL-6
IL-6 R-alpha
IL-7
IL-8/CXCL8
IL-10
IL-11
IL-12-p70
IL-13
IL-15
IL-16
IL-17A
IL-18
IL-18 BPa
IL-19
IL-22
IL-33
I-TAC
Insulin
LAG-3
LBP
LDLR *Discovery*
Leptin
LIGHT
Lipocalin-2/NGAL
M-CSF
MAdCAM-1
MBL *Discovery*
MCP-2/CCL8
MCP-3/CCL7
MDC/CCL22
Mesothelin
MICA
MIF
MIP-1 alpha/CCL3
MIP-1 beta/CCL4
MIP-3 alpha/CCL20
MMP-7
MMP-8
MMP-9
MMP-9, Total *Discovery*
Myeloperoxidase (MPO)
Nectin-2
Nectin-4 *Discovery*
Neprilysin/CD10
Neurofilament Light (NF-L)
Neurofilament Heavy (NF-H) *Discovery*
Neurogranin
Neutrophil Elastase/ELA2
NGF R/TNFRSF16
NGF-beta
Osteopontin (OPN)
Osteoprotegrin/OPG
OX40/TNFRSF4 *Discovery*
Pancreatic Polypeptide/PP *Discovery*
Park 7 *Discovery*
PCSK9/PC9
PD-1 *Discovery*
PD-L1/B7-H1
PD-L2 *Discovery*
PDGF-AB *Discovery*
PDGF-BB
Pentraxin 3/TSG-14
Perforin
Periostin/OSF-2
Placenta Growth Factor (PIGF)
Pro-Gastrin-releasing Peptide
Procalcitonin
P-Selectin/CD62P
RAGE
Reg-3a
Relaxin-2 *Discovery*
Resistin
SCF
Serpins A1
Serpins A4/Kallistatin
Serpins E1/PAI-1

Human Analytes

Siglec-9
SP-D
Spinesin *Discovery*
ST2/IL-1 R4 (IL-33 R)
Syndecan-4
S100B
TACI/TNFRSF13B
Tau, Phospho 217
Tau, Total
TFF3
TFPI *Discovery*
TfR *Discovery*
TGF-alpha *Discovery*
TGF-beta 1
Thrombomodulin *Discovery*
Thrombospondin-1 (THBS1)
Thrombospondin-2 *Discovery*
Tie-2
TIM-1/KIM-1/HAVCR
TIM-3
TIMP-1
TIMP-2
TNF alpha
TNF R1
TNF RII
TRAIL
Trappin-2/Elafin
TREM-1
TREM-2
TSLP *Discovery*
UCH-L1
uPAR
u-Plasminogen/uPA-1
Uteroglobin *Discovery*
VAP-1 *Discovery*
VCAM-1
VEGF (VEGF-A)
VEGF R1/Flt-1
VEGF R2/KDR
VEGF-B
VEGF-C
Vitamin D

Rat Analytes

Angiopoietin-2
beta-Synuclein
CD25/IL-2 R alpha
Chitinase 3-like 1
Cystatin C
FGF-21
Flt-3 Ligand
GDF-8
IL-2 R alpha
Lipocalin 2
MCP-1/CCL2
M-CSF
MMP-9, Total *Discovery*
Neurofilament Light (NF-L)
Osteopontin (OPN)
Park 7 *Discovery*
RANTES/CCL5
SP-D
S100B
TGF-beta 1
TIM-1
TREM-2
VCAM-1/CD106
VEGF-A
VEGF-D

Mouse Analytes

Adiponectin
AFP
Angiopoietin-2
Angiopoietin-like 3 (ANGPTL3)
BAFF
beta-Synuclein
CCL3/MIP-alpha
CCL4
CD25/IL-2 R alpha
Chitinase 3-like 1
C-Reactive Protein/CRP
CXCL1/KC
CXCL15
CXCL2/MIP-2
CXCL13/BLC/BCA-1
Cystatin C
Erythropoietin/EPO
FGF-21
Flt-3 Ligand
G-CSF
GDF-8
GDF-15
GM-CSF
Granzyme B
IFN-gamma
IL-1-beta
IL-17
IL-1ra/IL-1F3
IL-2
IL-2 R alpha
IL-4
IL-5
IL-6
IL-7
IP-10/CXCL10
IL-12 p70
Lipocalin-2/NGAL
MCP-1/CCL2
M-CSF
Neurofilament Light (NF-L)
Osteopontin (OPN)
P-Selectin
Park 7 *Discovery*
RANTES/CCL5
SDF-1 alpha/CXCL12
SP-D
ST2/IL-33R/IL-1R4
S100B
TGF-beta 1
TIM-1/KIM-1
TNF-alpha
TREM-2
VCAM-1/CD106
VEGF-A

Non-Human Primate Analytes

Angiopoietin-1

Angiopoietin-like 3 (ANGPTL3)

BDNF, free

CCL19/MIP-3 β

CCL2/MCP-1

CCL3/MIP-1 α

CCL4/MIP-1 β

CCL5/RANTES

CD163

C-Reactive Protein (CRP)

CTLA-4

CXCL10/IP-10

CXCL13/BLC/BCA-1

FAS Ligand

Fibronectin

FLT-3 Ligand

GM-CSF

Granzyme A

Granzyme B

IFN-alpha

IFN-gamma

IGFBP-rp1/IGFBP-7

IL-1b

IL-2

IL-4

IL-5

IL-6

IL-6 R α

IL-8/CXCL8

IL-13

IL-15

IL-17A

IL-18

Insulin

Nectin-2/CD112

Neurofilament Light (NF-L)

PDGF-BB

PD-L1/B7-H1

Perforin

P-Selectin/CD62P

TIMP-2

TNF R1

TNF RII

TNF α

uPAR



**Automated ELISA For
Single and Multiplex Detection**

Scan the QR Code or Visit:
bio-techne.com/simple-plex

Contact Us

Global info@bio-techne.com, bio-techne.com/find-us/distributors

North America TEL 800 343 7475

Europe // Middle East // Africa TEL +44 (0)1235 529449

China info.cn@bio-techne.com, TEL 400.821.3475

For research use or manufacturing purposes only. Trademarks and registered trademarks are the property of their respective owners.
8747214299_0425

bio-techne[®]

Global Developer, Manufacturer, and Supplier of High-Quality Reagents,
Analytical Instruments, and Precision Diagnostics.

INCLUDES R&D Systems™ Novus Biologicals™ Tocris Bioscience™ ProteinSimple™ ACD™ ExosomeDx™ Asuragen™ Lunaphore™