

# Quanterix SP-X™ Multiplex Biomarker Assay Menu

Unleash the power of next-generation Simoa® Planar Array technology for robust multiplex circulating biomarker detection at the earliest stages of disease progression - even at healthy baseline levels.

With a variety of species, multiplex and single plex options as well as our HomeBrew offerings, there is a wide array of options for your experiments. Also offered is our Custom assay design with our Accelerator Laboratory.

### Simoa CorPlex

Fully validated, reliable multiplex and single plex assays. Included are standard curve ranges, Limit of Detection (LOD), Lower Limit of Quantification (LLOQ), precision, spike and recovery, along with additional levels of validation.

### Simoa HomeBrew

Homebrew Assay Starter Kit has been configured with flexibility in mind, providing materials sufficient for making up to 25 plates worth of reagents to design your own assay with your antibody pair selection.

### Simoa Multi-Plex Assay

Available in Human, Mouse, Rat, and NHP options. Allows the ability to multiplex within a single assay. Subplexes of the established multiplexes are also available.

### Simoa Developer Single & Multi – Plex

Removes the time-consuming and expensive step of antibody pair screening because it has been done for you. Assays are available for Human, Mouse, Rat, NHP, and Porcine.

## SIMOA CORPlex PLANAR ARRAY ASSAYS - HUMAN

Assay	Analytes	LoD (pg/mL)	LoQ (pg/mL)	Dynamic Range (pg/mL)	Median Endogenous (pg/mL)	Sample Volume	Sample Type	Catalog Number
<b>CorPlex Cytokine Panel 1 (10-Plex)</b>	IFN- $\gamma$	0.011	0.195	0-200	0.77	12.5 $\mu$ L	E, S	85-0329
	IL-1 $\beta$	0.024	0.195	0-800	0.19			
	IL-4	0.014	0.098	0-400	0.08			
	IL-5	0.012	0.391	0-800	0.46			
	IL-6	0.038	0.445	0-456	5.18			
	IL-8	0.364	3.125	0-1,600	32.25			
	IL-10	0.032	0.195	0-800	2.11			
	IL-12p70	0.044	0.293	0-2,400	1.45			
	IL-22	0.01	0.156	0-400	1.66			
TNF- $\alpha$	0.055	1.563	0-1,600	0.93				
Assay	Analyte	LoD (pg/mL)	LoQ (pg/mL)	Dynamic Range (pg/mL)	Median Endogenous (pg/mL)	Sample Volume	Sample Type	Catalog Number
<b>CorPlex IFN-<math>\gamma</math></b>	IFN- $\gamma$	0.011	0.195	0-200	0.77	12.5 $\mu$ l	E,S	85-0415
<b>CorPlex IL-1<math>\beta</math></b>	IL-1 $\beta$	0.024	0.195	0-800	0.19			85-0348
<b>CorPlex IL-4</b>	IL-4	0.014	0.098	0-400	0.08			85-0416
<b>CorPlex IL-5</b>	IL-5	0.012	0.391	0-800	0.46			85-0406
<b>CorPlex IL-6</b>	IL-6	0.038	0.445	0-456	5.18			85-0341
<b>CorPlex IL-8</b>	IL-8	0.364	3.125	0-1,600	32.25			85-0417
<b>CorPlex IL-10</b>	IL-10	0.032	0.195	0-800	2.11			85-0418
<b>CorPlex IL-12p70</b>	IL-12p70	0.044	0.293	0-2,400	1.45			85-0419
<b>CorPlex IL-22</b>	IL-22	0.01	0.156	0-400	1.66			85-0420
<b>CorPlex TNF-<math>\alpha</math></b>	TNF- $\alpha$	0.055	1.563	0-1,600	0.93	85-0409		

<sup>†</sup>E = EDTA plasma, S = serum

## SIMOA MULTIPLEX PLANAR ARRAY ASSAYS - HUMAN

Assay	Analytes	LoD (pg/mL)	LoQ (pg/mL)	Dynamic Range (pg/mL)	Median Endogenous (pg/mL)	Sample Volume	Sample Type	Catalog Number
Chemokine Panel 1 (4-plex)	IP-10	0.19	0.58	0-300	108	25 µL	E, S	88-5002
	ITAC	0.015	0.195	0-100	7.3			
	MCP-1	0.56	1.17	0-600	440.5			
	MIP-3b	0.005	0.78	0-400	66.3			
Angiogenesis Panel 1 (7-plex)	ANG-2	5.5	43.8	0-5,600	285	25 µL	E, S	85-0004
	HBEGF	0.014	0.2	0-125	5.6			
	VEGF-C	23.3	52.1	0-16,000	253			
	PIGF	0.2	0.5	0-500	7.7			
	FGFb	0.9	2.6	0-2,000	ND			
	HGF	1	6.3	0-6,300	134			
Biomarker Panel 1 (2-plex)	RANTES / CCL5	0.09	7.81	0-800	3,617	25 µL	E, S	85-0006
	TIMP-1	0.22	4.88	0-5,000	50,430			

## SIMOA MULTIPLEX PLANAR ARRAY DEVELOPER ASSAYS - NON-HUMAN

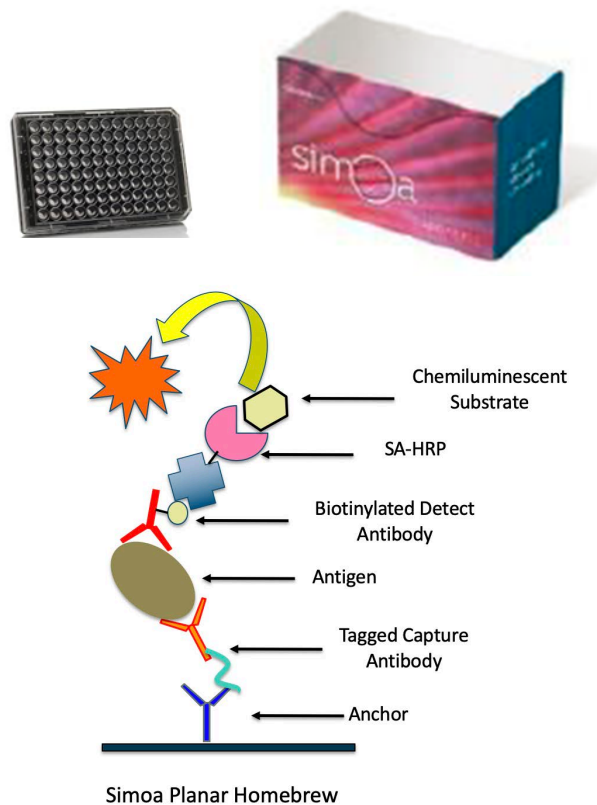
Species	Kit Name	Analytes	Catalog Number
Mouse	Mouse Cytokine Panel 1 (3-plex)	IL-10, IL-17, TNF-α	85-0450
Rat	Rat Cytokine Panel 1 (7-plex)	IFN-γ, IL-1β, IL-2, IL-6, IL-10, KC, TNF-α	85-0015
Rat	Rat Chemokine Panel 1 (4-plex)	MCP-1, MIP-1α, MIP-2, MIP-3α	85-0016
Non-Human Primate	NHP Cytokine Panel 1 (4-plex)	IFN-γ, IL-1β, IL-6, IL-8	85-0413

# SIMOA PLANAR HOMEBREW ASSAY DEVELOPMENT KITS

The Simoa Planar Array Homebrew Assay Starter kit is designed for customers developing Simoa homebrew planar array assays and contains all the components required for the development of a homebrew assay.

This kit includes sufficient materials to screen up to six antibody pairs (using one plate) and perform assay optimization (using two additional plates).

Item # 100-0461



The individual components of the Simoa Planar Array Homebrew Assay are designed for customers to scale up their Simoa homebrew planar array assays after initial HomeBrew Starter Kit has been tested. Some items included in the HomeBrew Starter kit can be purchased by a third-party vendor.

Planar Array Homebrew Plate (1-spot)	Simoa Planar Array homebrew plate (1-spot). 96 well plate for 1-plex homebrew assays. Qty 1.	197-0461
Simoa Planar Array Homebrew Tag 1	One vial of Simoa Planar Array Homebrew Tag 1 sufficient for preparing up to 100 plates	18910024
SuperSignal® Luminol	One bottle of SuperSignal Luminol reagent for processing one 96 well plate	1858089
SuperSignal® Peroxide	One bottle of SuperSignal Peroxide reagent for processing one 96 well plate	1858090
Simoa Planar Array Wash Buffer (25X)	25X Concentrate of Simoa Planar Array Wash Buffer (50 ml bottle)	1860055
Simoa Planar Array Homebrew Diluent A	Simoa Planar Array Homebrew Diluent A (25 ml bottle)	1880090

## SIMOA Developer Assay Kits

Developer assays take you beyond the HomeBrew to ready to use assay kits. Quanterix has screened several antibody pairs for each analyte and identified a working antibody pair and calibrator antigen. Standard curve can be generated and samples will read off the curve.

SP-X Assay Menu

### SIMOA PLANAR ARRAY DEVELOPER ASSAYS

Human Developer Assay				Rat, Mouse, & Porcine Assay			
Simoa Alpha-2 Macroglobulin Developer Kit	100-0429	Simoa MIG Developer Kit	100-0156	Simoa OPG Developer Kit	100-0025	Simoa porcine IL-1 $\beta$ Developer Kit	85-0393
Simoa Ang-1 Developer Kit	100-0067	Simoa MIP-1 $\beta$ Developer Kit	100-0351	Simoa PD-1 Developer Kit	100-0057	Simoa porcine IL-2 Developer Kit	100-0141
Simoa Amphiregulin Developer Kit	85-0369	Simoa Osteopontin Developer Kit	100-0002	Simoa PDGFAA Developer Kit	100-0394	Simoa porcine IL-6 Developer Kit	85-0394
Simoa BCA-1 (CXCL13) Developer Kit	100-0388	Simoa PD-L1 Developer Kit	100-0214	Simoa PDGFAB Developer Kit	100-0202	Simoa porcine IL-8 Developer Kit	85-0395
Simoa BMP-9 Developer Kit	100-0010	Simoa RAGE Developer Kit	100-0171	Simoa RANKL Developer Kit	100-0253		
Simoa C-Peptide Developer Kit	100-0292	Simoa TGF $\beta$ -1 Developer Kit	100-0021	Simoa SAA Developer Kit	100-0132		
Simoa CD14 Developer Kit	100-0001	Simoa TGF $\beta$ -2 Developer Kit	100-0105	Simoa SCF Developer Kit	100-0489		
Simoa CRP Developer Kit	100-0018	Simoa Tie-2 Developer Kit	100-0114	Simoa SDF-1 Developer Kit	100-0109		
Simoa EGF Developer Kit	85-0124	Simoa TSP-2 Developer Kit	100-0486	Simoa TGF $\alpha$ Developer Kit	85-0188		
Simoa Endoglin Developer Kit	100-0106	Simoa VEGF-D Developer Kit	100-0112	Simoa TIMP-2 Developer Kit	100-0189		
Simoa Eotaxin Developer Kit	100-0391	Simoa VEGF-R1 Developer Kit	100-0110	Simoa TNF-RI Developer Kit	100-0477		
Simoa Epregrulin Developer Kit	85-0366	Simoa VEGF-R2 Developer Kit	100-0030				
Simoa FasL Developer Kit	100-0332	Simoa VEGF-R3 Developer Kit	100-0111	Rat, Mouse, & Porcine Assay			
Simoa Fibrinogen Developer Kit	100-0012	Simoa MMP-9 Developer Kit	85-0021	Simoa rat MCP-1 Developer Kit	85-0370		
Simoa G-CSF Developer Kit	100-0304	Simoa Ang-2 Developer Kit	85-0108	Simoa rat MIP-1 $\alpha$ Developer Kit	85-0372		
Simoa gp130 Developer Kit	100-0386	Simoa FGFb Developer Kit	85-0280	Simoa rat MIP-2 Developer Kit	85-0373		
Simoa Granzyme B Developer Kit	100-0308	Simoa HB-EGF Developer Kit	85-0066	Simoa rat MIP-3 $\alpha$ Developer Kit	85-0374		
Simoa HGH Developer Kit	100-0293	Simoa HGF Developer Kit	85-0056	Simoa rat IFN $\gamma$ Developer Kit	85-0384		
Simoa I309 (CCL1) Developer Kit	100-0432	Simoa PDGFBB Developer Kit	85-0253	Simoa rat IL-10 Developer Kit	85-0385		
Simoa IgE Developer Kit	100-0457	Simoa PLGF Developer Kit	85-0352	Simoa rat IL-1 $\beta$ Developer Kit	85-0386		
Simoa IL-13 Developer Kit	100-0326	Simoa VEGF Developer Kit	85-0033	Simoa rat IL-2 Developer Kit	85-0387		
Simoa IL-15 Developer Kit	100-0485	Simoa VEGF-C Developer Kit	85-8049	Simoa rat KC/GRO $\alpha$ Developer Kit	85-0389		
Simoa IL-18 Developer Kit	100-0425	Simoa RANTES Developer Kit	85-0074	Simoa mouse MMP-9 Developer Kit	85-0096		
Simoa IL-1RA Developer Kit	100-0006	Simoa TIMP-1 Developer Kit	85-0161	Simoa mouse IL-10 Developer Kit	85-0310		
Simoa IL-1 $\alpha$ Developer Kit	100-0242	Simoa IP-10 Developer Kit	85-0018	Simoa mouse IL-17A Developer Kit	85-0196		
Simoa IL-2RG Developer Kit	100-0411	Simoa ITAC Developer Kit	85-0371	Simoa mouse VEGF Developer Kit	100-0325		
Simoa IL-3 Developer Kit	100-0331	Simoa MCP-1 Developer Kit	85-0039	Simoa mouse TNF $\alpha$ Developer Kit	85-0378		
Simoa IL-6R Developer Kit	100-0387	Simoa MIP-3 $\beta$ Developer Kit	85-0153	Simoa mouse ICAM-1 Developer Kit	85-0391		
Simoa IL-7 Developer Kit	100-0446	Simoa Insulin Developer Kit	100-0294	Simoa mouse IP-10 Developer Kit	85-0224		
Simoa KGF Developer Kit	100-0483	Simoa MDC Developer Kit	100-0488	Simoa mouse KC Developer Kit	100-0097		
Simoa M-CSF Developer Kit	100-0168	Simoa MIP-4 (PARC) Developer Kit	100-0066	Simoa mouse VCAM-1 Developer Kit	85-0392		
Simoa MIG Developer Kit	100-0156	Simoa MMP-2 Developer Kit	100-0011	Simoa porcine IFN $\alpha$ Developer Kit	85-0098		

### Custom Order Planar Array Assays

Custom singleplex and multiplex assay development services are provided by the Simoa Accelerator Laboratory. Please contact your Quanterix representative or email [info@quanterix.com](mailto:info@quanterix.com) to discuss custom assay development projects or sample testing services.

Quanterix Corporation  
900 Middlesex Turnpike, Billerica, MA 01821

**Quanterix™**  
The Science of Precision Health

©2021 Quanterix, Inc.  
Simoa™ and SP-X™ are registered trademarks of Quanterix, Inc.  
For research use only. Not for use in diagnostic procedures.

quanterix.com

472021 Rev1