



SP-X<sup>TM</sup> Ultrasensitive Imaging and Analysis System

Quanterix

# Ultrasensitive Biomarker Detection

Unleashing the power of next generation Simoa® planar array technology for robust multiplex circulating biomarker detection at the earliest stages of disease progression–even before recognizable symptoms begin

#### The value of The SP-X<sup>™</sup> System

The Quanterix SP-X<sup>™</sup> Imaging and Analysis System is a complete benchtop system that offers true multiplex detection at both acute and baseline levels in serum and plasma. Now oncology and immuno-oncology researchers and others who rely on multiplexing capabilities have an easy-to-use platform to help optimize workflows, speed up their research, and ultimately accelerate drug approvals.

# State of the art accuracy, sensitivity, precision, and reproducibility with chemiluminescent Simoa planar arrays



#### SP-X<sup>®</sup> unique capabilities

- · Easy to use and flexible data analysis software
- Open platform for custom ultra-sensitive assay development
- Combining Simoa ultra-sensitivity with 10-plex multiplexing
- Compact instrument footprint requiring only 11" (28 cm) of benchtop space
- Onboard high-resolution camera and custom lens for optimal light collection without user adjustments
- Optimized image analysis with proprietary acquisition algorithms and machine learning maximize dynamic range and S/N
- Touch screen tablet interface
- 21CFR Part 11 compatible workflow
- Imaging time in less than 10 minutes
- Low maintenance

### State of the Art Imaging and Data Analysis



Multiplex measurement of normal healthy baseline levels using the

Simoa CorPlex<sup>®</sup> Human Cytokine Panel 1on the SP-X system

### SP-X<sup>™</sup> Integrated Data Intelligence Program

1

Import panel and calibrator/sample dilution setup

Quanterix SPX	Quanterix SPX	Quanterix SPX	🤓 Quanterix SPX
Reports Tools Help	File Reports Tools Help	File Reports Tools Help	File Reports Tools Help
MONITOR LOAD LAYOUT SAVE LAYOUT TEN Catalog # 51-324-1-AB		SAVE LAYOUT SAVE LAYOUT SAVE LAYOUT	
Kit Catalog Custom Design	Catalog # 51-924-1-AB	Catalog # 51-924-1-AB	Catalog # 51-924-1-AB
51-924-1-AB Lookup Import A	Kit Catalog Cust	tom Design Kit Catalog Custom Design	Kit Catalog Custom Design
Empty B Empty	51-924-1-AB Lookup Impor	t <u>51-924-1-AB</u> Lookup Import A	51-924-1-AB Lookup Import
(12) °	Assay Units High C	Calibrator E Calibrators	Samples Import Map
Empty D	1 pg/ml 🗸	0 Dilution Factor C	Sample Name Dilution Factor C
9 Catalog # 51-924-1-AB 3 Empty	2 pg/ml Y		✓ 2
Species: Urknown			✓ 4
Empty	4 pg/ml Y		<ul> <li>✓ 10</li> </ul>
6 F	5 pg/ml 🗸		
G			
Plate: 96 Wells	7 pg/ml v		✓ 100 F
Plate: 96 Wells H	8 pg/ml V	256	✓ 1000
		600.0 S12 G	
	10 pg/ml Y	0 ●	
	11 pg/ml Y	H Map Dilutions to Analytes Select	Refresh & Purge
Plex Layout Assay Calibrators Samples	12 hMIP3b pg/ml 🗸	400.0	Remove Selected
rs'nehubiriOneDrive - QuanterivDesitlopi24-Feb-2022 Plate 1 Homebrew Plate 1 197			Vertical ↓ Horizontal →     Refresh Menu
	Plex Layout Assay Calibrators Sa	Plex Layout Assay Calibrators Samples	Piex Layout Assay Calibrators Samples
			2:\Users\nshubin\OneDrive - Quanterix\Desktop\24-Feb-2022 Plate 1 Homet



# 2 Build a plate map or load it from excel file

		SAVE LAYOUT	É TEMP										SP Quar	-X iterix
Catalog # 51-924-1-J														
Kit Catalog		Custom Design	1 Calbrator	2 Calibrator	3 Calibrator	4 Control	0 Control	6 Empty	7 Empty	8 Empty	9 Empty	10 Empty	11 Empty	12 Empty
51-924-1- <del>4</del> 8	Lookup	Import A												
			Calbrator	Calibrator	Calbrator	Control	Control							
							1:50							
			Calbrator	Calibrator	Calbrator	Sample 1	Sample 1							
	(12)					134	1.4							
							Sample 1							
9		3				14	14							
<b>.</b>		• •					Sample 1							
			1:128	1:128	1:128	14	14							
	6													
	•		1:255 Calibrator	1:255 Calibrator	1256 Celbreter	14	14							
			Calerator	Calorator			Sample 1							
			1:512 Cellbrator	1:512 Calibrator	1.512 Celbreter	1:4 Sample 6	1:4 Sample 1	Empty	Empty	Enoty	Evely	Empty	Empty	- Courter
Pla	ste: 96 Wells													
			1.0	1.0	1.0	14	1.4							_
Plex Layout Assa	y Calibrators	Samples												
rs'inshubiniOneOrive - Qu	anterixiDesktopi24-Fe	b-2022 Plate 1 Homebrew	Plate 1 197-04	61 3 19374P-05	2124SP0047 220	1224 145858 kit								

### **3** Open a .kit file

SAVELAYOUT											SF Quar	-X
Custom Design												
	Celbretor	Calbrator	Calibrator	Centrol	Cantrol							Em
Import												
	Calbrator	Calibrator	Calibrator	Certrol	Control							En
	Calbrator	Calibrator	Calibrator	Sample 1	Sample 1							
	Calbrator	Calibrator	Calibrator	Sample 2	Sample 1							
3	Calbrator	Calibrator	Calibrator	Sample 3	Sample 1							
	Calbrator	Calibrator	Calibrator	Sample 4	Sample 1							
	Calbrator	Calibrator	Calibrator	Sample 5	Sample 1							
	Calibrator	Calibrator	Calibrator	Sample 6	Sample 1							1
	SAVE LAYOUT Custom Design Import A Custom Design Custom Design Custom Design Custom Cu	SAVE LANGUT TERM Cottom Derign I I Control Derign I Control I I Control I I Control I I Control I I I Control I I I I Control I I I I I I I I I I I I I	SAVE LAVOUR         TUMALATES           Contrain Design         1           Opport         1           Contrain Design         1	SACE LAND?         TEMPLATE         PRIVEL AND/F           Outro Desire)         -         -         -           Concor         Salame         Concor         -           Import         -         1         -         -           Concor         Concor         Concor         Concor         Concor           10         10         10         10         Concor         Concor           13         10         Concor         T	Output         TEMPLATE         PRINTLANDOF         CERT           Caston Deepin	OWNER LINEOUT         TEMPLATES         INITE LINEOUT         CONTENTIONAL           Conton Design moort         1 <td>AVELAUXIT         TEMPLATE         PRINTLATION         DECRETIONAGE           CREAD Deprint         0         <td< td=""><td>South Langer         Time Arts         Initial Linear         Constrained         Official Linear         Officiar         Officiar         Offic</td><td>OWNER         COUNTY         TENDARTS         PRINTLANDAT         DECONTONIAGE         VETONIA           CARLIN DEATS         Carlier         Calabra         Carlier         Calabra         Carlier         Carlier</td><td>OWNER         TEMPLATE         PRIOR LANGET         DECRETIONALEE         PET FORM         MADE STATE           Contro Dealer mont        </td><td>Source Local         Titer.etts         Period         Description         Period         Period</td><td>OWNER LANGE         TEAL AND THE AND T</td></td<></td>	AVELAUXIT         TEMPLATE         PRINTLATION         DECRETIONAGE           CREAD Deprint         0 <td< td=""><td>South Langer         Time Arts         Initial Linear         Constrained         Official Linear         Officiar         Officiar         Offic</td><td>OWNER         COUNTY         TENDARTS         PRINTLANDAT         DECONTONIAGE         VETONIA           CARLIN DEATS         Carlier         Calabra         Carlier         Calabra         Carlier         Carlier</td><td>OWNER         TEMPLATE         PRIOR LANGET         DECRETIONALEE         PET FORM         MADE STATE           Contro Dealer mont        </td><td>Source Local         Titer.etts         Period         Description         Period         Period</td><td>OWNER LANGE         TEAL AND THE AND T</td></td<>	South Langer         Time Arts         Initial Linear         Constrained         Official Linear         Officiar         Officiar         Offic	OWNER         COUNTY         TENDARTS         PRINTLANDAT         DECONTONIAGE         VETONIA           CARLIN DEATS         Carlier         Calabra         Carlier         Calabra         Carlier         Carlier	OWNER         TEMPLATE         PRIOR LANGET         DECRETIONALEE         PET FORM         MADE STATE           Contro Dealer mont	Source Local         Titer.etts         Period         Description         Period         Period	OWNER LANGE         TEAL AND THE AND T

#### 4 Click "Process Kit"

File Reports Tools Help									1	$\sim$				
MONITOR Catalog # 51-924-1-	LOAD LAYOUT	SAVE LAYOUT	темр		PRINT LAYOUT				KITFOM	P	ROCESS KIT		SP	<b>-X</b> nterix:
Kit Catalog		Custom Design				4	5							
							Control							
51-924-1-AB	Lookup	Import				1;4	14							
							Control							
							1:50							
			Calibrator	Calibrator	Calbrator	Sample 1	Sample 1							
	12						14							
			Calbrator	Calbrator	Calbrator	Sample 2	Sample 1							
							14							
9	Catalog # 51-924-1-49	3	Calbrator	Calibrator	Celbrator	Sample 3	1.4 Sample 1	Empty	Emply	Empty	Empty	Empty	Empty	Empty
			1:125 Calbrator	1:125 Calibrator	1:125 Celbrator	1:4 Sample 4	1.4 Sample 1	Empty	Empty	Empty	Emply	Emply	Empty	Empty
	6		Calerator	Calorator	Celevator		Sample (	Chipty	CHEY	CHEVY	CHEV	CHUY		empty
	•		1.256	1256		1.6	194							
							Sample 1							
						1.6	14							
Р	late: 96 Wells						Sample 1							Enoty
							1.4							
Plex Layout Assa	ay Calibrators	Samples												

# 5 Review data and output report

ew Plate Layout pen a SPX Image File pen Recent pen a SPX File from an Imager sAVE LAYOUT		TEMPLATES		PRINT LAYOUT				R KIT FORM					SP-X Quanterix		
Save 76-SBX Kit File Close SPX Kit File															
1 UC	Custom Design	1 Celbreter	2 Calibrator	3 Celibrator	4 Centrol	5 Cantrol	6 Empty	7 Empty	8 Empty	9 Empty	10 Enpty	11 Empty	12 Empte		
51-924-1-AB Lookup	Import					1.4									
		Celbrator	Calibrator	Calibrator	Centrol	Control	Empty						Empl		
						150									
		Calbrator	Calibrator	Calibrator	Sample 1	Sample 1	Empty						Emp		
12						1.4									
						Sample 1	Empty						Eng		
Catalog #	3				1:4	1.4									
51824-148 Species:	- 👻 💡					Sample 1	Empty						Eng		
Unknown		1:128 Celbretor	1:128 Calibrator	1:128 Calibrator	1:4 Sample 4	1.4 Sample 1	Empty	Empty	Empty	Empty	Enpty	Empty	Emp		
6							0404								
		1.258 Calbrator	1:258 Calibrator	1:258 Calibrator	1:4 Sample 5	1.4 Sample 1	Empty	Empty	Empty	Empty	Empty	Empty	Emp		
						1.4									
Plate: 96 Wells		Calbretor	Calibrator	Calibrator	Sample 6	Sample 1	Empty						Eng		
						14									



# Simoa SP-X<sup>™</sup> workflow: simple, scalable, flexible.



Add reagents and samples to assay plate



Incubate and shake assay plate



Automated Wash of assay plate



Load assay plate onto SP-X<sup>™</sup> for quick imaging

### How the SP-X<sup>™</sup> Achieves Ultrasensitivity

The Simoa planar immunoassay technology is a revolutionary new digital biomarker solution, with features that provide researchers an incredibly simple, flexible, robust, and sensitive multiplexing platform.

#### Proprietary high-precision digital nanofluidic antibody deposition technology

Provides unprecedented surface chemistry optimization, minimizing non-specific binding and resulting in low background noise and excellent assay precision and accuracy.

#### Unique spot design

Antibodies are deposited into discrete spots in a circular pattern around the perimeter of each round microtiter plate well. Each spot contains the capture antibody for one target analyte. With enough space for 10 unique spots in a single well, up to a 10-plex assay can be performed saving sample volume, time, and cost without sacrificing assay performance characteristics.

# Vortex interaction between analyte molecules and capture antibodies

The Simoa planar approach of using capture antibodies spotted on a circular pattern combined with the fluid dynamics of vortex mixing, maximizes sensitivity by driving binding reactions to equilibrium and avoiding depletion regions in the center of the well

### Imaging occurs through bottom of translucent wells

The Simoa surface of the planar array plates are manufactured using translucent materials enabling imaging of the array through the bottom of the plate, eliminating the potential for optical signal deformations and for reduced sensitivity of the reading due to the meniscus of the liquid in the well.

#### Proprietary algorithm to calculate optimal exposure times for each individual plate

Proprietary algorithm allows the SP-X<sup>™</sup> Imager to automatically optimize exposure time and number of images to maximize sensitivity and dynamic range for each experiment.number of images to maximize sensitivity and dynamic range for each experiment.

Visit: quanterix.com/SP-X for more information

