

Product Number:	101643
Lot Number:	503988
Expiration:	18-Jul-2024
Platform(s):	HD-X



Data below represents results generated on the Simoa® HD-X

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	101645	327605	N/A
Detector Reagent	101644	327104	N/A
SBG Reagent (HD-1)	101646	327105	N/A
Sample Diluent	102200	319920	319920A
Calibrator Diluent	101648	324512	N/A
Calibrator Concentrate	100670	315305	315305A
SBG Diluent (SR-X) †	103158	N/A ¹	N/A
RGP Reagent	103159	N/A ¹	N/A

¹ These items are not Kit Lot Specific

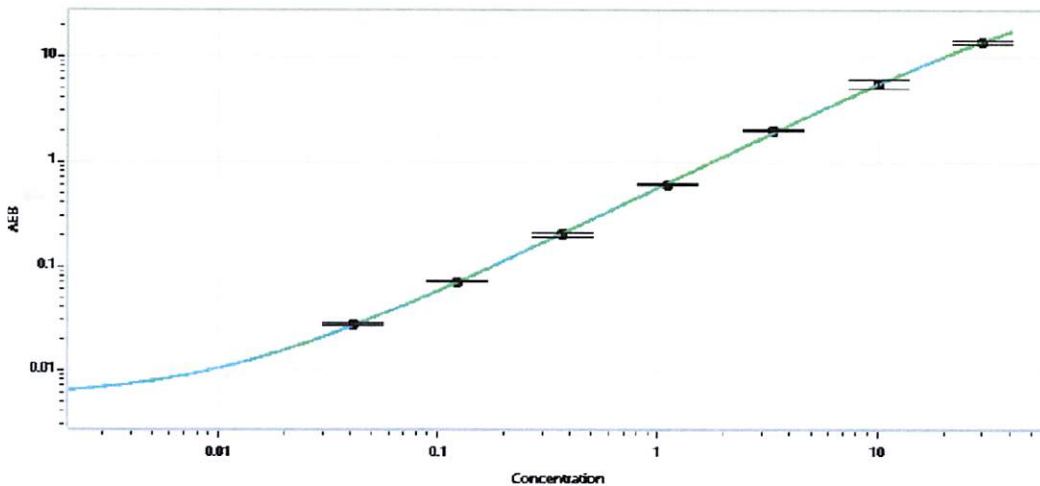
† Specific for use on the SR-X platform - refer to the Simoa™ Human IL-10 Advantage SR-X Kit Instructions (KI-0036).

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	315305	25.6	18.0 - 30.0
Control 2	315305	1.76	1.20 - 2.00
Panel 1	215202	0.465	0.213 - 0.561
Panel 2	215202	0.655	0.425 - 0.738
Panel 3	306507	63.5	47.1 - 71.5

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	603
---	-----

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.0411
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Sonia Khillan
Name

QC

Title

Sonia Khillan 06 OCT 2023
Signature/Date