

Product Number:	101643
Lot Number:	503533
Expiration:	21-Jun-2023
Platform(s):	HD-1/HD-X and SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101645	218813
Detector Reagent	101644	218901
SBG Reagent (HD-1)	101646	218917
Sample Diluent	102200	217218
Calibrator Diluent	101648	218919
Calibrator Concentrate	100670	218916
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ These items are not Kit Lot Specific

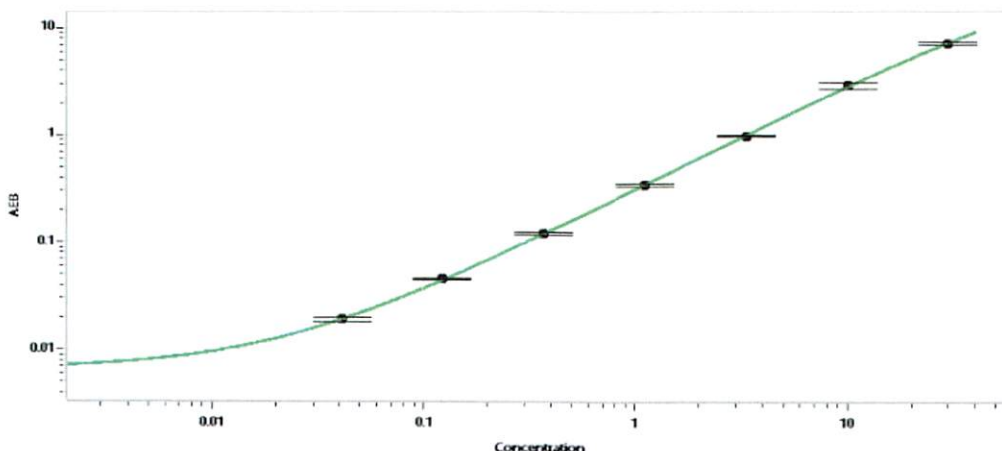
† Specific for use on the SR-X platform - refer to the Simoa™ Human IL-10 Advantage SR-X Kit Instructions (KI-0036).

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	308004	1.62	1.20 - 2.00
Control 2	308004	25.1	18.0 - 30.0
Panel 1	215202	0.428	0.260 - 0.553
Panel 2	215202	0.684	0.432 - 0.748
Panel 3	306507	58.3	48.4 - 72.6

* 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	0.041
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

Soni 12/16 11 APR 2023
Signature/Date

Signature/Date