

Product Number:	101643
Lot Number:	503850
Expiration:	13-Jan-2024
Platform(s):	HD-1/HD-X and SR-X



Component	Part Number	Lot Number
Bead Reagent	101645	307208
Detector Reagent	101644	306608
SBG Reagent (HD-1)	101646	306211
Sample Diluent	102200	306117
Calibrator Diluent	101648	304412
Calibrator Concentrate	100670	308004
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).

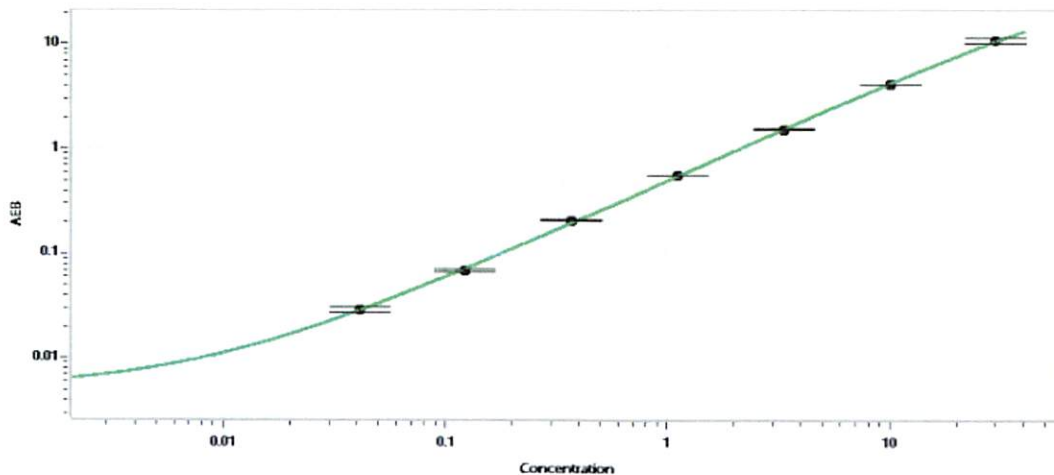
Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	308004	1.51	1.20 - 2.00
Control 2	308004	24.7	18.0 - 30.0
Panel 1	215202	0.497	0.260 - 0.553
Panel 2	215202	0.607	0.432 - 0.748
Panel 3	306507	62.6	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	0.0411
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 12/11/23 11 APR 2023
Signature/Date