

Product Number:	101643
Lot Number:	503146
Expiration:	1-Jun-2022
Platform(s):	HD-1/HD-x and SR-X

Component	Part Number	Lot Number
Bead Reagent	101645	118501
Detector Reagent	101644	126801
SBG Reagent (HD-1)	101646	118603
Sample Diluent	102200	115222
Calibrator Diluent	101648	117710
Calibrator Concentrate	100670	118903
SBG Diluent (SRX) †	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).



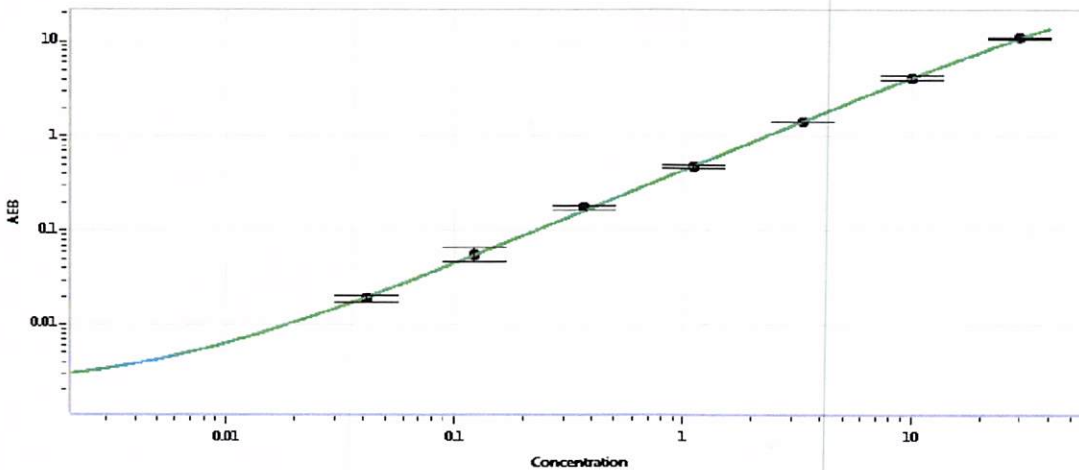
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	105704	1.48	0.830 - 1.61
Control 2	105704	26.8	15.6 - 31.5
Panel 1	029001	0.795	0.436 - 1.18
Panel 2	029001	1.19	0.685 - 1.49
Panel 3	029001	15.6	10.2 - 20.3

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

Calibrator Concentrate Value Assignment (pg/mL)	187
---	-----

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	Junpil Park	QC Supervisor	
	Name	Title	Signature/Date