

Product Number:	101643
Lot Number:	502832
Expiration:	31-Mar-2022
Platform(s):	HD-1/HD-x and SR-X

Component	Part Number	Lot Number
Bead Reagent	101645	109004
Detector Reagent	101644	109010
SBG Reagent (HD-1)	101646	109011
Sample Diluent	102200	109015
Calibrator Diluent	101648	109019
Calibrator Concentrate	100670	108904
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is 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).

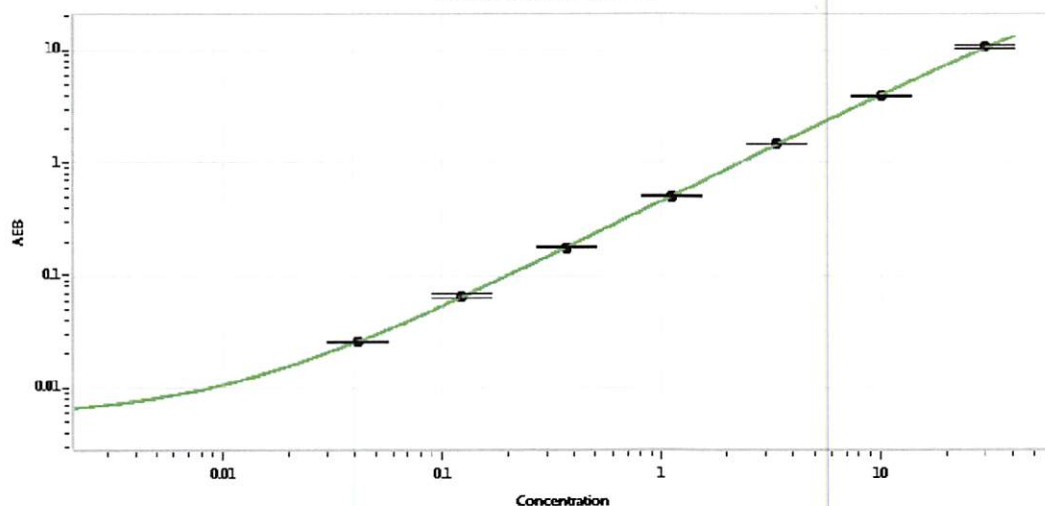


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	105704	1.01	0.805-1.65	
Control 2	105704	23.1	14.7-32.2	
Panel 1	029001	0.549	0.436-1.18	
Panel 2	029001	0.777	0.685-1.49	
Panel 3	029001	16.4	10.2-20.3	

Calibrator Concentrate Value Assignment (pg/mL)	506
---	-----

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
Name

QC Supervisor
Title

[Signature]
Signature/Date
03 May 2021