

Product Number:	101643
Lot Number:	503185
Expiration:	30-Jul-2022
Platform(s):	HD-1/HD-x and SR-X

Component	Part Number	Lot Number
Bead Reagent	101645	128605
Detector Reagent	101644	128705
SBG Reagent (HD-1)	101646	128706
Sample Diluent	102200	121115
Calibrator Diluent	101648	124010
Calibrator Concentrate	100670	128707
SBG Diluent (SRX) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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).

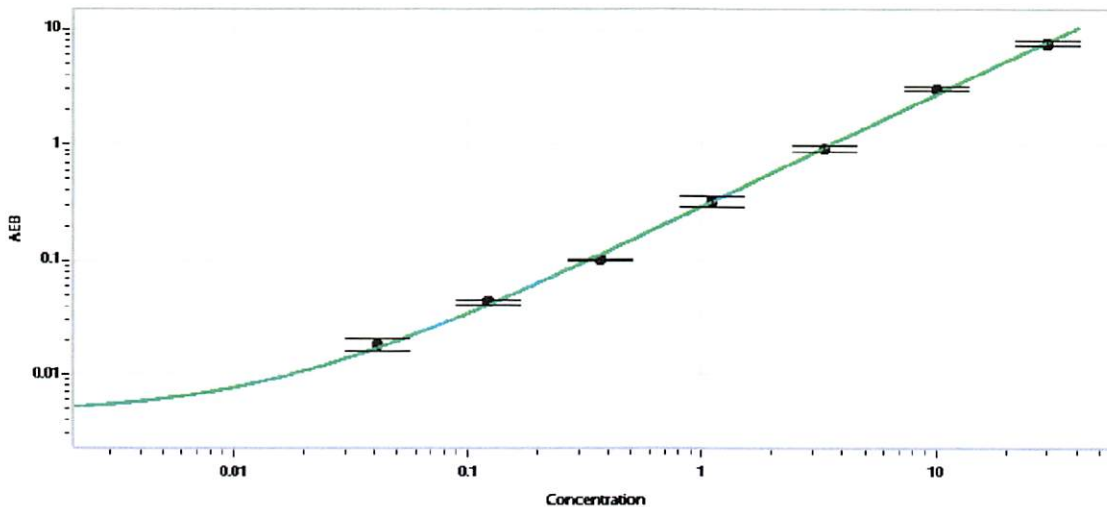


Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials*				* 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	1.32	0.830 - 1.61	
Control 2	105704	29.0	15.6 - 31.5	
Panel 1	029001	0.572	0.436 - 1.18	
Panel 2	029001	0.774	0.685 - 1.49	
Panel 3	029001	11.3	10.2 - 20.3	

Calibrator Concentrate Value Assignment (pg/mL)	300
---	-----

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]* 01 NOV 2021  
Signature/Date