

Product Number:	101643
Lot Number:	502757
Expiration:	11-Jan-2022
Platform(s):	HD-1/HD-x and SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101645	104611
Detector Reagent	101644	104603
SBG Reagent (HD-1)	101646	106419
Sample Diluent	102200	104616
Calibrator Diluent	101648	102922
Calibrator Concentrate	100670	105603
SBG Diluent (SRX) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

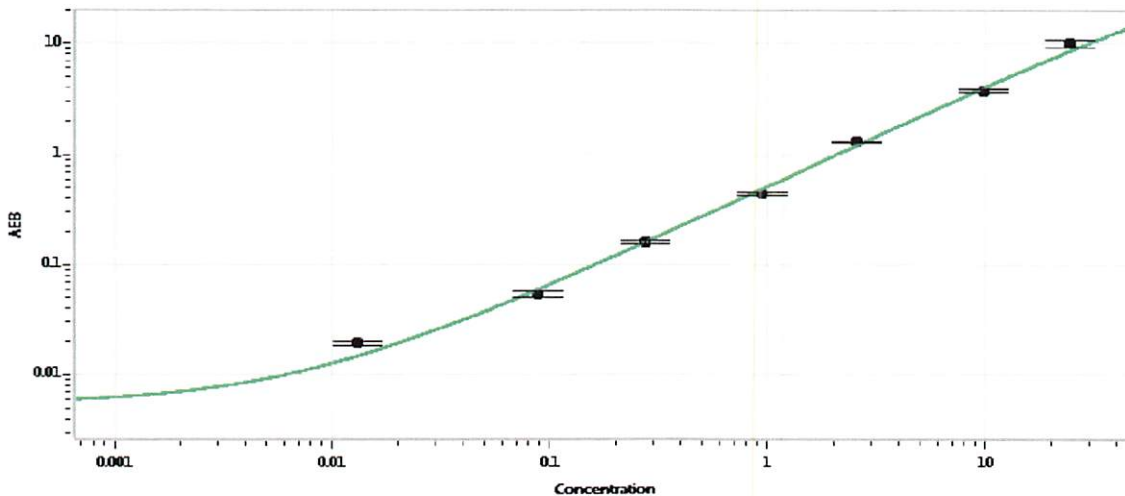
¹ RGP is not Kit Lot Specific

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

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	1.40	0.879-1.61	
Control 2	105704	25.6	12.1-33.7	
Panel 1	029001	0.725	0.611-1.11	
Panel 2	029001	1.06	0.917-1.38	
Panel 3	029001	11.4	9.69-20.4	

Calibrator Concentrate Value Assignment (pg/mL)	468
---	-----

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/Date
22 Mar 2021