

Product Number:	101643
Lot Number:	503017
Expiration:	3-Feb-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	117404
Detector Reagent	101644	117702
SBG Reagent (HD-1)	101646	117706
Sample Diluent	102200	110014
Calibrator Diluent	101648	115922
Calibrator Concentrate	100670	003413
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

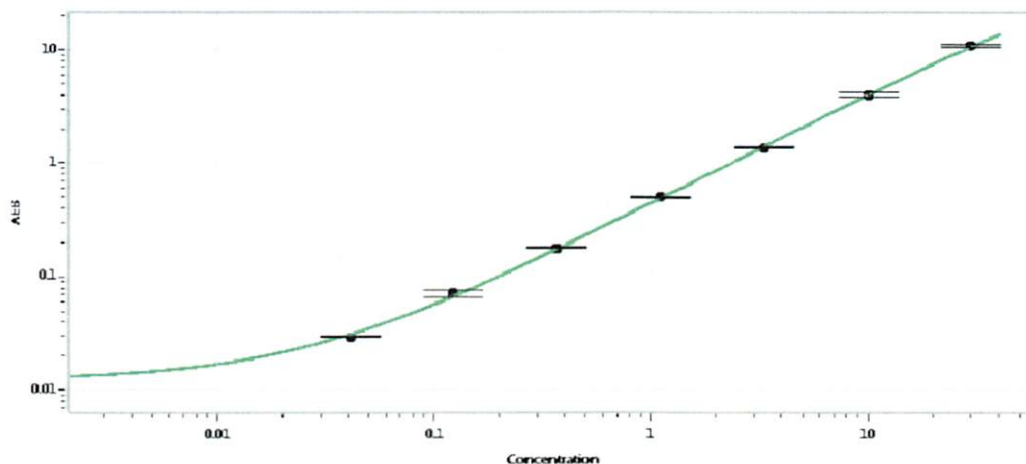
† Specific for use on the SR-X platform - refer to the Simoa™ Human IL-10 Advantage SR-X Kit Instructions (KI-0036).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	105704	1.18	0.830-1.61
Control 2	105704	23.9	15.6-31.5
Panel 1	029001	0.876	0.436-1.18
Panel 2	029001	1.15	0.685-1.49
Panel 3	029001	17.3	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)	817
---	-----

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

 08 Jul 2021  
Signature/Date