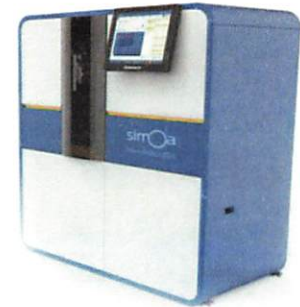


Product Number:	101643
Lot Number:	503167
Expiration:	28-Jun-2022
Platform(s):	HD-1/HD-x and SR-X

Component	Part Number	Lot Number
Bead Reagent	101645	121901
Detector Reagent	101644	127404
SBG Reagent (HD-1)	101646	122002
Sample Diluent	102200	117922
Calibrator Diluent	101648	122007
Calibrator Concentrate	100670	122003
SBG Diluent (SRX) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>



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

<sup>1</sup> These items are not Kit Lot Specific

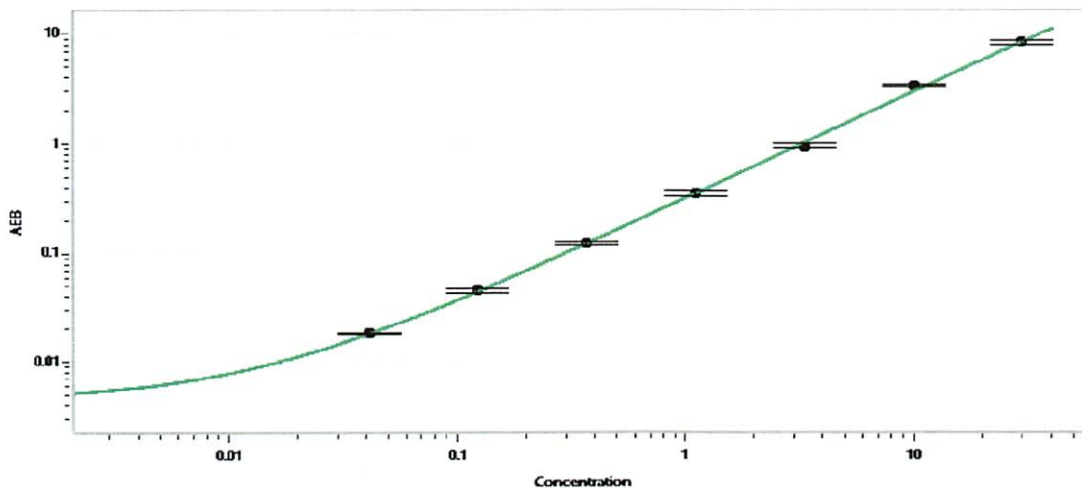
† Specific for use on the SR-X platform - refer to the Simoa™ Human IL-10 Advantage SR-X Kit Instructions (KI-0036).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	105704	1.38	0.830 - 1.61
Control 2	105704	26.6	15.6 - 31.5
Panel 1	029001	0.610	0.436 - 1.18
Panel 2	029001	0.852	0.685 - 1.49
Panel 3	029001	12.9	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)	278
---	-----

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

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel (20 Oct 2021)  
Signature/Date