

Product Number:	101643
Lot Number:	502715
Expiration:	14-Dec-2021
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	102701
Detector Reagent	101644	102704
SBG Reagent (HD-1)	101646	102907
Sample Diluent	102200	034915
Calibrator Diluent	101648	100704
Calibrator Concentrate	100670	003413
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

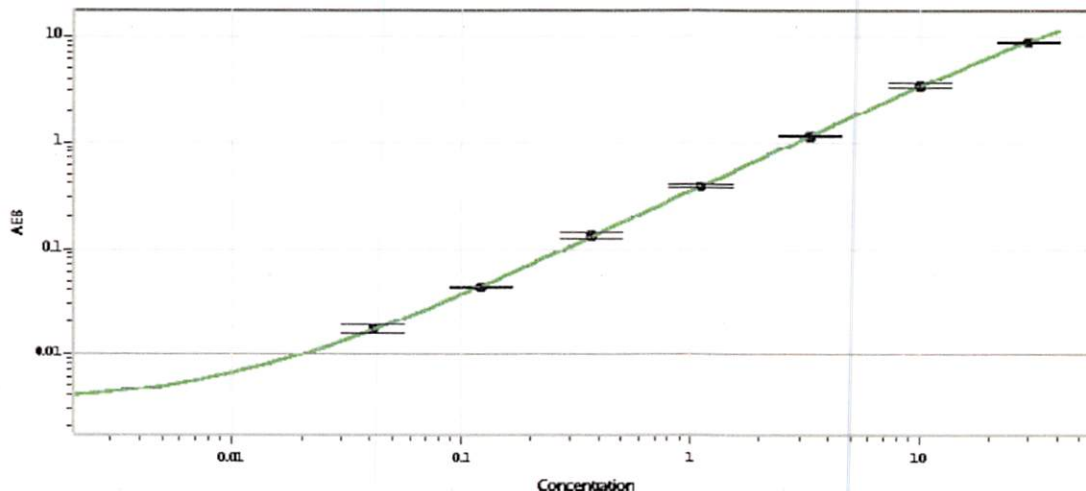
† Specific for use on the SR-X platform - refer to the Simoa™ Human IL-10 Advantage SR-X Kit Instructions (KI-0036).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-060518-03	1.29	0.740-1.52
Control 2	MF-060518-03	27.1	11.1-32.3
Panel 1	029001	1.00	0.701-1.09
Panel 2	029001	1.21	0.934-1.40
Panel 3	029001	17.2	12.8-19.2

* 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

Shivani Goel
Name

QC Manager
Title

Shivani Goel 26 Feb 2021
Signature/Date