

Product Number:	101643
Lot Number:	501757
Expiration:	21-Dec-2019
Platform(s):	HD-1 and SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101645	905812
Detector Reagent	101644	905808
SBG Reagent (HD-1)	101646	905809
Sample Diluent	102200	902931
Calibrator Diluent	101648	835505
Calibrator Concentrate	100670	909504
SBG Diluent (SRX) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

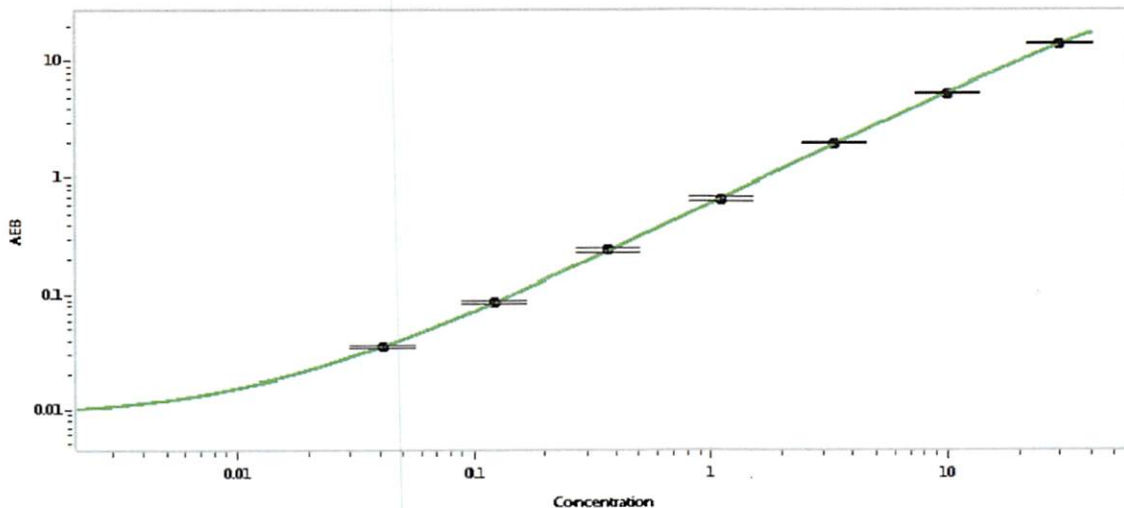
¹ Reagents are not Kit Lot Specific

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

Release Materials				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	MF-060518-03	1.10	0.740-1.52	
Control 2	MF-060518-03	23.2	11.1-32.3	
Panel 1	MF-060718-01	0.378	0.188-0.563	
Panel 2	MF-060718-02	0.999	0.536-1.32	
Panel 3	MF-060718-03	61.2	49.8-89.0	

Calibrator Concentrate Value Assignment
1057 pg/mL

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	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	10 Apr 2019
	Name	Title	Signature/Date	