

Product Number:	101643
Lot Number:	501508
Expiration:	19-Sep-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	832003
Detector Reagent	101644	832004
SBG Reagent (HD-1)	101646	831814
Sample Diluent	102200	826218
Calibrator Diluent	101648	827413
Calibrator Concentrate	100670	833805
SBG Diluent (SRX) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

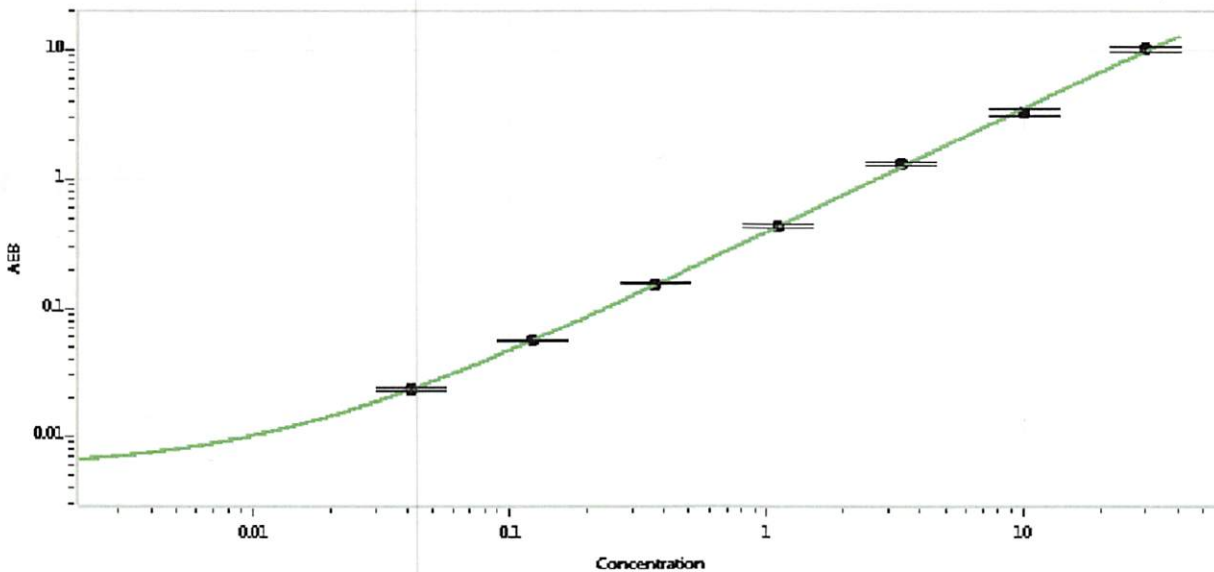
<sup>1</sup> Reagents are not Kit Lot Specific

† 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	1.21	0.753-1.31	
Control 2	MF-060518-03	23.9	11.8-25.9	
Panel 1	MF-060718-01	0.361	0.220-0.436	
Panel 2	MF-060718-02	0.917	0.610-1.04	
Panel 3	MF-060718-03	74.5	48.7-84.6	

**Calibrator Concentrate Value Assignment**  
209 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

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 17DEC2018  
Signature/Date