

Product Number:	101643
Lot Number:	501311
Expiration:	10-Apr-2019
Platform(s):	HD-1 and SR-X



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

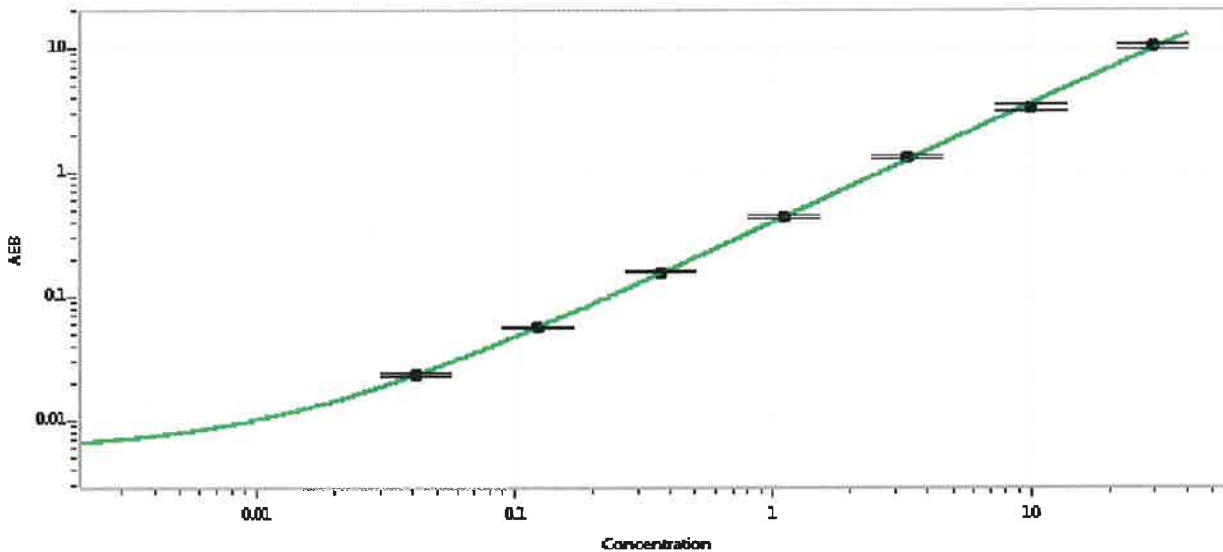
Component	Part Number	Lot Number
Bead Reagent	101645	818408
Detector Reagent	101644	81787
SBG Reagent (HD-1)†	101646	81792
Sample Diluent	102200	81007
Calibrator Diluent	101648	81298
Calibrator Concentrate	100670	73482
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

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	52232	1.64	0.741-1.88	
Control 2	52232	27.5	12.7-29.6	
Panel 1	MF-011916-08	0.557	0.235-0.621	
Panel 2	MF-011916-09	2.18	1.21-2.32	
Panel 3	MF-011916-10	58.1	24.9-63.3	

Calibrator Concentrate Value Assignment
589 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, S. Murugan 13/4/2018 Signature/Date