

Product Number:	101643
Lot Number:	501695
Expiration:	19-Sep-2019
Platform(s):	HD-1 and SR-X



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

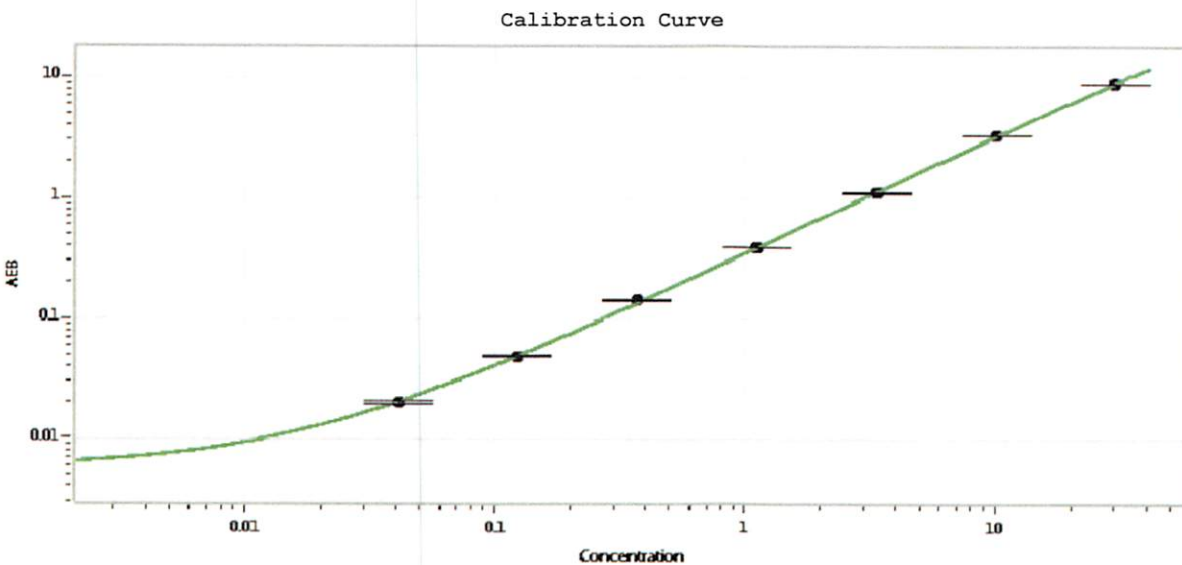
Component	Part Number	Lot Number
Bead Reagent	101645	907107
Detector Reagent	101644	906302
SBG Reagent (HD-1)	101646	833712
Sample Diluent	102200	826222
Calibrator Diluent	101648	827414
Calibrator Concentrate	100670	833805
SBG Diluent (SRX) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

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	1.25	0.740-1.52	
Control 2	MF-060518-03	23.7	11.1-32.3	
Panel 1	MF-060718-01	0.522	0.188-0.563	
Panel 2	MF-060718-02	1.23	0.536-1.32	
Panel 3	MF-060718-03	64.8	49.8-89.0	

Calibrator Concentrate Value Assignment
209 pg/mL



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

P. Stadler 15 Nov 2019

Name

Title

Signature/Date