

Product Number:	101643
Lot Number:	501801
Expiration:	21-Dec-2019
Platform(s):	HD-1 and SR-X

Component	Part Number	Lot Number
Bead Reagent	101645	911911
Detector Reagent	101644	912106
SBG Reagent (HD-1)	101646	9127107
Sample Diluent	102200	908908
Calibrator Diluent	101648	835510
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).



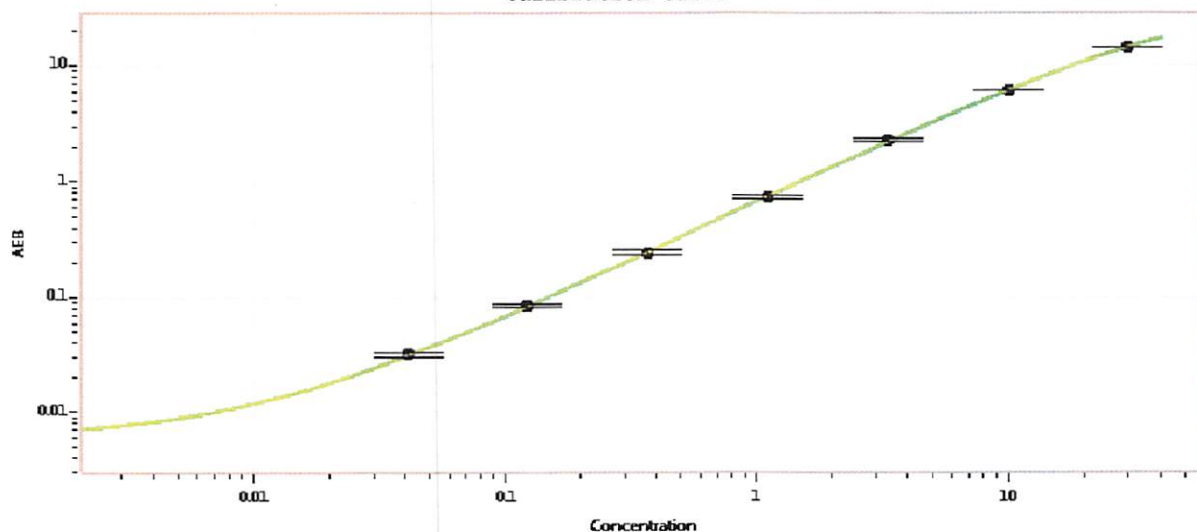
Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	1.16	0.740-1.52	
Control 2	MF-060518-03	22.2	11.1-32.3	
Panel 1	MF-060718-01	0.436	0.188-0.563	
Panel 2	MF-060718-02	1.04	0.536-1.32	
Panel 3	MF-060718-03	79.5	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

Name

Title

Signature/Date

P. Stadler

09 May 2019