

Product Number:	101643
Lot Number:	502126
Expiration:	29-Apr-2020
Platform(s):	HD-1/HD-x and SR-X

Component	Part Number	Lot Number
Bead Reagent	101645	911911
Detector Reagent	101644	912106
SBG Reagent (HD-1)	101646	912107
Sample Diluent	102200	928023
Calibrator Diluent	101648	932212
Calibrator Concentrate	100670	909504
SBG Diluent (SRX)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is 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).



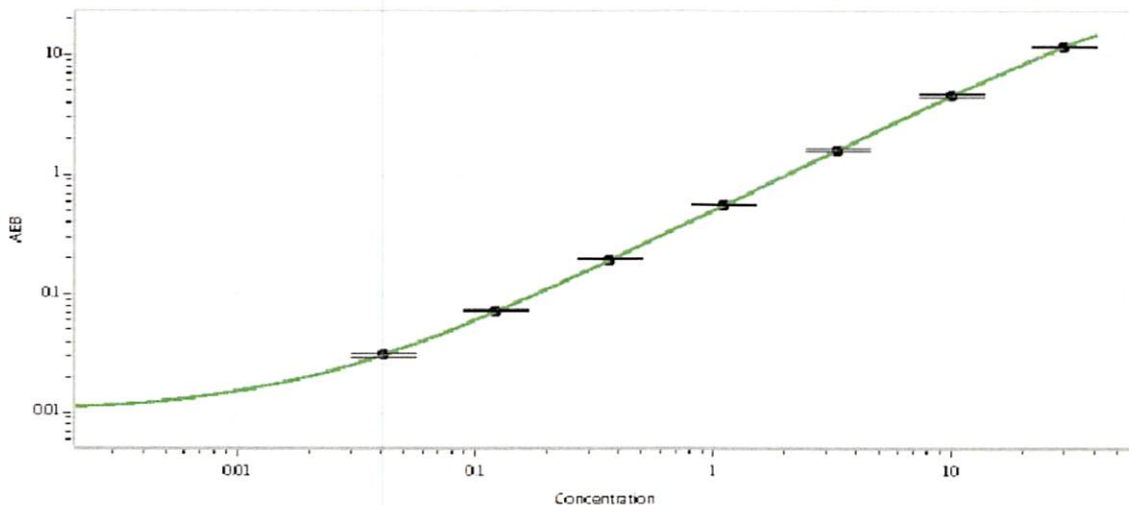
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-060518-03	1.30	0.740-1.52
Control 2	MF-060518-03	24.8	11.1-32.3
Panel 1	MF-060718-01	0.386	0.188-0.563
Panel 2	MF-060718-02	1.000	0.536-1.32
Panel 3	MF-060718-03	74.4	49.8-89.0

\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	1057
---	------

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

*P. Stadler* 09 Dec 2019

Name

Title

Signature/Date