

Product Number:	101643
Lot Number:	501181
Expiration:	30-Jan-2019
Platform(s):	HD-1 and SR-X

Component	Part Number	Lot Number
Bead Reagent	101645	80751
Detector Reagent	101644	80680
SBG Reagent (HD-1)†	101646	80652
Sample Diluent	102200	80653
Calibrator Diluent	101648	80654
Calibrator Concentrate	100670	73482
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific



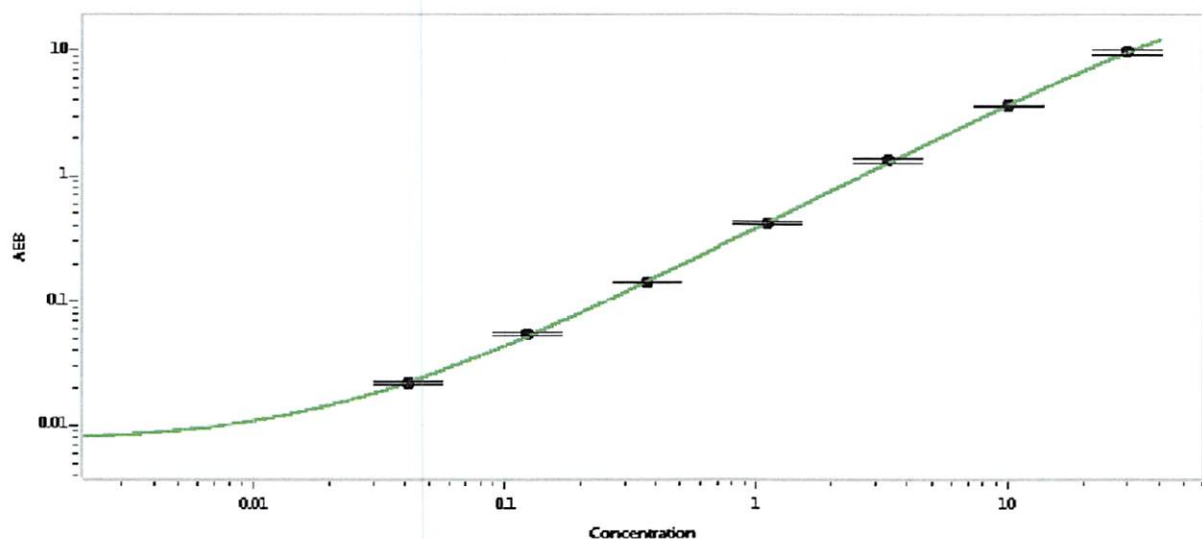
Data below represents results generated on the Simoa™ HD-1 Analyzer.

† Specific for use on the SRX platform – refer to the Simoa™ Human IL-10 Advantage SRX 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	1.50	0.753-1.78	
Control 2	52232	21.6	14.1-26.6	
Panel 1	MF-011916-08	0.332	0.251-0.585	
Panel 2	MF-011916-09	1.54	1.26-2.21	
Panel 3	MF-011916-10	43.4	25.7-59.8	

Calibrator Concentrate Value Assignment (pg/mL)	589
---	-----

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
Signature/Date 30 MAR 2018