

Product Number:	100198
Lot Number:	501254
Expiration:	23-Oct-2018
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	100815	81151
Detector Reagent	100816	81152
SBG Reagent	100817	81353
Sample Diluent	103204	81133
Calibrator Diluent	100829	81134
Calibrator Concentrate	100828	73380
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

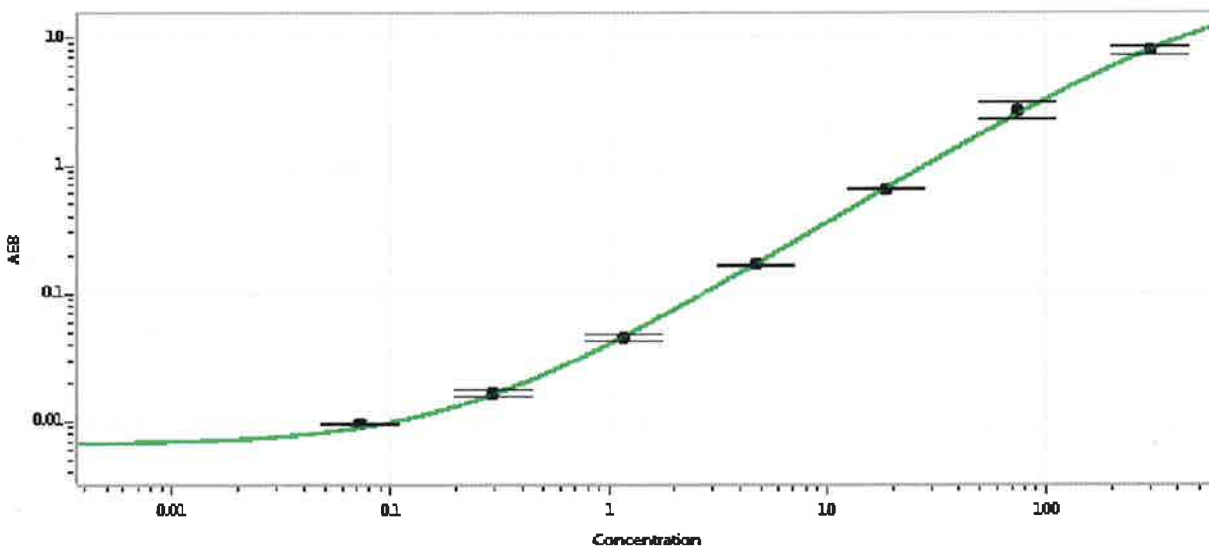
† Specific for use on the SRX platform - refer to the Simoa™ IL-8 SR-X Kit Instructions (KI-0092).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-091217-02	10.1	7.76-11.6
Control 2	MF-091217-02	172	147-220
Panel 1	MF-061617-01	5.94	3.46-6.45
Panel 2	MF-061617-02	52.1	42.0-62.9
Panel 3	MF-061617-03	339	279-459

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

Calibrator Concentrate Value Assignment (pg/mL)	15,328
---	--------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.073
C	0.293
D	1.17
E	4.69
F	18.8
G	75.0
H	300

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 20/10/2018
Signature/Date