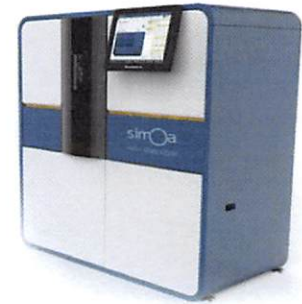


Product Number:	100198
Lot Number:	501432
Expiration:	17-Mar-2019
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	100815	827102
Detector Reagent	100816	826007
SBG Reagent	100817	826002
Sample Diluent	103204	826003
Calibrator Diluent	100829	826004
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

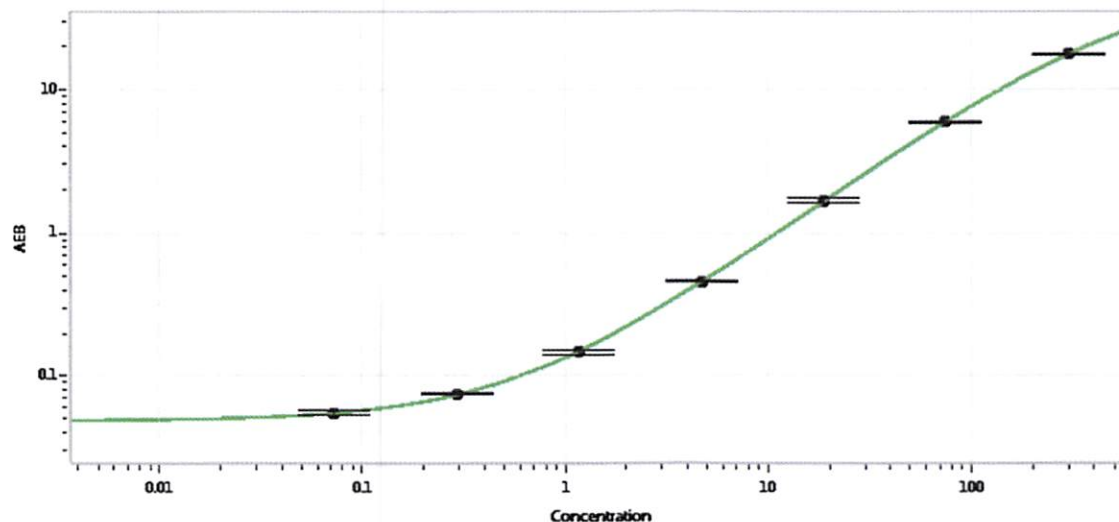
† Specific for use on the SRX platform - refer to the Simoa™ IL-8 SR-X Kit Instructions (KI-0092).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-091217-02	9.74	7.76-11.6
Control 2	MF-091217-02	166	145-217
Panel 1	MF-061617-01	4.43	3.41-6.37
Panel 2	MF-061617-02	50.1	41.7-62.6
Panel 3	MF-061617-03	309	261-459

* 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/Date

[Handwritten Signature] 05 OCT 2018