

Product Number:	100198
Lot Number:	502627
Expiration:	17-Jun-2021
Platform(s):	HD-1/HD-x & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	100815	032401
Detector Reagent	100816	032503
SBG Reagent	100817	032505
Sample Diluent	103204	032506
Calibrator Diluent	100829	032501
Calibrator Concentrate	100828	016905
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

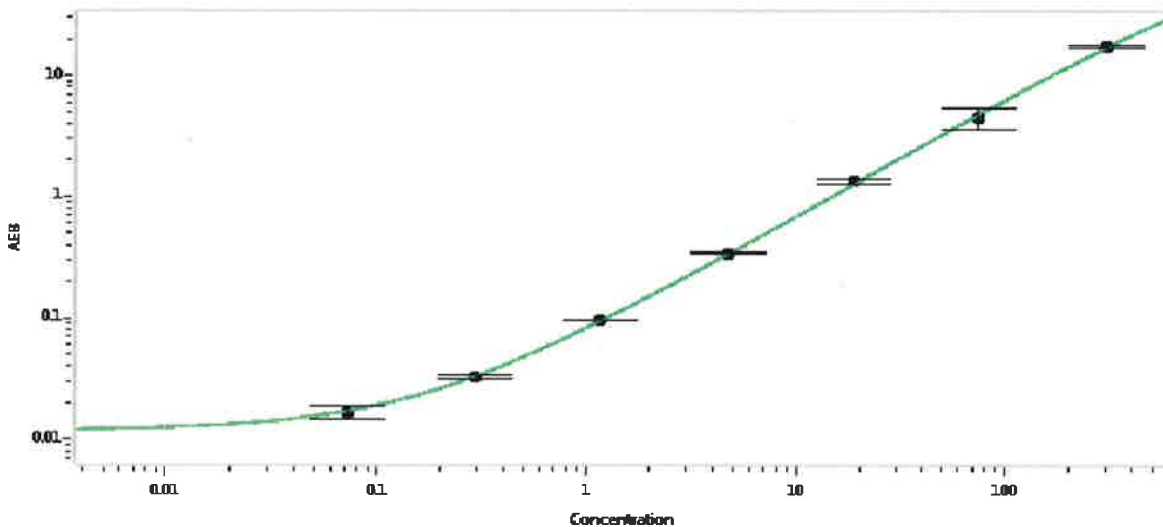
† Specific for use on the SRX platform - refer to the Simoa™ IL-8 SR-X Kit Instructions (KI-0092).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-091217-02	12.0	7.54-12.5
Control 2	MF-091217-02	219	137-234
Panel 1	032901	4.76	3.17-6.50
Panel 2	032901	38.5	27.2-56.5
Panel 3	033005	484	358-660

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

Calibrator Concentrate Value Assignment (pg/mL)	5,050
---	-------

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	Junpil Park	QC Supervisor	
	Name	Title	Signature/Date

[Handwritten Signature] 28 Dec 2020