

Product Number:	100198
Lot Number:	502521
Expiration:	3-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	016810
Detector Reagent	100816	016805
SBG Reagent	100817	016804
Sample Diluent	103204	016807
Calibrator Diluent	100829	016803
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

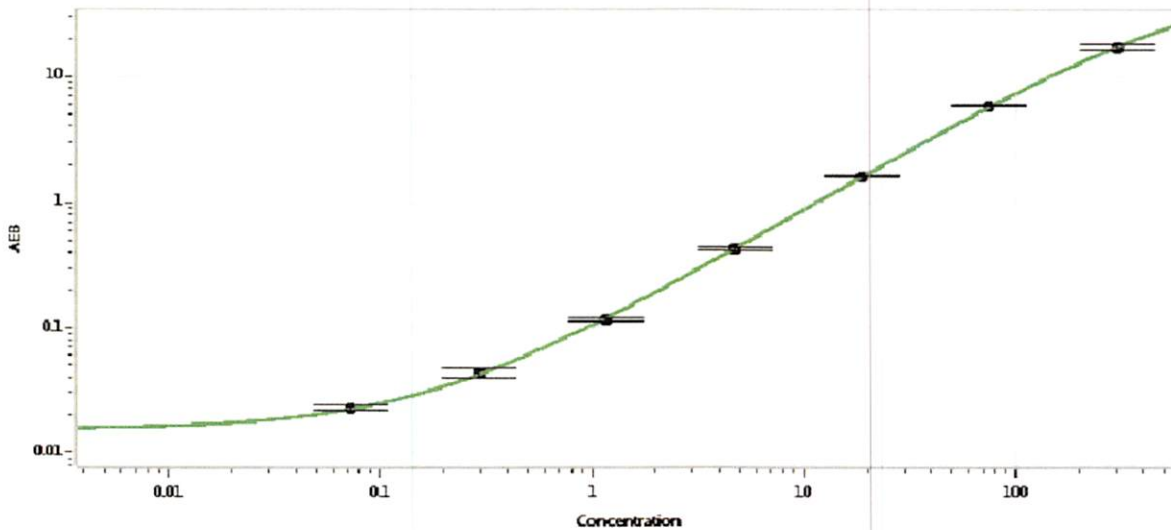
† Specific for use on the SRX platform - refer to the Simoa™ IL-8 SR-X Kit Instructions (KI-0092).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-091217-02	9.97	7.70-11.6
Control 2	MF-091217-02	186	143-220
Panel 1	MF-061617-01	5.05	3.17-6.91
Panel 2	MF-061617-02	54.5	42.6-63.9
Panel 3	MF-061617-03	311	267-459

* 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	Shivani Goel	QC Manager	Shivani Goel 08 Sep 2020
	Name	Title	Signature/Date