

Product Number:	100198
Lot Number:	502227
Expiration:	15-Jan-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	007606
Detector Reagent	100816	007603
SBG Reagent	100817	007701
Sample Diluent	103204	007602
Calibrator Diluent	100829	007601
Calibrator Concentrate	100828	007903
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

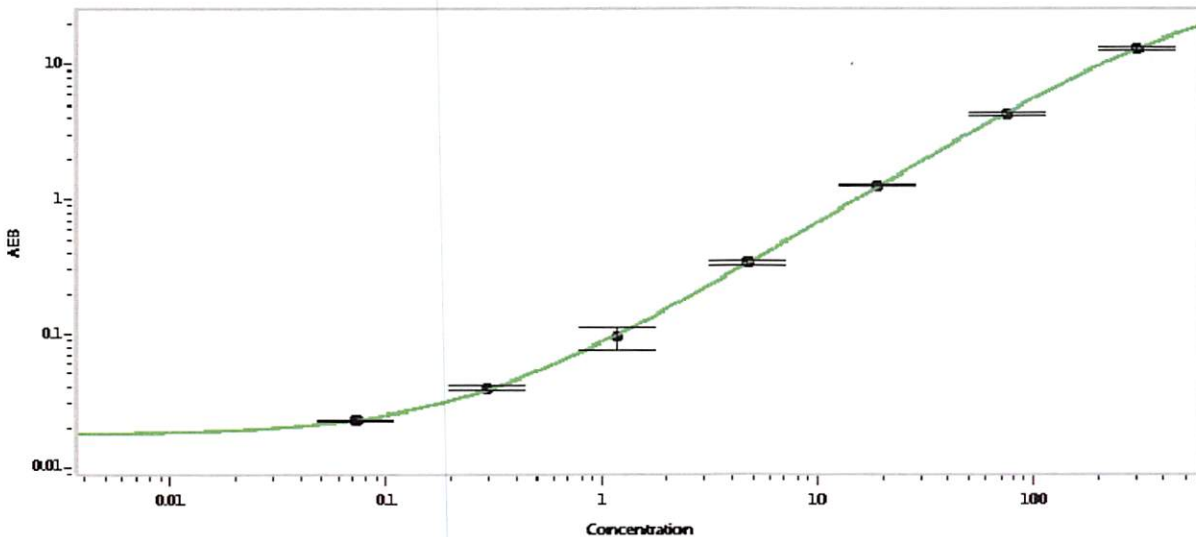
† Specific for use on the SRX platform - refer to the Simoa™ IL-8 SR-X Kit Instructions (KI-0092).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-091217-02	8.73	7.70-11.6
Control 2	MF-091217-02	159	143-220
Panel 1	MF-061617-01	4.61	3.17-6.91
Panel 2	MF-061617-02	51.3	42.6-63.9
Panel 3	MF-061617-03	313	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)	5,293
---	-------

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

Patrizia Stadler Assoc. Director, QC
Name Title

P. Stadler 23/10/2020
Signature/Date