

Product Number:	100198
Lot Number:	502103
Expiration:	29-Oct-2020
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	100819	931819
Detector Reagent	100816	930807
SBG Reagent	100817	930805
Sample Diluent	100818	930804
Calibrator Diluent	100829	930806
Calibrator Concentrate	100828	928804
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

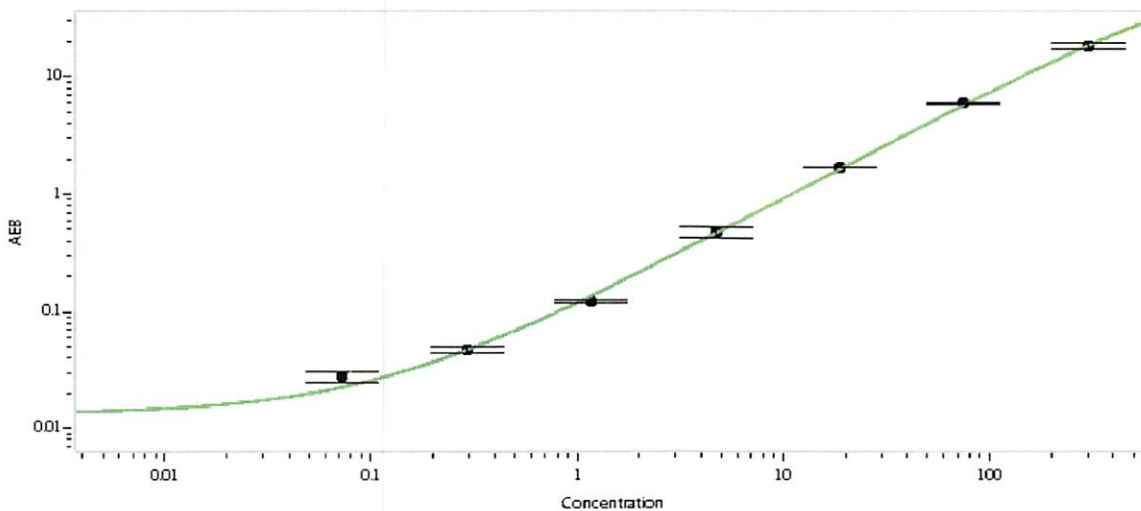
† Specific for use on the SRX platform - refer to the Simoa™ IL-8 SR-X Kit Instructions (KI-0092).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-091217-02	11.5	7.70-11.6
Control 2	MF-091217-02	205	143-220
Panel 1	MF-061617-01	6.02	3.17-6.91
Panel 2	MF-061617-02	53.8	42.6-63.9
Panel 3	MF-061617-03	298	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)	4,574
---	-------

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	<i>P. Stadler</i>	25 NOV 2019
	Name	Title	Signature/Date	