

Product Number:	100198
Lot Number:	502066
Expiration:	9-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	100815	928202
Detector Reagent	100816	928306
SBG Reagent	100817	928305
Sample Diluent	103204	928304
Calibrator Diluent	100829	928206
Calibrator Concentrate	100828	928804
SBG Diluent (SR-X) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

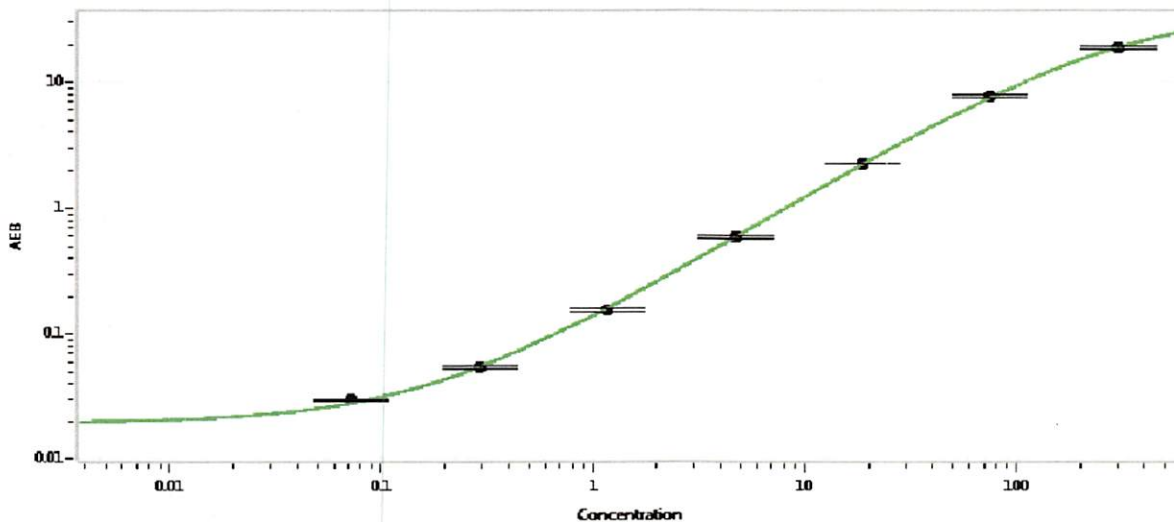
† Specific for use on the SRX platform - refer to the Simoa™ IL-8 SR-X Kit Instructions (KI-0092).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-091217-02	10.7	7.70-11.6
Control 2	MF-091217-02	196	143-220
Panel 1	MF-061617-01	5.92	3.17-6.91
Panel 2	MF-061617-02	55.1	42.6-63.9
Panel 3	MF-061617-03	332	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>	07 NOV 2019
	Name	Title	Signature/Date	