

Product Number:	100198
Lot Number:	501708
Expiration:	25-Feb-2020
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	100819	905705
Detector Reagent	100816	905606
SBG Reagent	100817	906406
Sample Diluent	103204	905605
Calibrator Diluent	100829	905608
Calibrator Concentrate	100828	905708
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

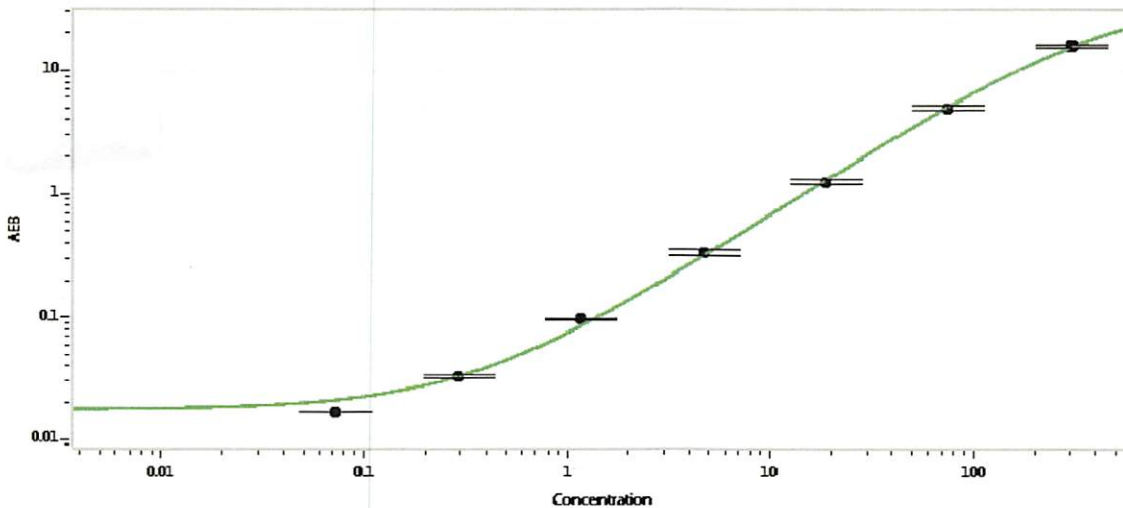
¹ Reagents are not Kit Lot Specific

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

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	MF-091217-02	10.3	7.70-11.6	
Control 2	MF-091217-02	198	143-220	
Panel 1	MF-061617-01	6.47	3.17-6.91	
Panel 2	MF-061617-02	63.1	42.2-63.3	
Panel 3	MF-061617-03	380	267-459	

Calibrator Concentrate Value Assignment (pg/mL)	20,638
---	--------

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 04 Mar 2019
Signature/Date