

Product Number:	100198
Lot Number:	501864
Expiration:	13-Jun-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	916806
Detector Reagent	100816	916813
SBG Reagent	100817	916812
Sample Diluent	103204	916811
Calibrator Diluent	100829	916810
Calibrator Concentrate	100828	905708
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

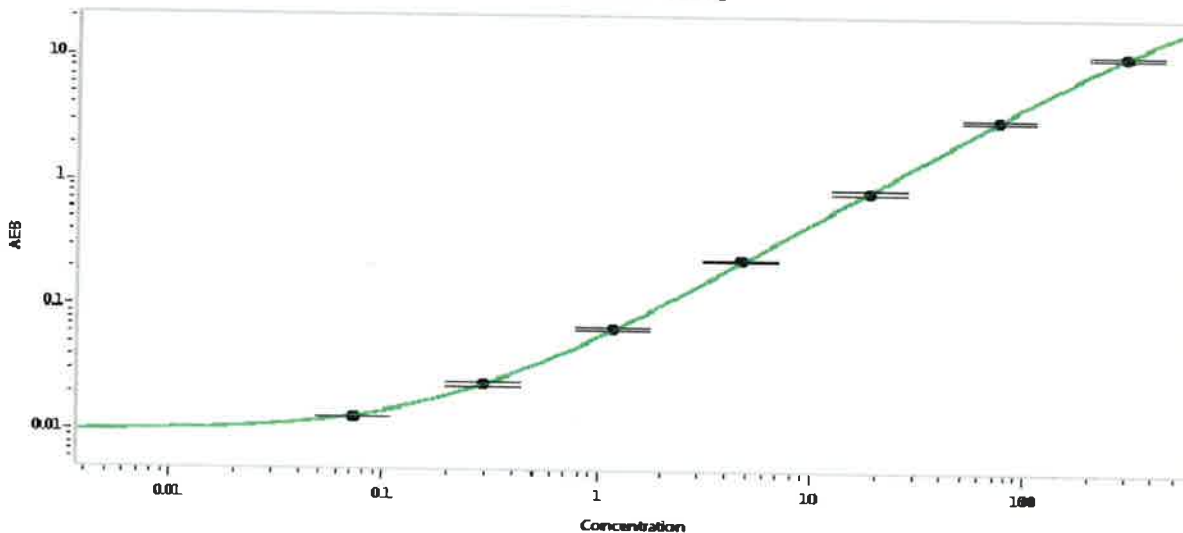
† Specific for use on the SRX platform - refer to the Simoa™ IL-8 SR-X Kit Instructions (KI-0092).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-091217-02	10.5	7.70-11.6
Control 2	MF-091217-02	195	143-220
Panel 1	MF-061617-01	5.96	3.17-6.91
Panel 2	MF-061617-02	63.6	42.6-63.9
Panel 3	MF-061617-03	388	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)	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	<i>P. Stadler</i>	26 JUN 2019
	Name	Title	Signature/Date	