

Product Number:	100198
Lot Number:	502874
Expiration:	21-Mar-2022
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	110321
Detector Reagent	100816	110322
SBG Reagent	100817	110323
Sample Diluent	103204	110328
Calibrator Diluent	100829	110329
Calibrator Concentrate	100828	110507
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

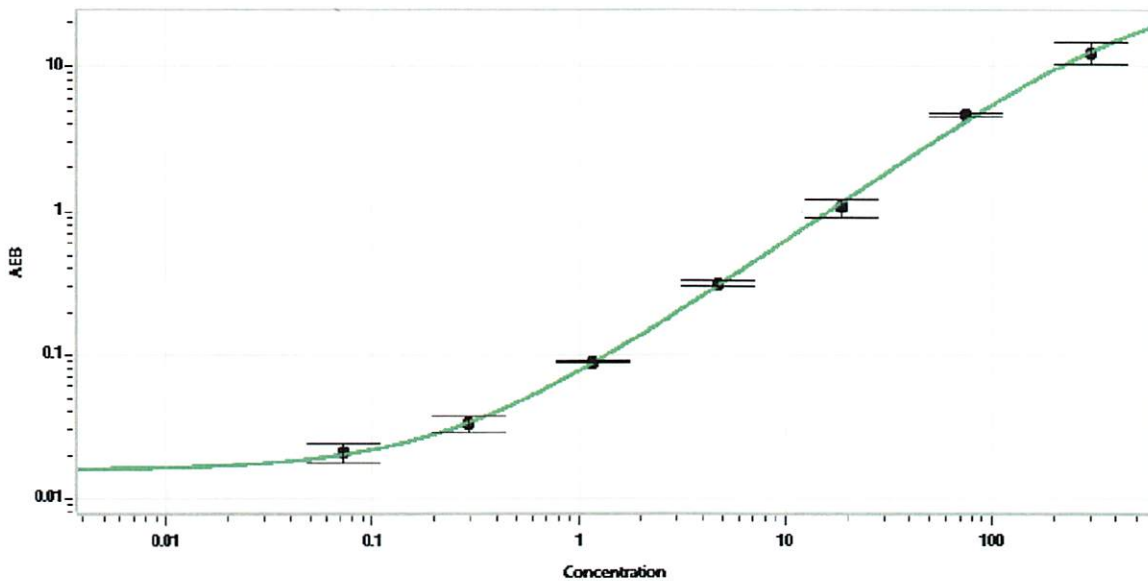
† Specific for use on the SRX platform - refer to the Simoa™ IL-8 SR-X Kit Instructions (KI-0092).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	110507	14.5	11.3-18.8
Control 2	110507	258	188-313
Panel 1	032901	5.69	3.17-6.50
Panel 2	032901	45.9	27.2-56.5
Panel 3	033005	504	358-660

\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	5,753
---	-------

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: Junpil Park, QC Supervisor, Signature: [Signature], Date: 13 May 2021