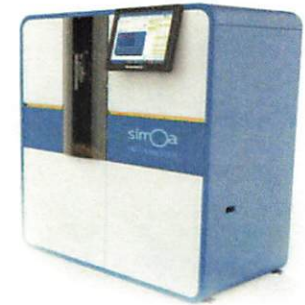


Product Number:	100198
Lot Number:	503159
Expiration:	15-Apr-2022
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	100815	114803
Detector Reagent	100816	114810
SBG Reagent	100817	115316
Sample Diluent	103204	114808
Calibrator Diluent	100829	114811
Calibrator Concentrate	100828	110507
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

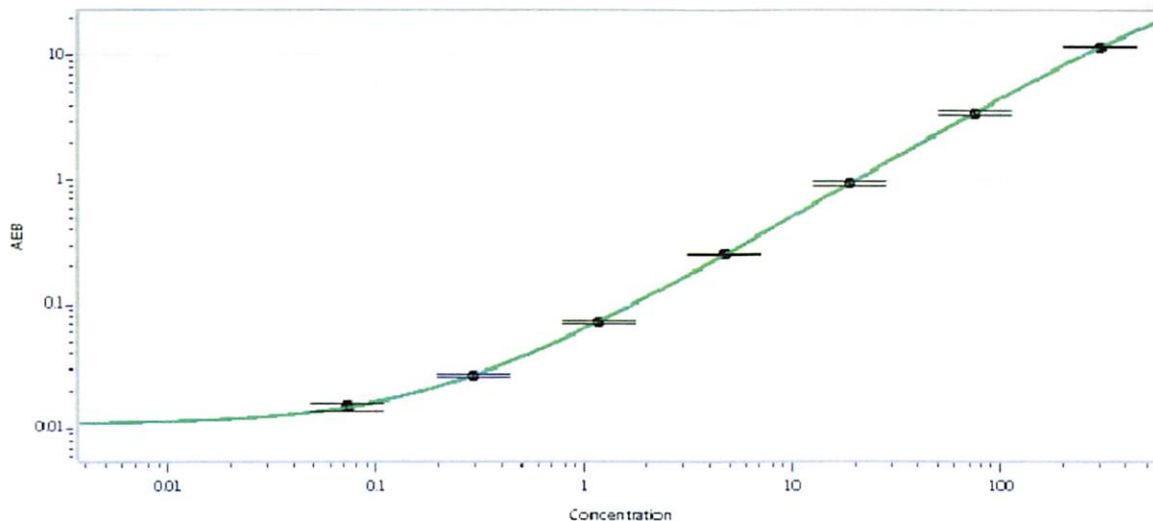
† Specific for use on the SRX platform - refer to the Simoa™ IL-8 SR-X Kit Instructions (KI-0092).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	110507	13.9	11.3-18.8
Control 2	110507	238	188-313
Panel 1	032901	5.23	3.34-6.49
Panel 2	032901	48.4	28.5-57.9
Panel 3	033005	543	363-670

* 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

Shivam Goel
Name

QC

Title

Shivam Goel 27 Sep 2021
Signature/Date