

Product Number:	100198
Lot Number:	503377
Expiration:	3-Dec-2022
Platform(s):	HD-1/HD-X & SR-X



Data below represents results generated on the Simoa® HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	100815	225805
Detector Reagent	100816	212701
SBG Reagent	100817	225901
Sample Diluent	103204	225902
Calibrator Diluent	100829	225903
Calibrator Concentrate	100828	133714
SBG Diluent (SR-X) †	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> These components are not Kit Lot Specific

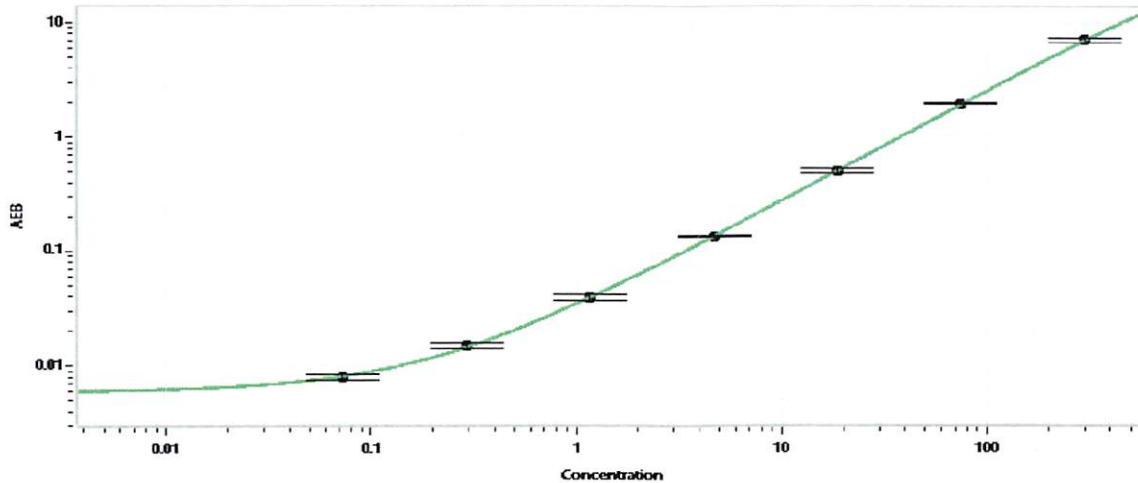
† Specific for use on the SRX platform - refer to the Simoa® IL-8 SR-X Kit Instructions (KI-0092).

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	133714	15.7	11.3-18.8
Control 2	133714	244	188-313
Panel 1	032901	4.15	2.17-6.98
Panel 2	032901	39.5	24.2-58.4
Panel 3	033005	459	355-671

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

Calibrator Concentrate Value Assignment (pg/mL)	4,871
---	-------

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: Sonia Khillan Name, QC Title, Sonia Khillan 12/07/2022 Signature/Date