

Product Number:	100198
Lot Number:	503243
Expiration:	10-Nov-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	133520
Detector Reagent	100816	133701
SBG Reagent	100817	133702
Sample Diluent	103204	133703
Calibrator Diluent	100829	133704
Calibrator Concentrate	100828	133714
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ These components are not Kit Lot Specific

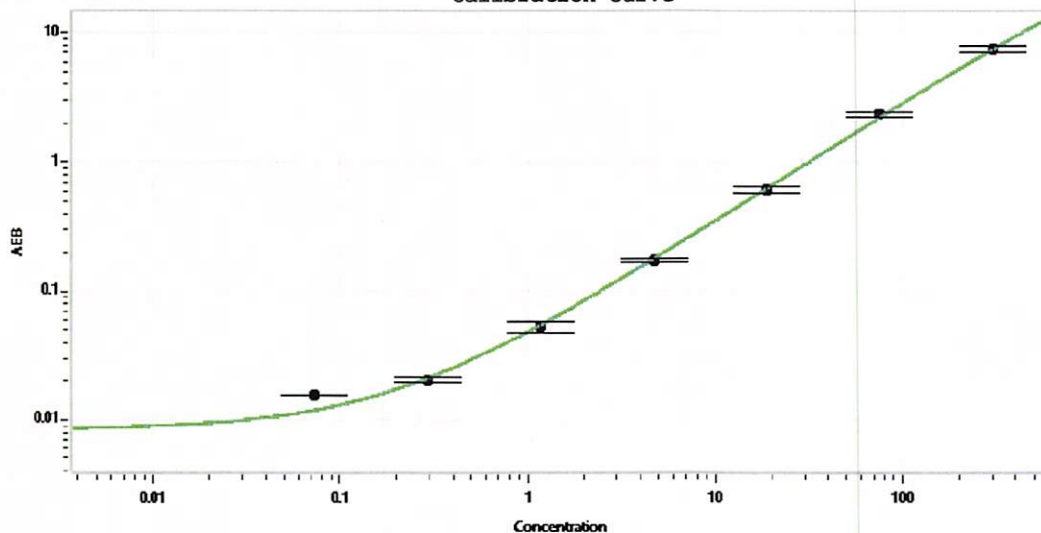
† Specific for use on the SRX platform - refer to the Simoa™ IL-8 SR-X Kit Instructions (KI-0092).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	115807	7.07	5.79 - 11.3
Control 2	115807	139	102 - 191
Panel 1	032901	2.87	2.17 - 6.98
Panel 2	032901	32.1	24.2 - 58.4
Panel 3	033005	497	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)	4,929
---	-------

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

Wenka Tsai

QC Supervisor

06 Jan 2022

Name

Title

Signature/Date