

Product Number:	100794
Lot Number:	503512
Expiration:	1-Jul-2023
Platform(s):	HD-1/HD-X & SR-X



Data below represents results generated on the Simoa® HD-1 / HD-X & SR-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	100781	218203
Detector Reagent	100782	218808
SBG Reagent	100783	218809
Sample Diluent	103202	227229
Calibrator Diluent	100793	231917
Calibrator Concentrate	100780	201216
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

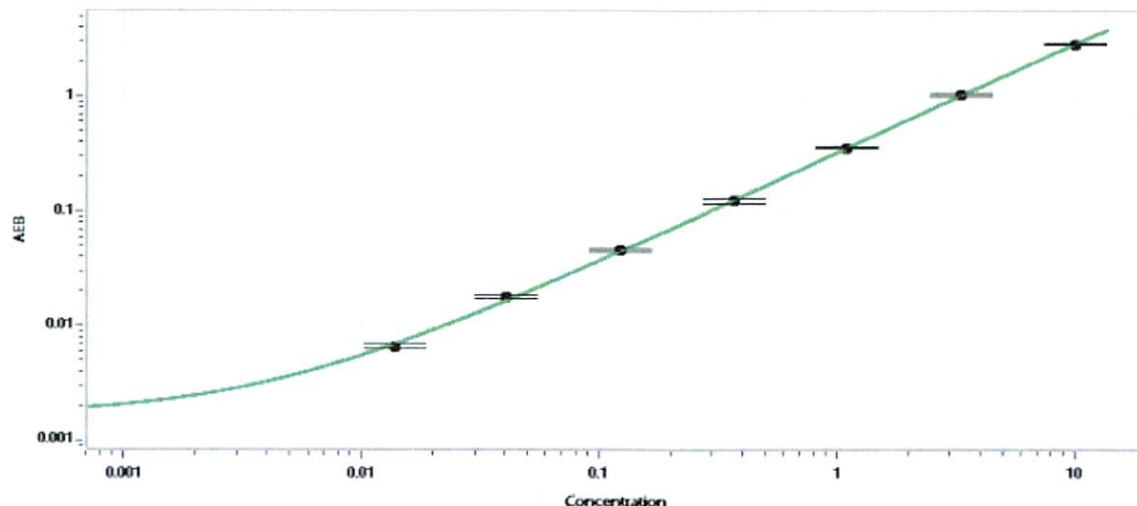
† Specific to use on the SR-X platform - refer to the SR-X Kit Instructions (KI-0085).

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	201216	0.467	0.375 - 0.625
Control 2	201216	5.90	4.50 - 7.50
Panel 1	100801	3.00	1.87 - 4.07
Panel 2	100801	20.3	12.5 - 25.0
Panel 3	100801	2.31	1.12 - 3.15

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

Calibrator Concentrate Value Assignment (pg/mL)	159
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.014
C	0.041
D	0.123
E	0.370
F	1.11
G	3.33
H	10.0

Review/Approval

Sonia Khillan
Name

QC

Title

Soni Kolla / 07MAR2023

Signature/Date