

Product Number:	100794
Lot Number:	503277
Expiration:	30-Jul-2022
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	135202
Detector Reagent	100782	135206
SBG Reagent	100783	135208
Sample Diluent	103202	121121
Calibrator Diluent	100793	134922
Calibrator Concentrate	100780	201216
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

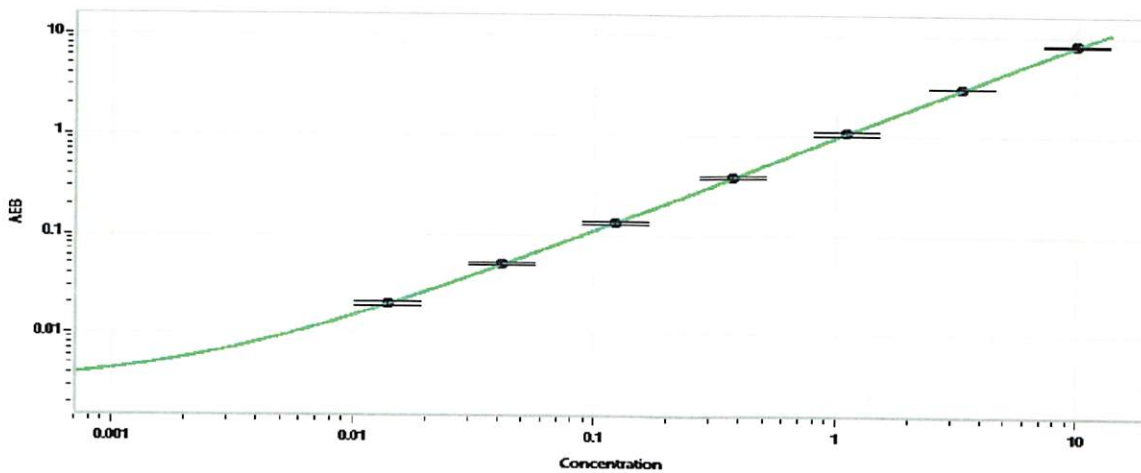
† Specific to use on the SR-X platform - refer to the SR-X Kit Instructions (KI-0085).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	203703	0.348	0.295-0.442
Control 2	203703	4.31	3.44-5.15
Panel 1	100801	2.73	2.32-3.49
Panel 2	100801	17.5	14.3-21.5
Panel 3	100801	1.91	1.71-2.56

* 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

Shivani Goel

QC Manager

Shivani Goel 28 Mar 2022

Name

Title

Signature/Date