

Product Number:	100794
Lot Number:	503916
Expiration:	23-Mar-2024
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	100781	318104	318104A
Detector Reagent	100782	318601	318601A
SBG Reagent	100783	318602	318602A
Sample Diluent	103202	308222	N/A
Calibrator Diluent	100793	313511	313511A
Calibrator Concentrate	100780	319205	319205A
SBG Diluent (SRX) †	103158	N/A ¹	N/A ¹
RGP Reagent	103159	N/A ¹	N/A ¹

¹ RGP is not Kit Lot Specific

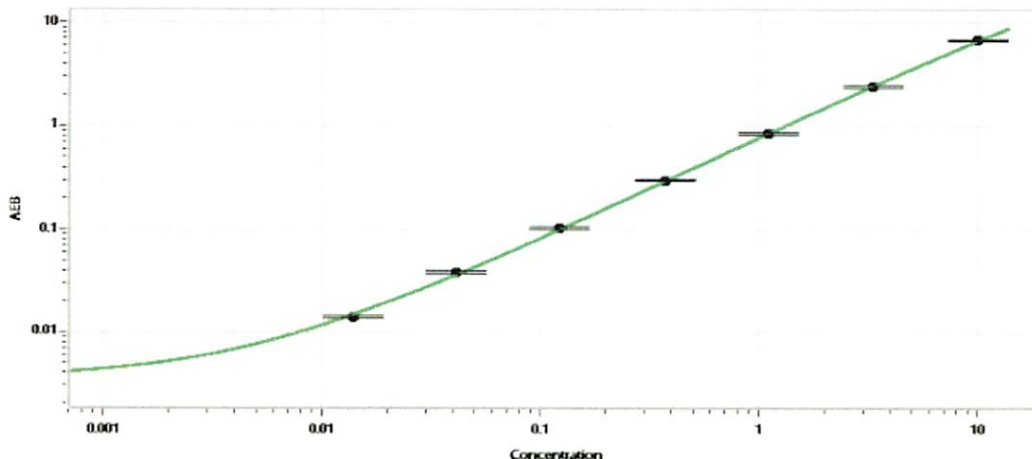
† Specific to use on the SR-X platform - refer to the SR-X Kit Instructions (KI-0085).

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	203703	0.310	0.281 - 0.441
Control 2	203703	3.81	3.062 - 5.706
Panel 1	100801	2.62	1.87 - 4.07
Panel 2	100801	15.4	12.5 - 25.0
Panel 3	100801	1.85	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)	1031
---	------

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 Khillan 20 JUL 2023
Signature/Date