

Product Number:	101599
Lot Number:	502580
Expiration:	17-Apr-2021
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101601	028704
Detector Reagent	102012	028902
SBG Reagent	101602	028712
Sample Diluent	101603	028802
Calibrator Diluent	102014	010815
Calibrator Concentrate	100778	001414
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



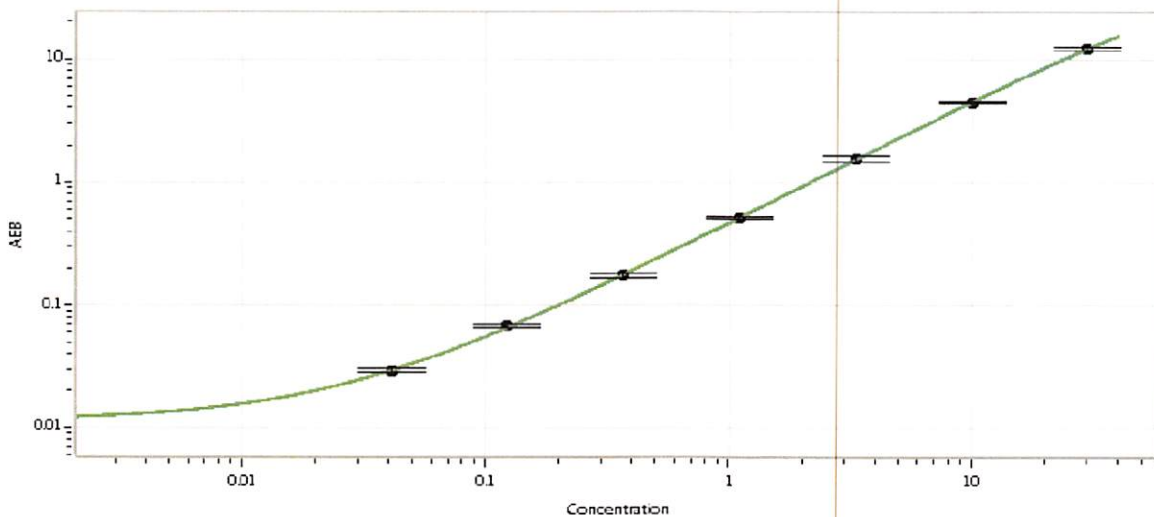
Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-012518-02	0.430	0.352-0.537
Control 2	MF-012518-02	22.2	14.5-28.3
Panel 3	900903	8.34	6.37-10.7
Panel 4	900904	31.3	21.1-38.7

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

Calibrator Concentrate Value Assignment (pg/mL)	706
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Shivani Goel

QC Manager

Shwam Goel 02Nov2020

Name

Title

Signature/Date