

Product Number:	101622
Lot Number:	502359
Expiration:	21-Feb-2021
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101624	006904
Detector Reagent	101623	007203
SBG Reagent	101625	006509
Sample Diluent	103203	005220
Calibrator Diluent	101627	006516
Calibrator Concentrate	100652	006408
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



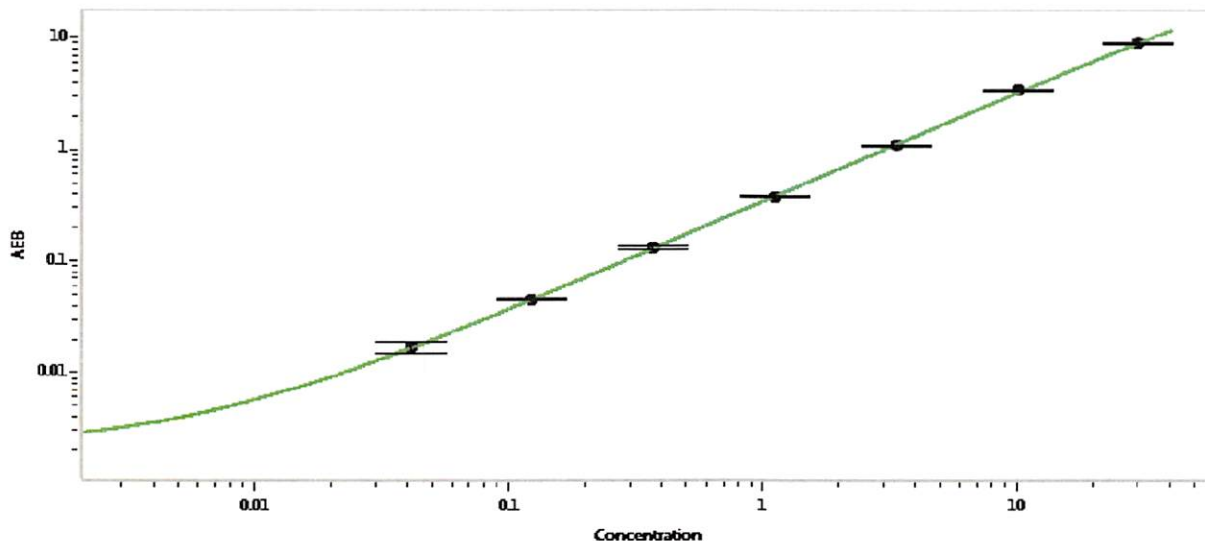
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	901702	1.04	0.944-1.42
Control 2	901702	15.5	15.4-23.1
Panel 1	MF-050918-01	46.9	41.6-67.5
Panel 2	MF-050918-02	6.73	5.03-8.84
Panel 3	MF-050918-03	0.938	0.644-1.24

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

Calibrator Concentrate Value Assignment (pg/mL)	675
---	-----

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

Shivani Goel 11 May 2020

Name

Title

Signature/Date