

Product Number:	101622
Lot Number:	502483
Expiration:	23-Jun-2021
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101624	021010
Detector Reagent	101623	021105
SBG Reagent	101625	021202
Sample Diluent	103203	021207
Calibrator Diluent	101627	020417
Calibrator Concentrate	100652	021203
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



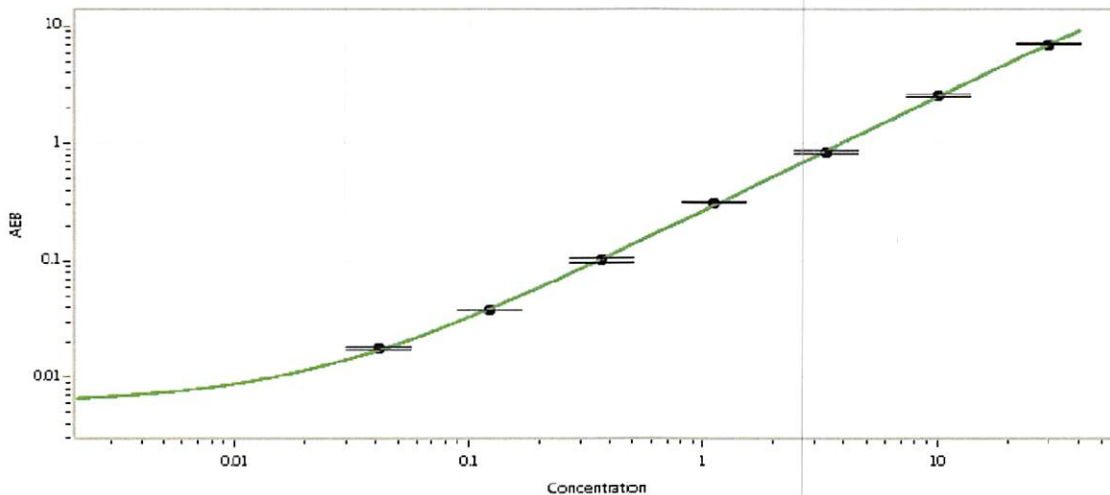
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	901702	1.08	0.944-1.42
Control 2	901702	16.3	15.4-23.1
Panel 1	MF-050918-01	50.2	41.6-67.5
Panel 2	MF-050918-02	6.78	5.03-8.84
Panel 3	MF-050918-03	0.827	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)	606
---	-----

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 13 Aug 2020

Name

Title

Signature/Date