

Product Number:	101635
Lot Number:	502374
Expiration:	29-Apr-2021
Platform(s):	HD-1/ HD-X



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

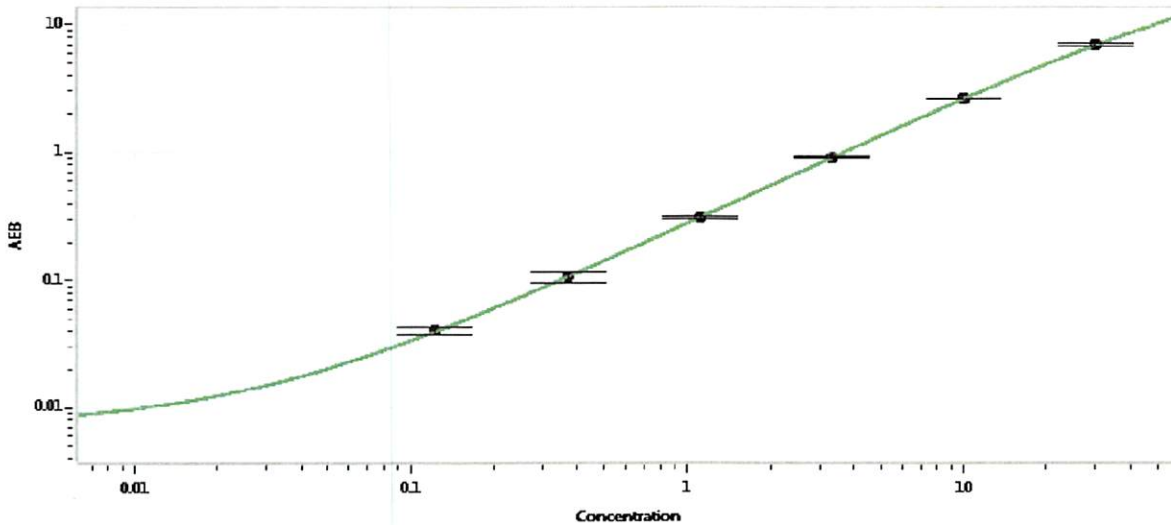
Component	Part Number	Lot Number
Bead Reagent	101637	013101
Detector Reagent	101636	013203
SBG Reagent	101638	013104
Sample Diluent	101639	013202
Calibrator Diluent	101640	013108
Calibrator Concentrate	100661	908402
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	904406	0.656	0.550-0.825	
Control 2	904406	67.9	54.1-81.2	
Panel 1	MF-040618-02	0.638	0.414-0.657	
Panel 2	MF-040618-03	5.50	4.18-6.26	
Panel 3	MF-040618-04	57.1	44.1-66.2	

Calibrator Concentrate Value Assignment (pg/mL)	1760
---	------

Calibration Curve



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

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 21 May 2020
	Name	Title	Signature/Date