

Product Number:	101622
Lot Number:	502414
Expiration:	13-May-2021
Platform(s):	HD-1 / HD-X & SR-X



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

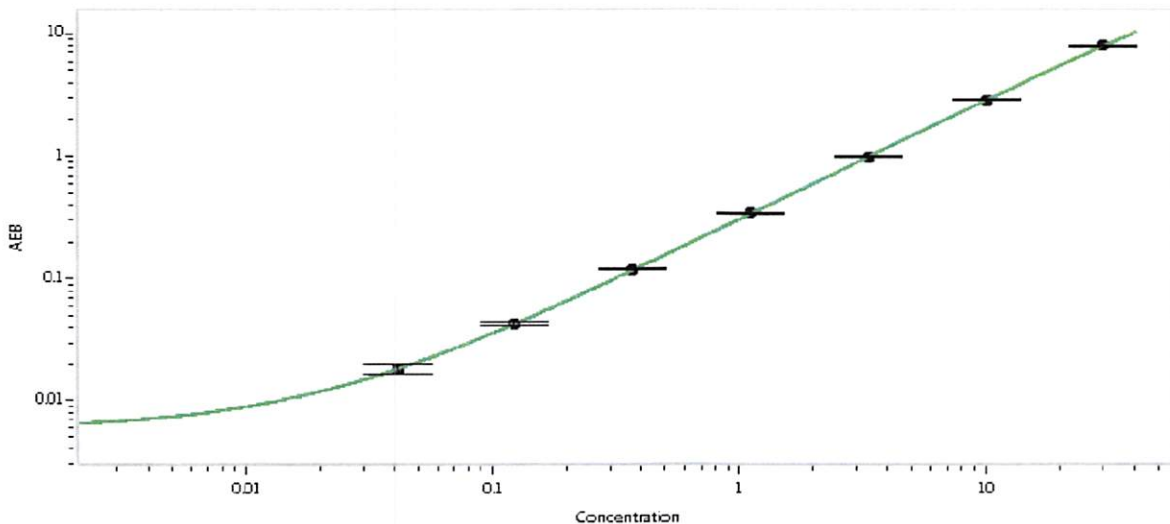
Component	Part Number	Lot Number
Bead Reagent	101624	016408
Detector Reagent	101623	016702
SBG Reagent	101625	015903
Sample Diluent	103203	015413
Calibrator Diluent	101627	015805
Calibrator Concentrate	100652	006408
RGP Reagent	103159	N/A ¹

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	901702	1.13	0.944-1.42	
Control 2	901702	17.3	15.4-23.1	
Panel 1	MF-050918-01	53.5	41.6-67.5	
Panel 2	MF-050918-02	6.97	5.03-8.84	
Panel 3	MF-050918-03	0.930	0.644-1.24	

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 23/6/2020
	Name	Title	Signature/Date