

Product Number:	101664
Lot Number:	503313
Expiration:	15-Jun-2022
Platform(s):	HD-1/HD-X & SR-X



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

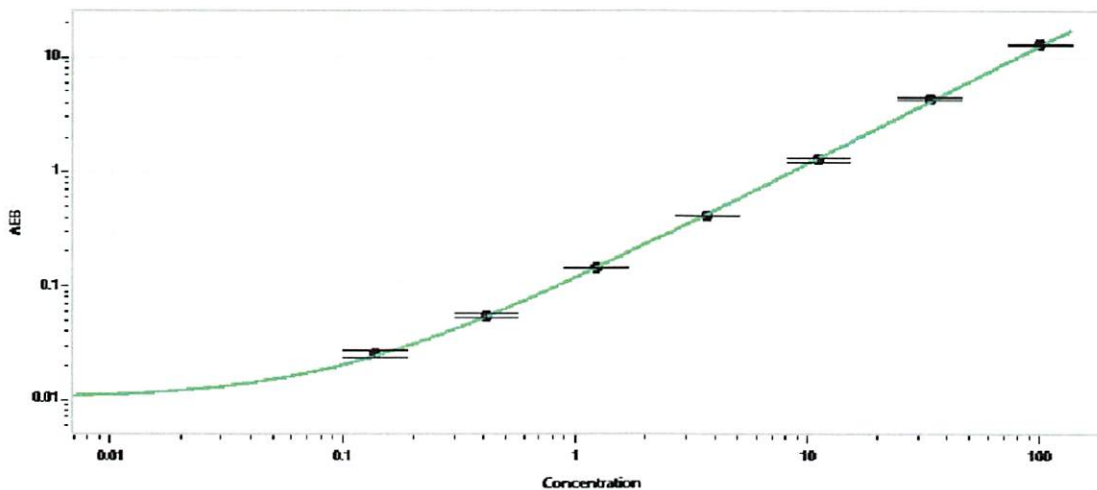
Component	Part Number	Lot Number
Bead Reagent	101666	135004
Detector Reagent	101665	202813
SBG Reagent	101667	202814
Sample Diluent	101668	201810
Calibrator Diluent	101669	203108
Calibrator Concentrate	103217	202410
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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	201802	1.11	0.690-1.36	
Control 2	201802	67.3	46.2-75.1	
Panel 1	202604	7.90	5.87-8.81	
Panel 2	202604	49.3	32.2-54.6	
Panel 3	202604	116	90.4-136	
Panel 4	N/A	N/A	N/A	

Calibrator Concentrate Value Assignment (pg/mL)	1417
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval	Shivani Goel	QC Manager	Shivani Goel	18Feb2022
	Name	Title	Signature/Date	