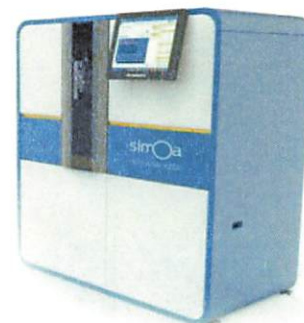


Product Number:	103337
Lot Number:	503298
Expiration:	29-Jun-2022
Platform(s):	HD-1/HD-X & SR-X



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

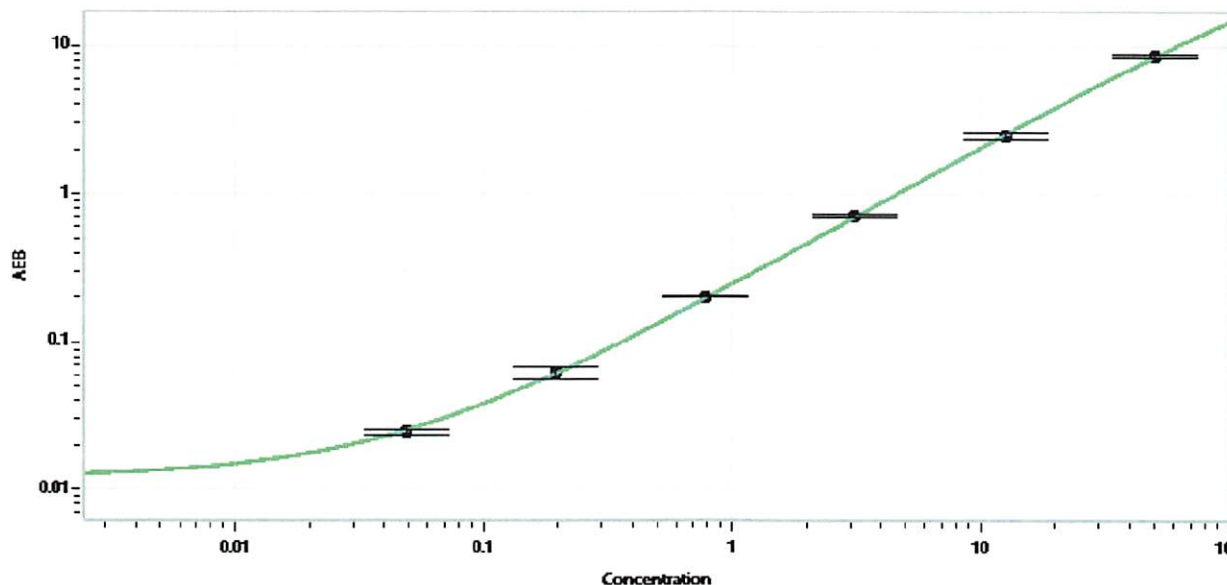
Component	Item Number	Lot Number
Bead Reagent	103334	118210
Detector Reagent	101091	118402
SBG Reagent	101101	201209
Sample Diluent	103336	118403
Calibrator Diluent	101095	118020
Calibrator Concentrate	101093	118213
RGP Reagent	103159	N/A ¹

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	118213	0.789	0.750-1.25	
Control 2	118213	16.2	15.0-25.0	
Panel 2	024004	3.19	2.26-4.95	
Panel 3	024004	45.31	32.5-57.3	

Calibrator Concentrate Value Assignment (pg/mL)	1,895
---	-------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.049
C	0.195
D	0.781
E	3.13
F	12.5
G	50.0

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 28 Jan 2022

Name

Title

Signature/Date