

Product Number:	103186
Lot Number:	503196
Expiration:	28-Jul-2022
Platform(s):	HD-1/ HD-X

Component	Part Number	Lot Number
Bead Reagent	102246	121015
Detector Reagent	102248	121028
SBG Reagent	102250	121030
Sample Diluent	102252	121031
Calibrators	102255	129303
Control 1	102256	129304
Control 2	102257	129304
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

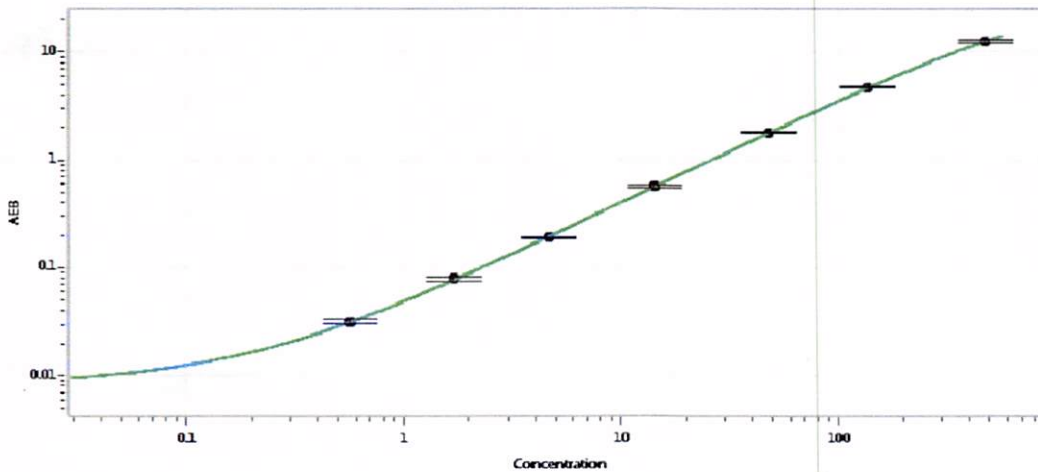


Data below represents results generated on the Simoa HD-1® Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	129304	4.40	3.34-5.02
Control 2	129304	288	179-398
Panel 1	008501	12.7	8.59-20.5
Panel 2	008501	10.0	7.39-15.1
Panel 3	008501	60.8	43.8-91.8
Panel 4	009401	199	128-302

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.569
C	1.71
D	4.61
E	14.2
F	47.8
G	137
H	473

Review/Approval: Junpil Park, QC Supervisor, Signature: [Signature], Date: 26 Oct 2021