

Product Number:	101132
Lot Number:	502998
Expiration:	28-Dec-2021
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	101119	116403
Detector Reagent	101121	105305
SBG Reagent	101131	106501
Sample Diluent	103319	036320
Calibrator Diluent	101125	105013
Calibrator Concentrate	101123	105511
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



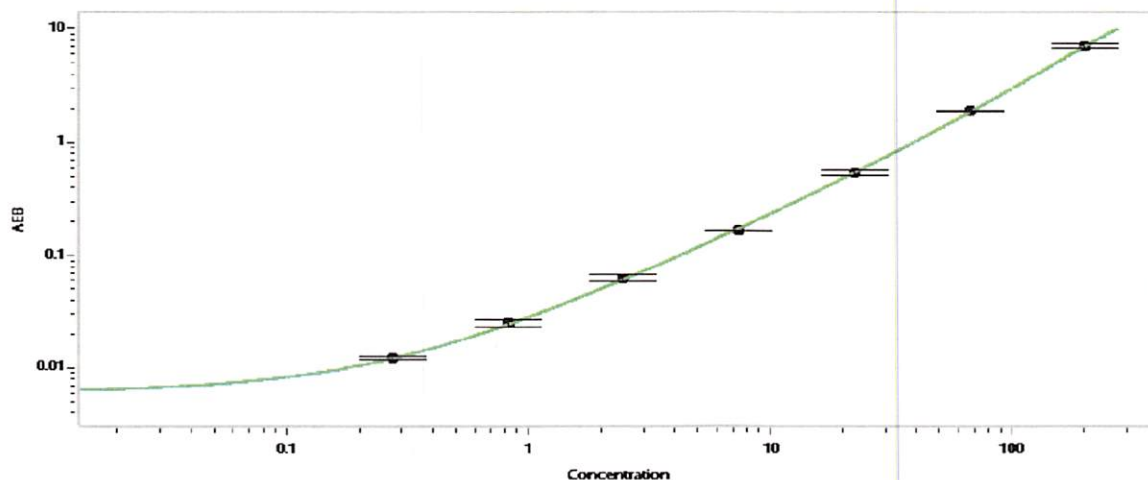
Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920702	2.80	1.68-3.11
Control 2	920702	288	122-349
Panel 1	106201	153	95.3-224
Panel 2	106201	81.2	66.4-121
Panel 3	106201	286	219-328
Panel 4	106201	251	195-292

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

Calibrator Concentrate Value Assignment (pg/mL)	9021
---	------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Junpil Park
Name

QC Supervisor
Title

 24 Jun 2021
Signature/Date