

Product Number:	103186
Lot Number:	503019
Expiration:	27-Jun-2022
Platform(s):	HD-1/ HD-X

Component	Part Number	Lot Number
Bead Reagent	102246	117801
Detector Reagent	102248	117803
SBG Reagent	102250	117808
Sample Diluent	102252	117914
Calibrators	102255	118706
Control 1	102256	118707
Control 2	102257	118707
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

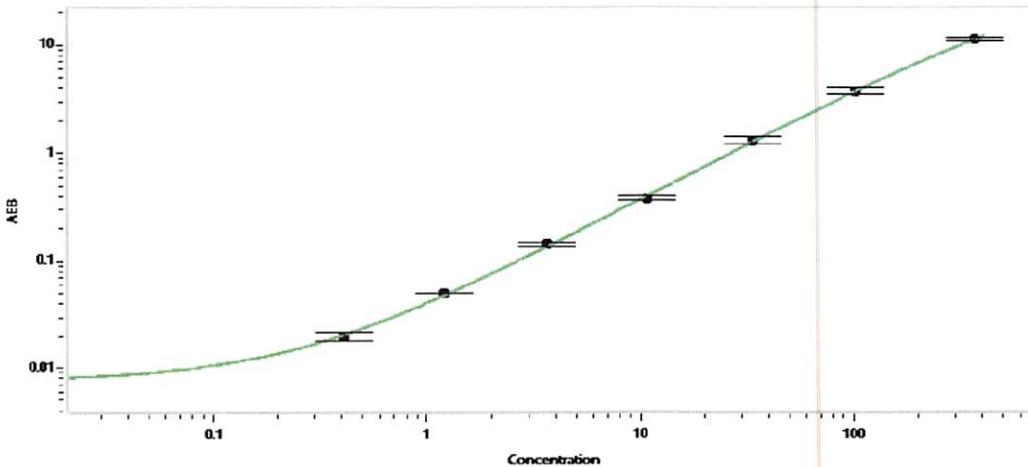


Data below represents results generated on the Simoa HD-X® Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	118707	3.43	2.62-3.93
Control 2	118707	186	138-214
Panel 1	008501	14.4	8.59-20.5
Panel 2	008501	11.4	7.39-15.1
Panel 3	008501	61.4	43.8-91.8
Panel 4	009401	189	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	0.410
C	1.21
D	3.67
E	10.6
F	33.6
G	102
H	369

Review/Approval

Junpil Park
Name

QC Supervisor
Title

Signature/Date

 19 Jul 2021