

Product Number:	103186
Lot Number:	502625
Expiration:	24-Sep-2021
Platform(s):	HD-1/ HD-X

Component	Part Number	Lot Number
Bead Reagent	102246	026804
Detector Reagent	102248	026904
SBG Reagent	102250	026908
Sample Diluent	102252	026905
Calibrators	102255	030009
Control 1	102256	030201
Control 2	102257	030201
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

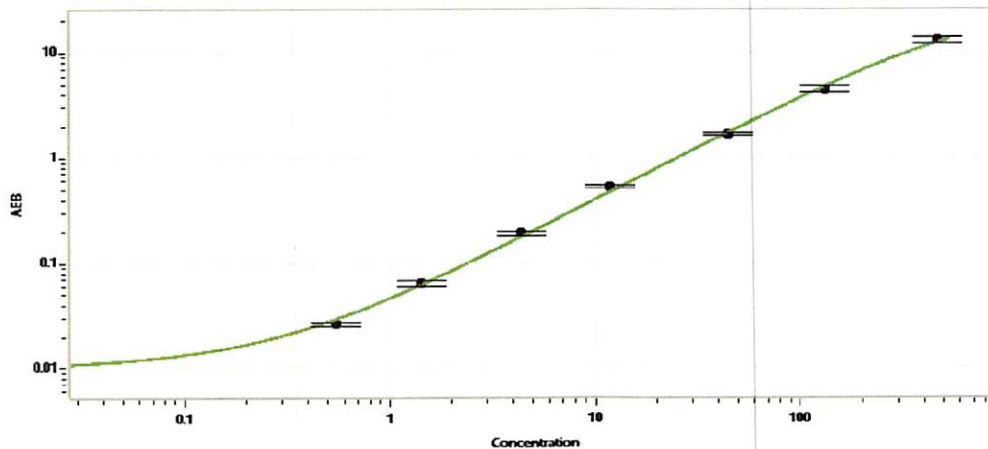


Data below represents results generated on the Simoa HD-X® Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	030201	4.08	2.93 - 5.21
Control 2	030201	159	120 - 203
Panel 1	922601	19.0	14.3 - 21.4
Panel 2	922602	6.67	4.97 - 7.45
Panel 3	922603	40.8	28.8 - 43.2
Panel 4	922604	175	133 - 201

* 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	0.544
C	1.43
D	4.41
E	11.8
F	44.5
G	133
H	471

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 04 Dec 2020
Signature/Date