

Product Number:	103186
Lot Number:	502607
Expiration:	14-Oct-2021
Platform(s):	HD-1/ HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	102246	028804
Detector Reagent	102248	029003
SBG Reagent	102250	031118
Sample Diluent	102252	029005
Calibrators	102255	030009
Control 1	102256	030201
Control 2	102257	030201
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

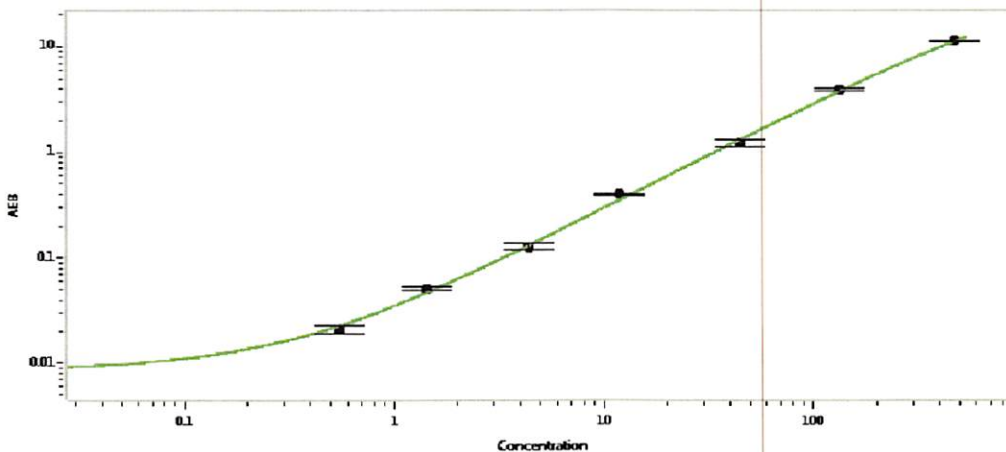


Data below represents results generated on the Simoa HD-1® Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	030201	4.46	2.84-5.30
Control 2	030201	153	118-207
Panel 1	926601	17.3	14.3-21.4
Panel 2	922602	6.22	4.97-7.45
Panel 3	922603	38.7	28.8-43.2
Panel 4	922604	156	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: Shivani Goel QC Manager
Name Title

Shivani Goel 13 Nov 2020
Signature/Date