

Product Number:	103186
Lot Number:	502323
Expiration:	13-Mar-2021
Platform(s):	HD-1/HD-x

Component	Part Number	Lot Number
Bead Reagent	102246	010001
Detector Reagent	102248	009701
SBG Reagent	102250	009707
Sample Diluent	102252	009702
Calibrators	102255	007001
Controls	102256 - 102257	007002
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

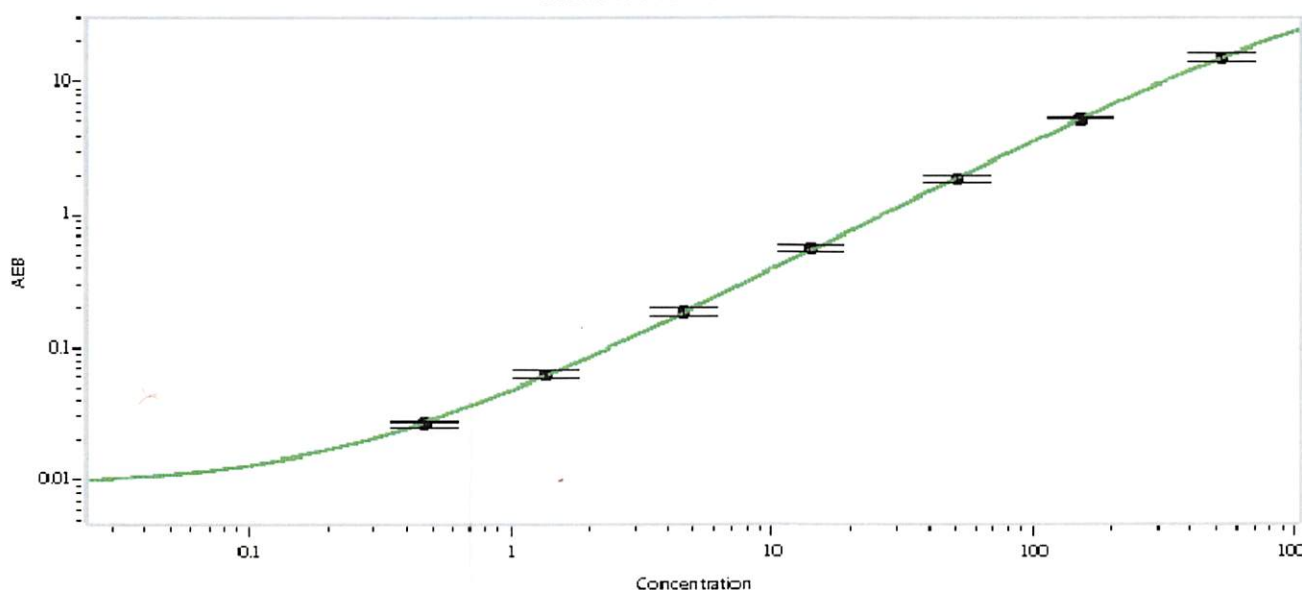


Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	007002	4.17	2.52-5.10
Control 2	007002	173	124-186
Panel 1	922601	18.8	14.1-21.1
Panel 2	922602	6.45	4.94-7.41
Panel 3	922603	41.0	28.4-42.6
Panel 4	922604	148	129-204

* 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	0.474
C	1.36
D	4.58
E	14.1
F	51.4
G	152
H	526

Review/Approval

Shivani Goel

Name

QC Manager

Title

Shivani Goel 16 Apr 2020

Signature/Date