

Product Number:	101132
Lot Number:	502174
Expiration:	26-Jun-2020
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	101119	936103
Detector Reagent	101121	936005
SBG Reagent	101131	936006
Sample Diluent	103319	933810
Calibrator Diluent	101125	932923
Calibrator Concentrate	101123	81771
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



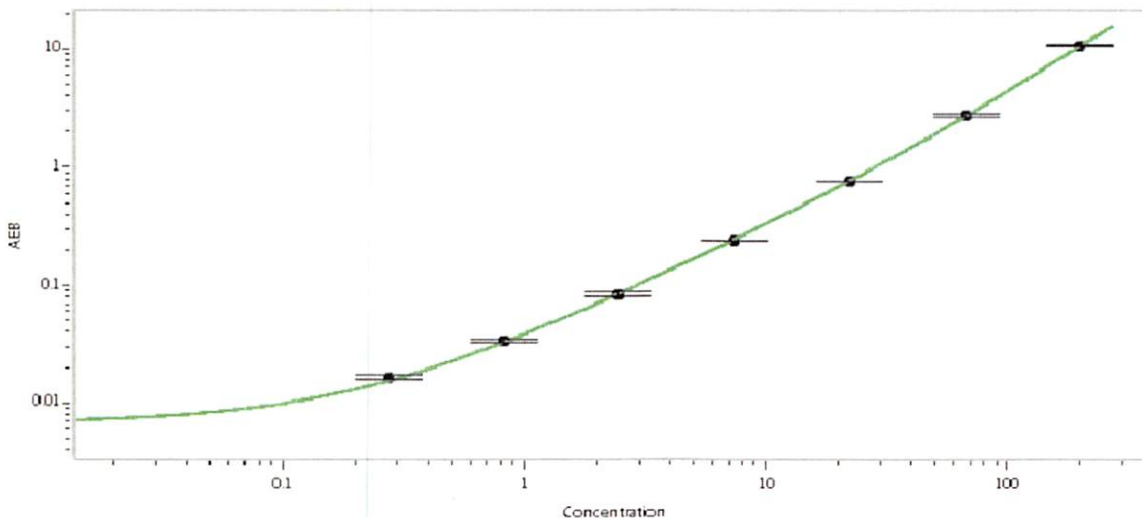
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-051716-04	2.14	1.39-3.39
Control 2	MF-051716-04	249	127-431
Panel 1	830204	100	79.0-119
Panel 2	830205	88.1	69.2-104
Panel 3	830206	246	193-289
Panel 4	830207	315	243-365

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

Calibrator Concentrate Value Assignment (pg/mL)	55360
---	-------

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

Patrizia Stadler Assoc. Director, QC

P. Stadler

13 Jun 2020

Name

Title

Signature/Date