

Product Number:	101552
Lot Number:	502694
Expiration:	25-Nov-2021
Platform(s):	HD-1/HD-x & SR-X



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

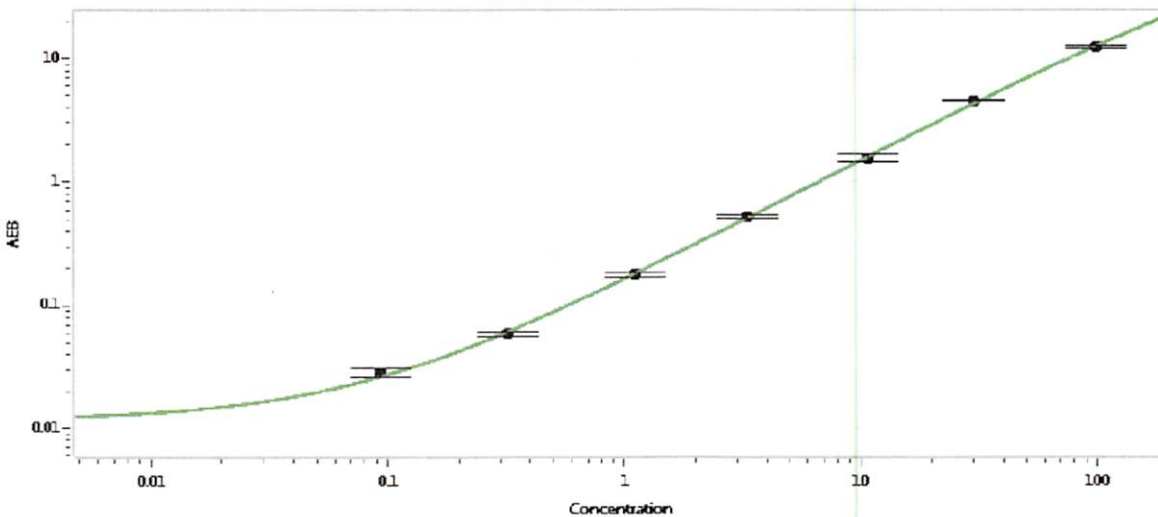
Component	Part Number	Lot Number
Bead Reagent	101554	033006
Detector Reagent	101553	105601
SBG Reagent	101555	033904
Sample Diluent	101556	033905
Calibrators	101559	034301
Control 1	101560	034302
Control 2	101561	034302
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	034302	0.985	0.714-1.22
Control 2	034302	66.1	37.4-76.5
Panel 2	916902	1.26	0.635-1.32
Panel 3	916903	31.1	13.0-36.2
Panel 4	916904	92.1	37.0-118

* 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	0.091
C	0.279
D	0.837
E	2.57
F	8.53
G	26.5
H	78.4

Review/Approval

Junpil Park
Name

QC Supervisor
Title

Signature/Date

25 Mar 2021