

Product Number:	101552
Lot Number:	503136
Expiration:	10-May-2022
Platform(s):	HD-1/HD-x & SR-X



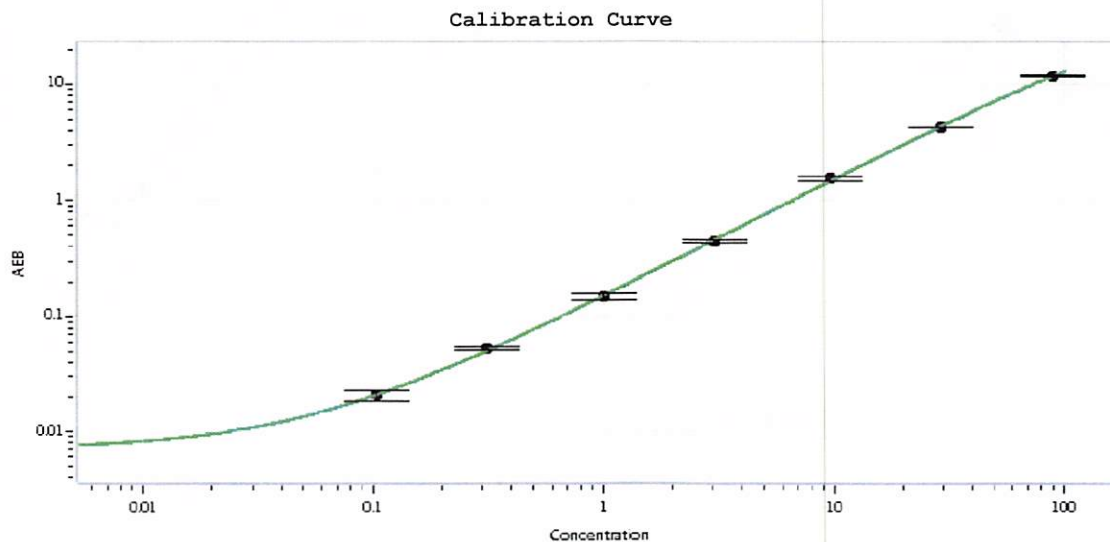
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101554	113902
Detector Reagent	101553	125604
SBG Reagent	101555	114111
Sample Diluent	101556	113019
Calibrators	101559	116005
Control 1	101560	116006
Control 2	101561	116006
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	116006	1.38	0.888 - 1.57
Control 2	116006	57.1	47.1 - 70.6
Panel 1	117205	4.04	3.38 - 5.07
Panel 2	117205	2.56	1.90 - 5.09
Panel 3	117205	24.0	17.7 - 26.5
Panel 4	119612	37.6	22.1 - 40.5

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



Calibrator Levels (pg/mL)

A	0.000
B	0.103
C	0.311
D	1.00
E	3.03
F	9.61
G	28.8
H	89.1

Review/Approval

Junpil Park
Name

QC Supervisor
Title

Signature/Date

23 Sep 2021