

Product Number:	101478
Lot Number:	502878
Expiration:	15-Apr-2022
Platform(s):	HD-1/HD-X & SR-X



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

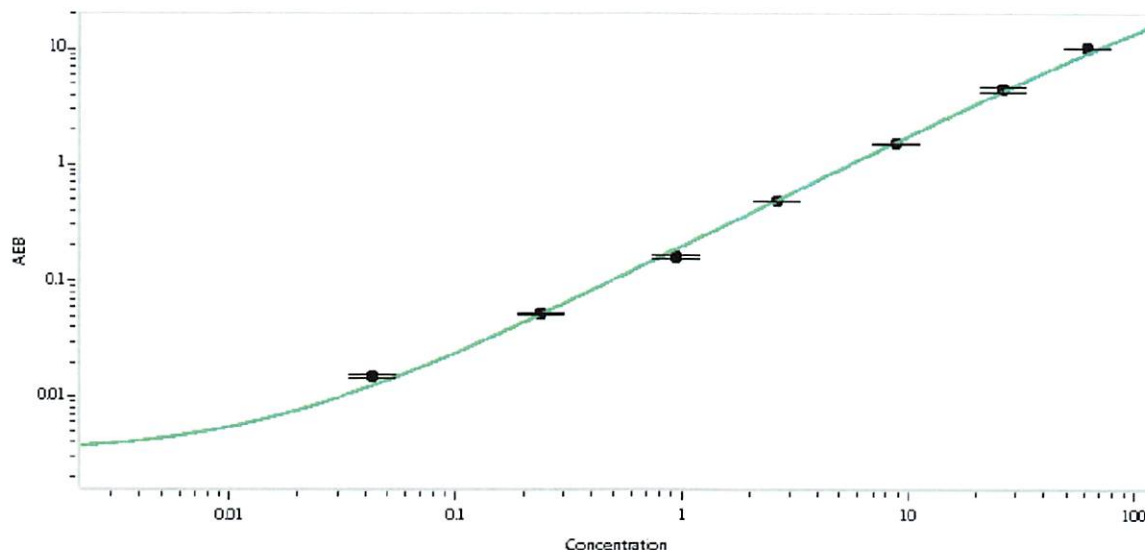
Component	Part Number	Lot Number
Bead Reagent	101480	110511
Detector Reagent	101481	110703
SBG Reagent	101482	110704
Sample Diluent	101483	110705
Calibrators	100758-100765	121603
Control 1	100768	121604
Control 2	100769	121604
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	121604	2.53	1.69 - 3.70
Control 2	121604	54.5	42.7 - 64.0
Panel 1	122912	0.238	0.178 - 0.297
Panel 2	122912	5.16	3.99 - 5.99
Panel 3	122912	50.4	44.6 - 67.0

* 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	0.043
C	0.237
D	0.937
E	2.62
F	8.76
G	26.2
H	61.6

Review/Approval

Junpil Park
Name

QC Supervisor
Title

Signature/Date

02 Sep 2021