

Product Number:	100860
Lot Number:	502675
Expiration:	23-Dec-2021
Platform(s):	HD-1/ HD-X



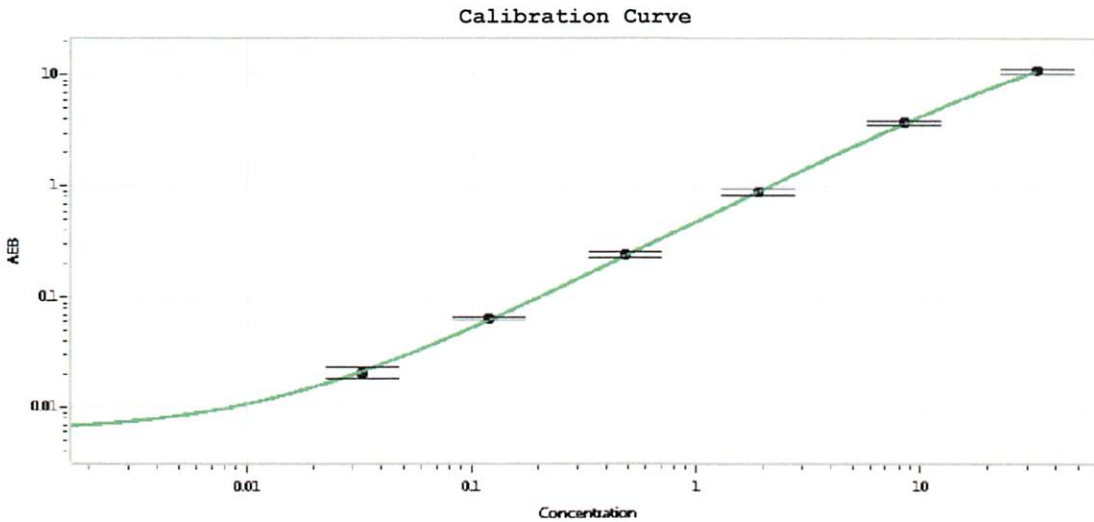
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	100865	100809
Detector Reagent	100862	100903
SBG Reagent	100868	106416
Sample Diluent	100870	035801
Calibrators	101111	101105
Control 1	100875	101106
Control 2	100876	101106
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	101106	6.46	4.94-7.42
Control 2	101106	0.110	0.082-0.123
Panel 1	908501	1.27	0.631-2.10
Panel 2	908502	8.73	4.38-12.5
Panel 3	908503	26.9	12.1-47.2

* 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	0.033
C	0.121
D	0.485
E	1.90
F	8.47
G	32.8

Review/Approval

Junpil Park
Name

QC Supervisor
Title

16 Mar 2021
Signature/Date