

Product Number:	100860
Lot Number:	502831
Expiration:	23-Mar-2022
Platform(s):	HD-1/ HD-X



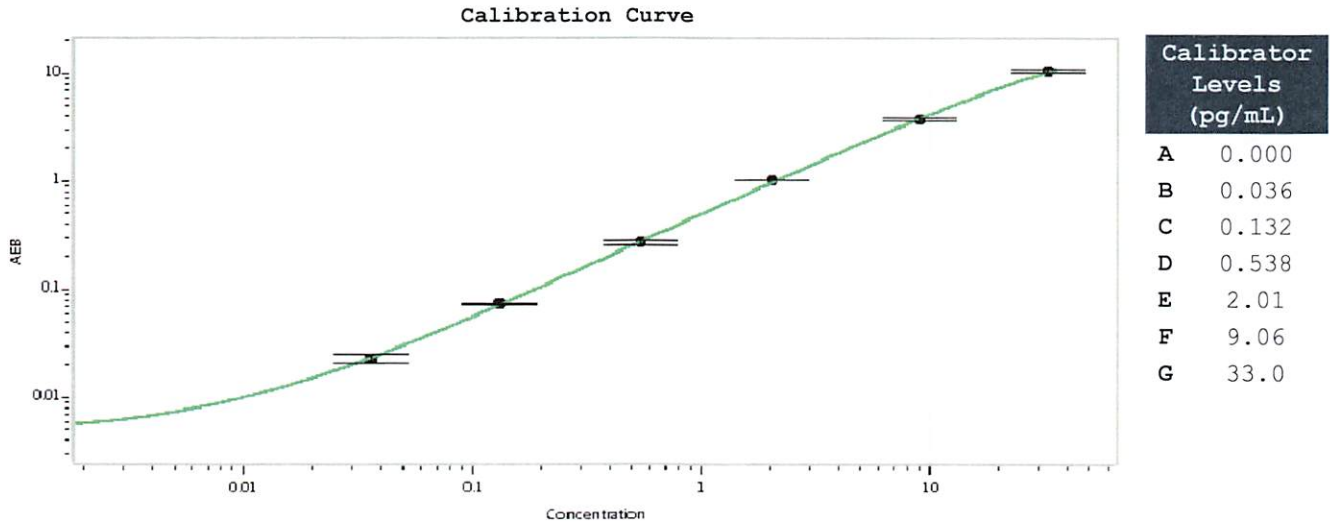
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	100865	108407
Detector Reagent	100862	108511
SBG Reagent	100868	108512
Sample Diluent	100870	108217
Calibrators	101111	108701
Control 1	100875	108702
Control 2	100876	108702
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	108702	6.53	5.47-8.20
Control 2	108702	0.114	0.097-0.146
Panel 1	908501	1.22	0.631-2.10
Panel 2	908502	8.35	4.38-12.5
Panel 3	908503	35.3	12.1-47.2

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



Review/Approval	Junpil Park <hr/> Name	QC Supervisor <hr/> Title	 <hr/> Signature/Date
			15 Apr 2021