

Product Number:	103714
Lot Number:	503008
Expiration:	18-Apr-2022
Platform(s):	HD-1 / HD-X & SR-X

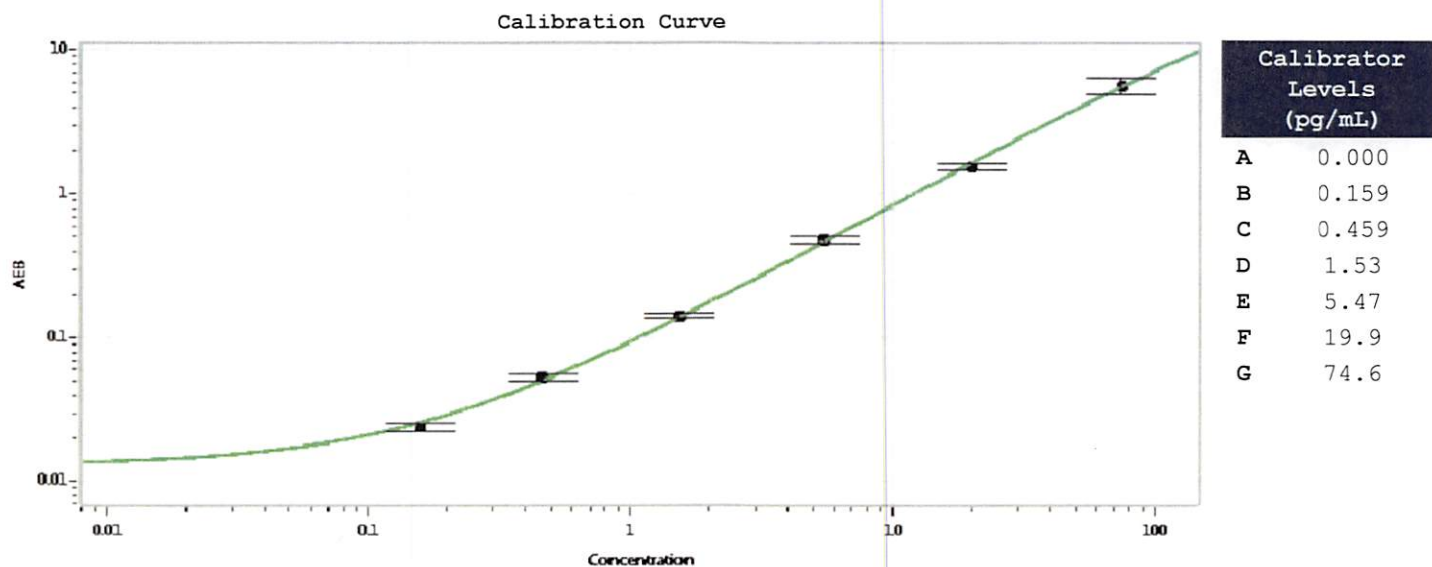
Component	Part Number	Lot Number
Bead Reagent	103698	116904
Detector Reagent	103700	117007
SBG Reagent	103702	116907
Sample Diluent	103704	117113
Reference Calibrator	103711	117303
Control 1	103712	117304
Control 2	103713	117304
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



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

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	117304	3.21	2.48-3.72	
Control 2	117304	124	92.2-138	
Panel 1	013903	0.528	0.258-1.11	
Panel 2	013903	1.92	0.841-2.44	
Panel 3	013903	2.43	1.79-4.31	



Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 08 Jul 2021
Signature/Date