

Product Number:	103714
Lot Number:	502793
Expiration:	8-Sep-2021
Platform(s):	HD-1 / HD-X & SR-X

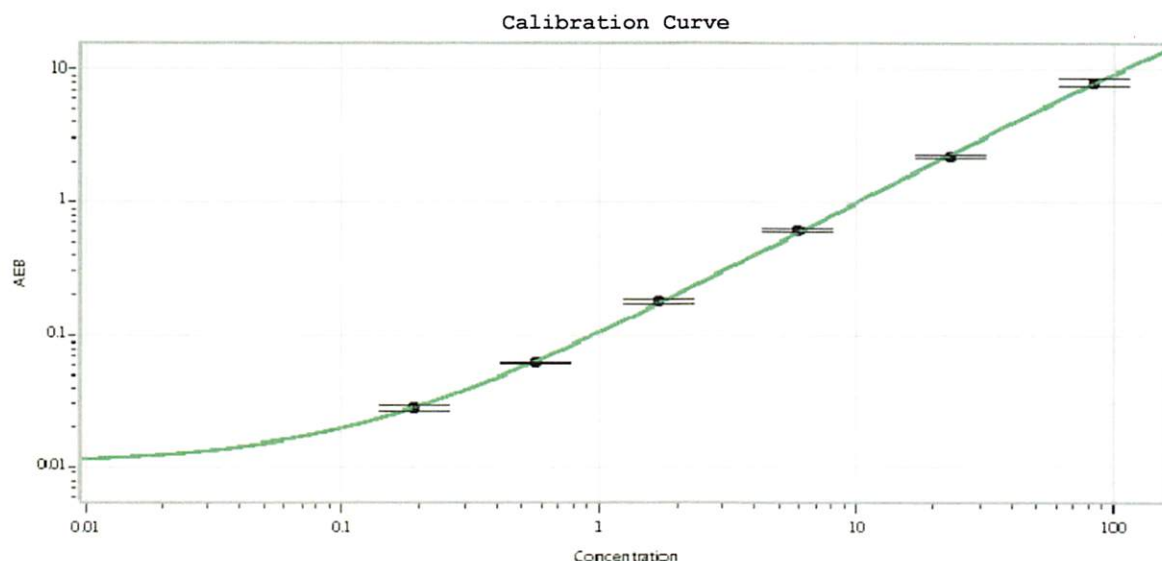
Component	Part Number	Lot Number
Bead Reagent	103698	106712
Detector Reagent	103700	106715
SBG Reagent	103702	106714
Sample Diluent	103704	107016
Reference Calibrator	103711	107004
Control 1	103712	107018
Control 2	103713	107018
RGP Reagent	103159	N/A ¹

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	107018	3.73	1.73-4.48	
Control 2	107018	133	89.4-158	
Panel 1	013903	0.821	0.258-1.11	
Panel 2	013903	2.06	0.841-2.44	
Panel 3	013903	3.47	1.79-4.31	



Calibrator Levels (pg/mL)	
A	0.000
B	0.191
C	0.560
D	1.70
E	5.93
F	23.1
G	83.8

Review/Approval

Junpil Park

QC Supervisor

Name

Title

Signature/Date

19 Mar 2021