

Product Number:	103714
Lot Number:	503410
Expiration:	3-Aug-2022
Platform(s):	HD-1 / HD-X & SR-X

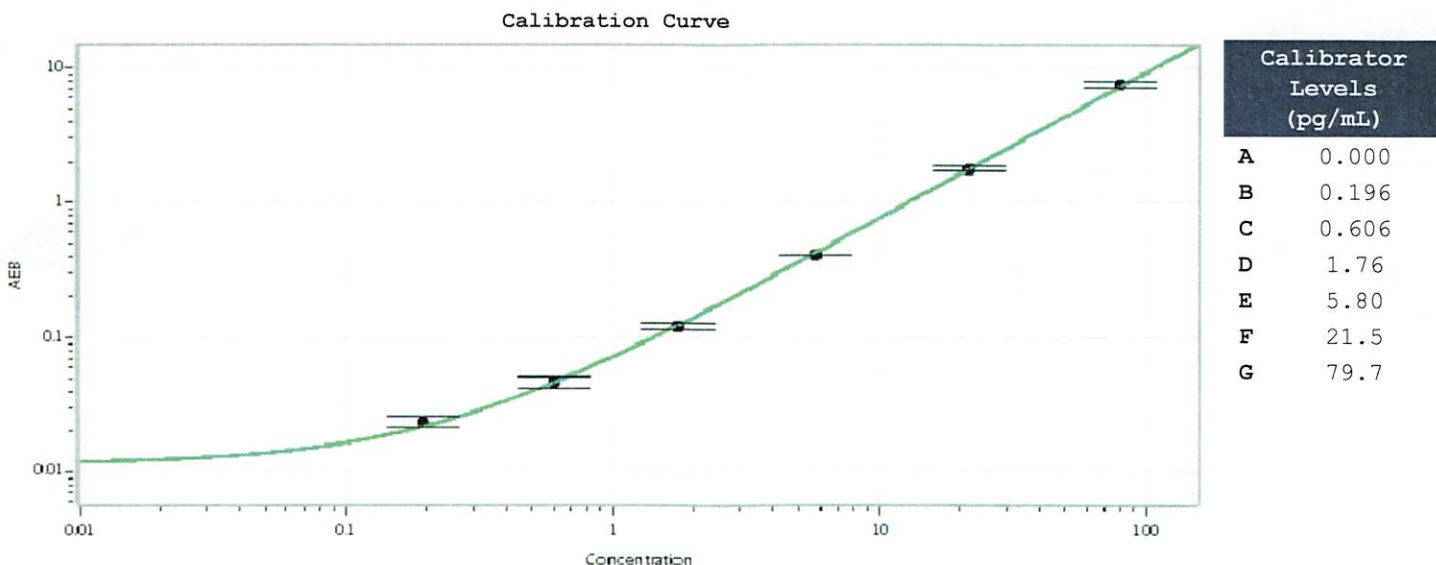


Data below represents results generated on the Simoa® HD-1 / HD-X & SR-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	103698	121508
Detector Reagent	103700	121509
SBG Reagent	103702	121605
Sample Diluent	103704	121511
Reference Calibrator	103711	123501
Control 1	103712	123502
Control 2	103713	123502
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

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	123502	4.14	2.88-4.77	
Control 2	123502	53.3	44.2-66.3	
Panel 1	013903	0.601	0.258-1.11	
Panel 2	013903	1.67	0.841-2.44	
Panel 3	013903	2.90	1.79-4.31	



Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 22 Mar 2022
Signature/Date