

Product Number:	103714
Lot Number:	502613
Expiration:	11-May-2021
Platform(s):	HD-1 / HD-X & SR-X

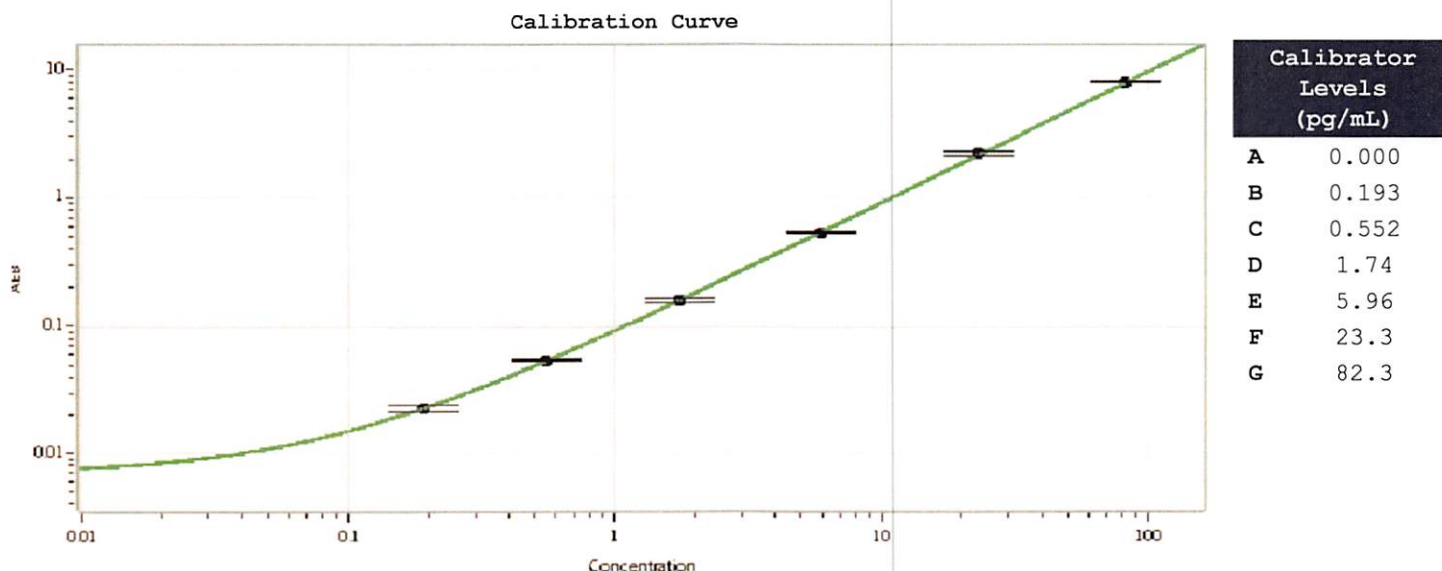


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	103698	031610
Detector Reagent	103700	031801
SBG Reagent	103702	031802
Sample Diluent	103704	031808
Reference Calibrator	103711	031601
Control 1	103712	031602
Control 2	103713	031602
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	031602	3.28	1.91-4.30	
Control 2	031602	34.5	20.8-62.0	
Panel 1	013903	0.355	0.230-1.08	
Panel 2	013903	1.09	0.855-2.26	
Panel 3	013903	2.50	1.74-4.18	



Review/Approval

Shivani Goel

QC Manager

Shivani Goel 03 Dec 2020

Name

Title

Signature/Date