

Product Number:	103714
Lot Number:	502508
Expiration:	6-Jul-2021
Platform(s):	HD-1 / HD-X & SR-X

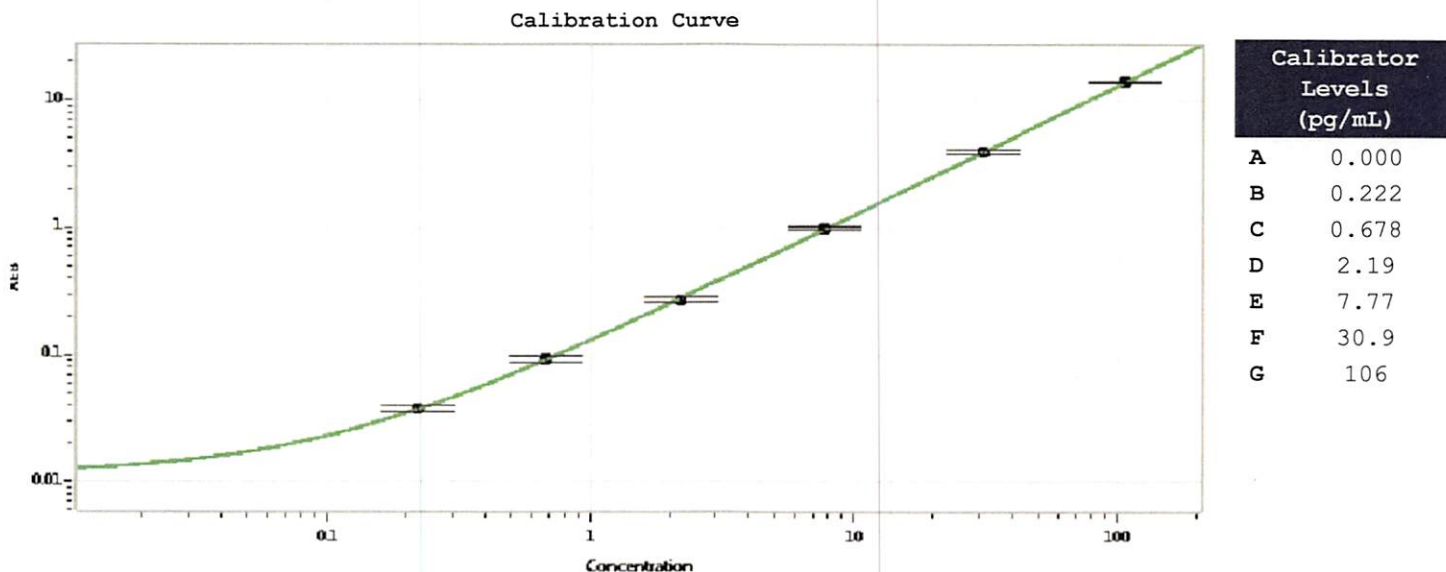


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	103698	018804
Detector Reagent	103700	018902
SBG Reagent	103702	018803
Sample Diluent	103704	019705
Reference Calibrator	103711	023201
Control 1	103712	021901
Control 2	103713	021901
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	021901	3.82	2.98-4.47	
Control 2	021901	62.2	48.8-75.1	
Panel 1	013903	0.769	0.575-0.916	
Panel 2	013903	1.68	1.35-2.04	
Panel 3	013903	2.68	1.73-4.89	



Review/Approval

Patrizia Stadler Assoc. Director, QC

Name

Title

P. Stadler 27 Aug 2020

Signature/Date