

Product Number: 100860
Lot Number: 502999
Expiration: 17-May-2022
Platform(s): HD-1/ HD-X

Component	Part Number	Lot Number
Bead Reagent	100865	115006
Detector Reagent	100862	115004
SBG Reagent	100868	115315
Sample Diluent	100870	114523
Calibrators	101111	116101
Control 1	100875	116102
Control 2	100876	116102
RGP Reagent	103159	N/A ¹

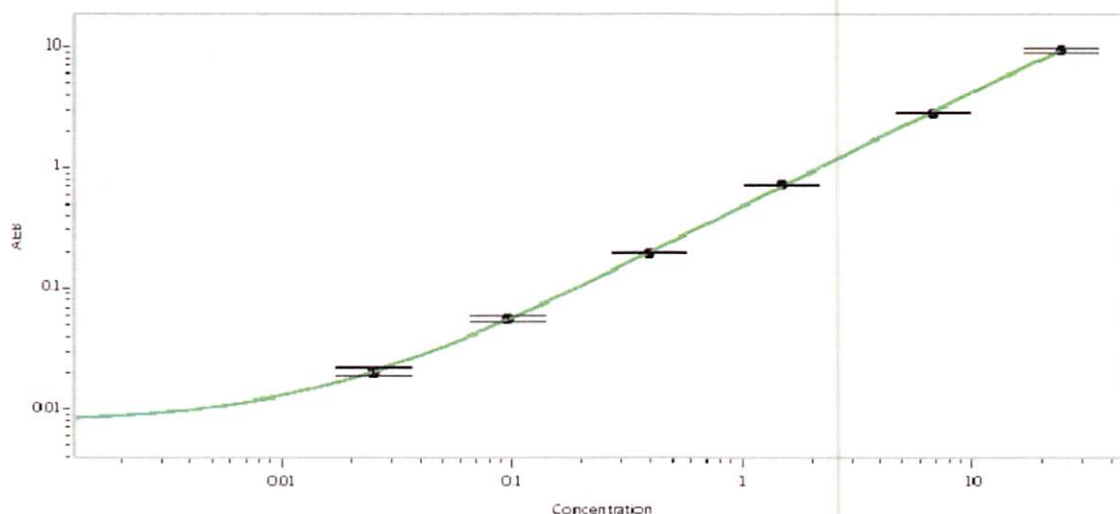
¹ RGP is not Kit Lot Specific



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	116102	4.58	3.48-5.50	
Control 2	116102	0.076	0.049-0.102	
Panel 1	115916	1.24	0.900-1.35	
Panel 2	115916	6.50	4.94-9.05	
Panel 3	115916	38.3	28.0-50.3	

Calibration Curve



Review/Approval: Shivani Goel, QC Manager, Signature/Date: Shrivam Goel 22 Jun 2021