

Product Number:	101599
Lot Number:	502253
Expiration:	26-Nov-2020
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	101601	005211
Detector Reagent	102012	005105
SBG Reagent	101602	005004
Sample Diluent	101603	004909
Calibrator Diluent	102014	003512
Calibrator Concentrate	100778	001414
RGP Reagent	103159	N/A ¹

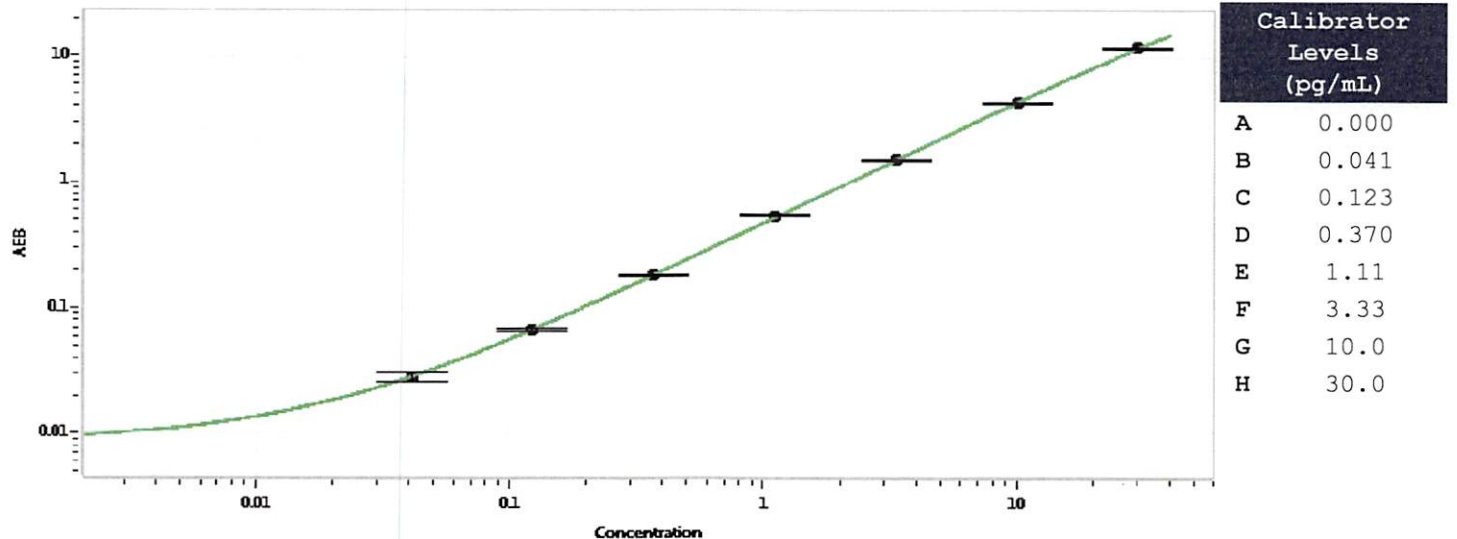
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-012518-02	0.443	0.352-0.537
Control 2	MF-012518-02	23.8	14.5-28.3
Panel 3	900903	8.92	6.49-9.73
Panel 4	900904	32.3	22.3-35.1

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	706
---	-----

Calibration Curve



Review/Approval

Patrizia Stadler Assoc. Director, QC

f. Stadler 28 March 20

Name

Title

Signature/Date