

Product Number:	101599
Lot Number:	502325
Expiration:	13-Mar-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	101601	010305
Detector Reagent	102012	007703
SBG Reagent	101602	007805
Sample Diluent	101603	007804
Calibrator Diluent	102014	007806
Calibrator Concentrate	100778	007810
RGP Reagent	103159	N/A ¹

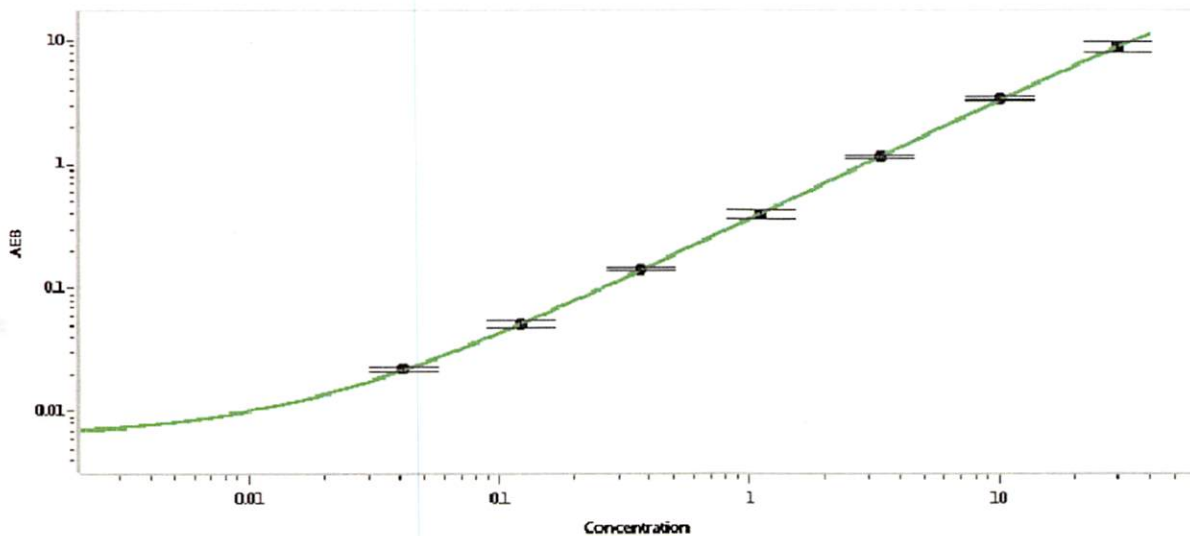
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-012518-02	0.410	0.352-0.537
Control 2	MF-012518-02	23.8	14.5-28.3
Panel 3	900903	8.68	6.55-9.82
Panel 4	900904	31.4	22.6-35.6

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

Calibrator Concentrate Value Assignment (pg/mL)	825
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval: Patrizia Stodler BC Associate Director 21 Apr 2020
Name Title Signature/Date