

Product Number:	101599
Lot Number:	502364
Expiration:	24-Mar-2021
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101601	012203
Detector Reagent	102012	012204
SBG Reagent	101602	012404
Sample Diluent	101603	012403
Calibrator Diluent	102014	010810
Calibrator Concentrate	100778	001414
RGP Reagent	103159	N/A ¹

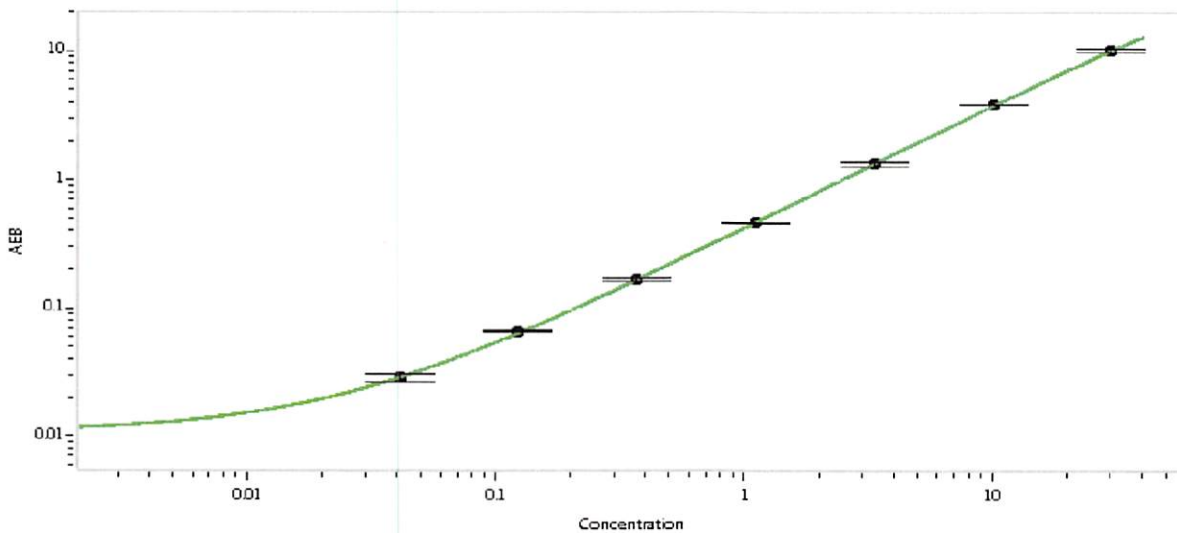
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-012518-02	0.506	0.352-0.537
Control 2	MF-012518-02	25.9	14.5-28.3
Panel 3	900903	10.2	6.37-10.7
Panel 4	900904	36.8	21.1-38.7

* 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



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

Shivani Goel
Name

QC Manager
Title

Shivani Goel 15 May 2020
Signature/Date