

Product Number:	100794
Lot Number:	502406
Expiration:	21-Feb-2021
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	100781	015703
Detector Reagent	100782	015105
SBG Reagent	100783	015012
Sample Diluent	103202	005235
Calibrator Diluent	100793	010814
Calibrator Concentrate	100780	932311
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹



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

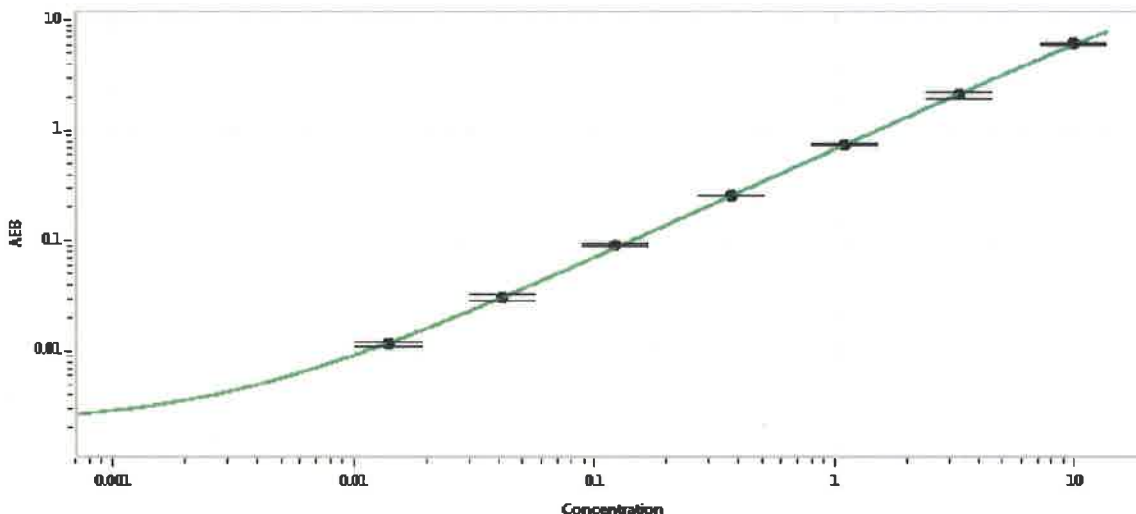
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-062316-02	9.63	7.17-10.8
Control 2	MF-062316-02	0.797	0.606-0.910
Panel 1	MF-110316-05	4.03	3.54-5.31
Panel 2	MF-110316-06	47.1	39.6-59.3
Panel 3	MF-110316-07	4.08	3.16-4.74

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

Calibrator Concentrate Value Assignment (pg/mL)	134
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.014
C	0.041
D	0.123
E	0.370
F	1.11
G	3.33
H	10.0

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 08 Sep 2020
Signature/Date

