

Product Number:	101605
Lot Number:	502314
Expiration:	13-Mar-2021
Platform(s):	HD-1/ HD-X



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

Component	Part Number	Lot Number
Bead Reagent	101607	009402
Detector Reagent	101606	009003
SBG Reagent	101608	009002
Sample Diluent	103200	009001
Calibrator Diluent	101610	009012
Calibrator Concentrate	100895	912609
RGP Reagent	103159	N/A ¹

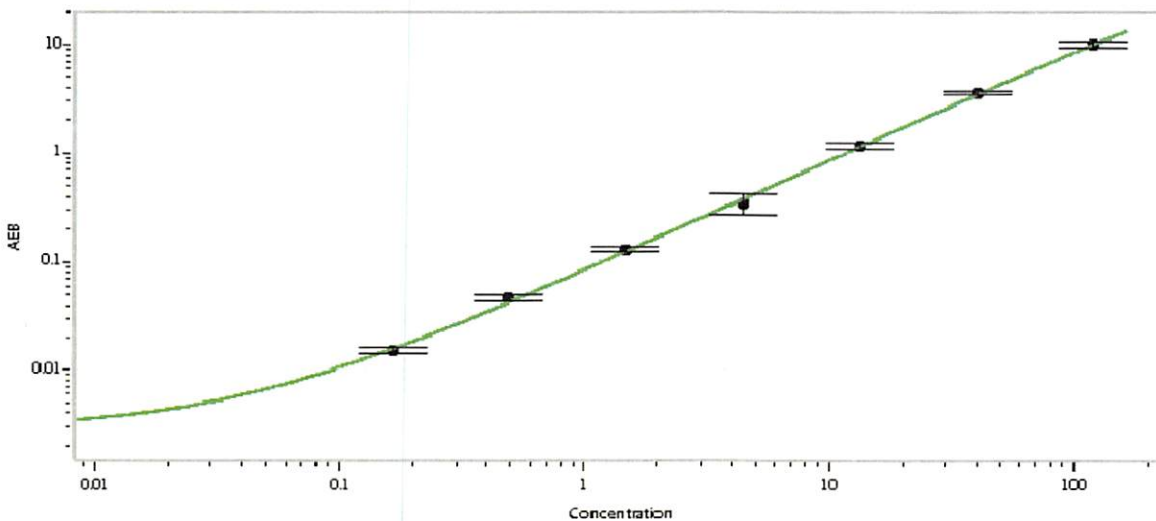
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	911406	3.29	1.00-3.71
Control 2	911406	186	87.0-203
Panel 1	829001	4.26	1.76-4.78
Panel 2	829002	8.00	3.80-8.88
Panel 3	829003	13.5	7.07-16.5
Panel 4	829004	59.2	38.5-65.6

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

Calibrator Concentrate Value Assignment (pg/mL)	6163
---	------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.165
C	0.494
D	1.48
E	4.44
F	13.3
G	40.0
H	120

Review/Approval

Patrizia Stadler

Name

Assoc. Director, QC

Title

P. Stadler

Signature/Date

08 Apr 2020