

Product Number:	102732
Lot Number:	501920
Expiration:	16-Jul-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	102722	919702
Detector Reagent	102724	920711
SBG Reagent	102726	920710
Sample Diluent	102728	920510
Calibrator Diluent	102731	920509
Calibrator Concentrate	100658	920712
RGP Reagent	103159	N/A ¹

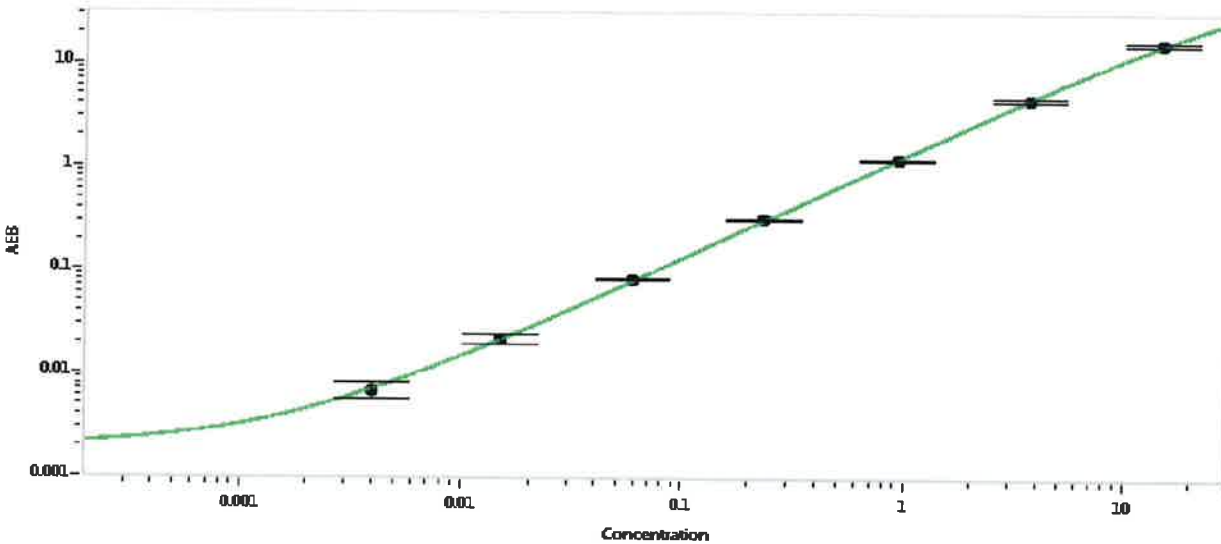
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	909502	0.135	0.097-0.158
Control 2	909502	3.67	3.03-4.54
Panel 1	911401	0.045	0.011-0.052
Panel 3	911403	1.59	1.30-2.00
Panel 4	911404	5.55	4.50-7.31

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

Calibrator Concentrate Value Assignment (pg/mL)	412
---	-----

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.004
C	0.015
D	0.059
E	0.234
F	0.938
G	3.75
H	15.0

Review/Approval: Patrizia Stadler Assoc. Director, QC *P. Stadler* 08 Aug 2019
Name Title Signature/Date