

Product Number:	102732
Lot Number:	501798
Expiration:	16-Apr-2020
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	102722	911205
Detector Reagent	102724	911605
SBG Reagent	102726	911606
Sample Diluent	102728	911607
Calibrator Diluent	102731	911312
Calibrator Concentrate	100658	908008
RGP Reagent	103159	N/A ¹

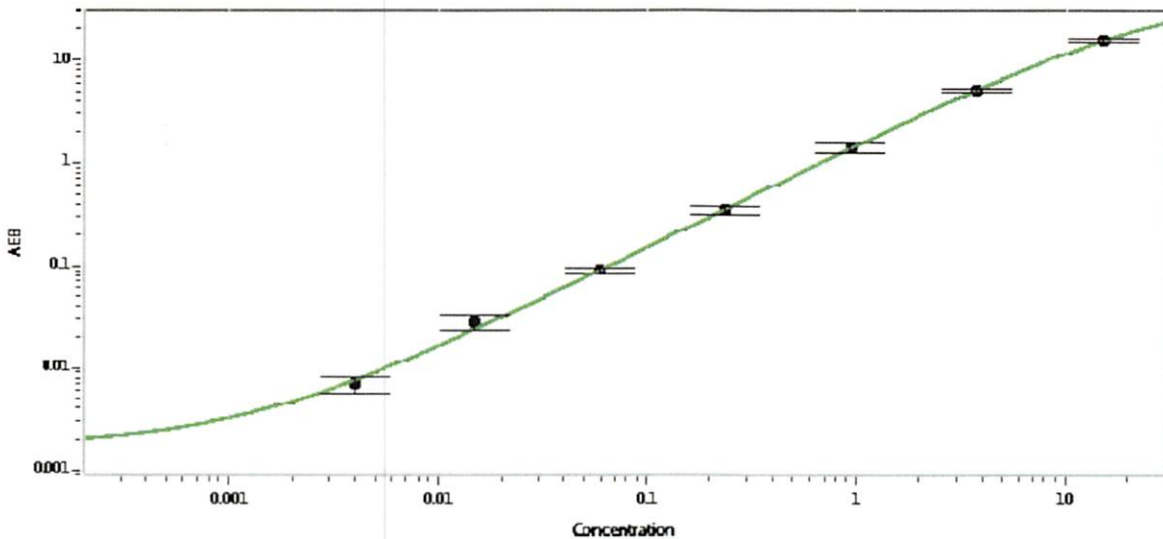
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	909502	0.125	0.097-0.156
Control 2	909502	3.21	3.04-4.56
Panel 1	911401	0.027	0.013-0.045
Panel 3	911403	1.36	1.31-2.02
Panel 4	911404	5.28	4.50-7.41

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

Calibrator Concentrate Value Assignment (pg/mL)	237
---	-----

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 Associate Director, QC
 Name: Patrizia Stadler Title: QC Signature/Date: P. Stadler 09 May 19