

Product Number:	102732
Lot Number:	502378
Expiration:	29-Jan-2021
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	012807
Detector Reagent	102724	012804
SBG Reagent	102726	012805
Sample Diluent	102728	012806
Calibrator Diluent	102731	010715
Calibrator Concentrate	100658	013605
RGP Reagent	103159	N/A ¹

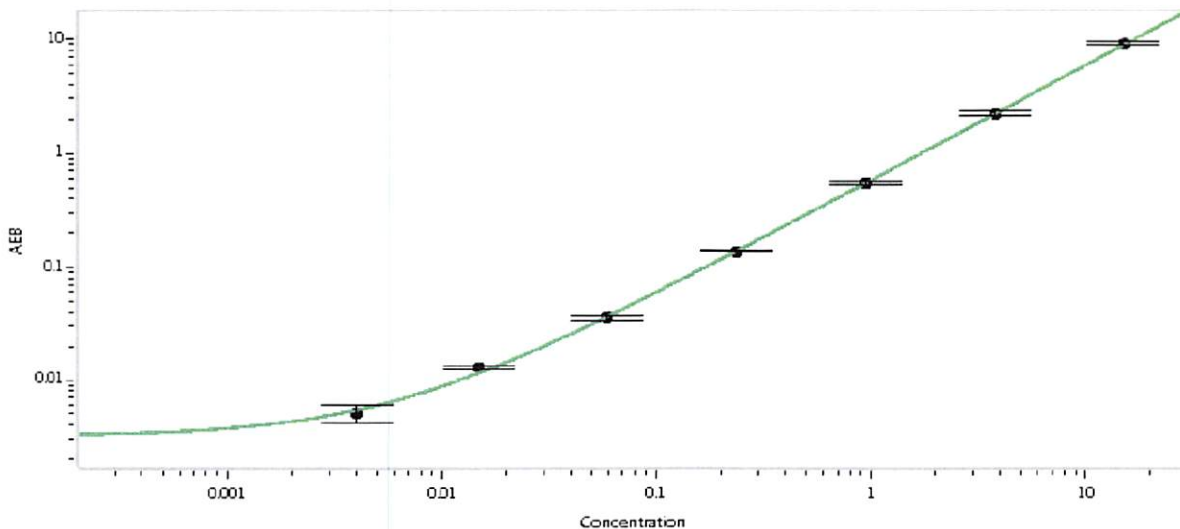
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	909502	0.124	0.095-0.162
Control 2	909502	3.25	2.88-4.52
Panel 1	911401	0.026	0.013-0.050
Panel 3	911403	1.53	1.22-2.02
Panel 4	911404	5.15	3.95-7.43

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

Calibrator Concentrate Value Assignment (pg/mL)	497
---	-----

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

Shivani Goel
Name

QC Manager
Title

Shivani Goel 27 May 2020
Signature/Date