

Product Number:	100988
Lot Number:	502775
Expiration:	2-Dec-2021
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	100974	105203
Detector Reagent	100976	105404
SBG Reagent	100987	105405
Sample Diluent	103734	103415
Calibrator Diluent	100980	033715
Calibrator Concentrate	100978	934402
RGP Reagent	103159	N/A ¹

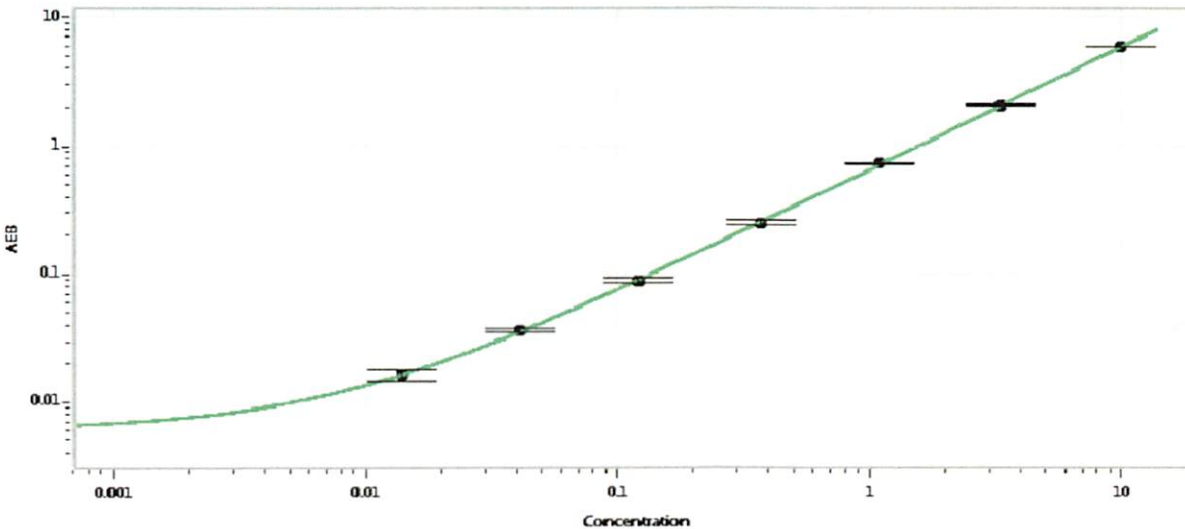
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108314	0.202	0.082-0.254
Control 2	108314	15.0	10.1-15.6
Panel 1	030001	0.215	0.134-0.227
Panel 3	030001	34.3	22.4-38.0
Panel 4	030001	0.184	0.129-0.210

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

Calibrator Concentrate Value Assignment (pg/mL)	225
---	-----

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	QC Manager	Shivani Goel (5Apr 2021)
	Name	Title	Signature/Date