

Product Number:	101622
Lot Number:	502564
Expiration:	3-Aug-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	101624	027909
Detector Reagent	101623	028002
SBG Reagent	101625	028108
Sample Diluent	103732	028106
Calibrator Diluent	101627	024716
Calibrator Concentrate	100652	021203
RGP Reagent	103159	N/A ¹

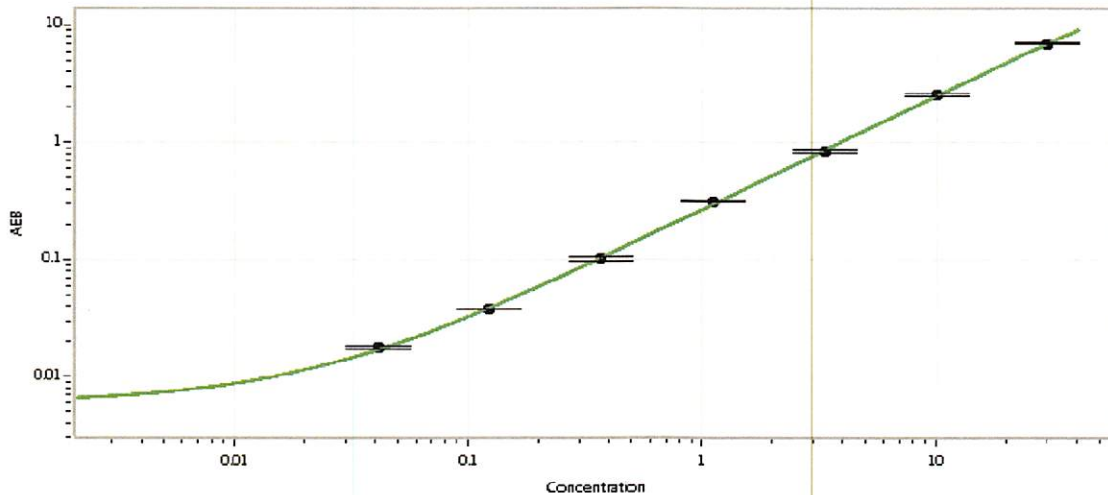
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	901702	1.10	0.944-1.42
Control 2	901702	16.9	15.4-23.1
Panel 1	029301	52.7	40.2-60.3
Panel 2	029401	5.25	4.33-6.50
Panel 3	029301	1.39	1.21-1.81

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

Calibrator Concentrate Value Assignment (pg/mL)	606
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Shivani Goel QC Manager
Patrizia Stadler Assoc. Director, QC
Name Title

Shivani Goel 04 Nov 2020
Signature/Date