

Product Number:	101622
Lot Number:	503273
Expiration:	15-May-2022
Platform(s):	HD-1 / HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101624	134905
Detector Reagent	101623	134915
SBG Reagent	101625	134916
Sample Diluent	103732	113510
Calibrator Diluent	101627	134920
Calibrator Concentrate	100652	126802
RGP Reagent	103159	N/A ¹

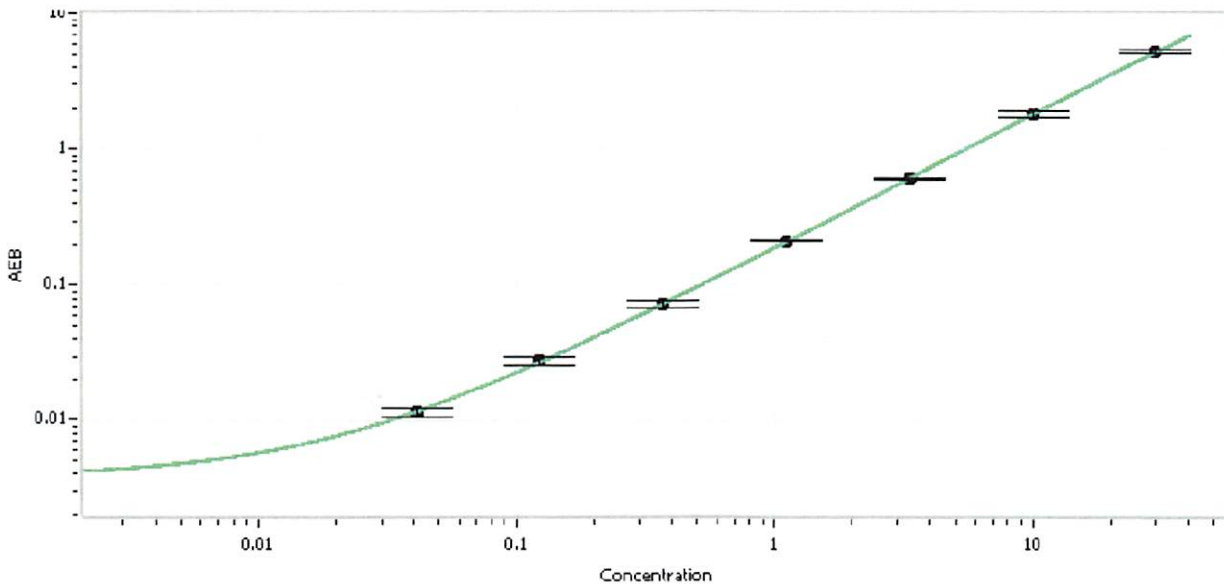
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	126802	2.08	1.50 - 2.50
Control 2	126802	19.5	15.0 - 25.0
Panel 1	029301	49.7	39.8 - 59.7
Panel 2	029401	5.91	4.35 - 6.53
Panel 3	029301	1.72	1.22 - 1.82

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

Calibrator Concentrate Value Assignment (pg/mL)	541
---	-----

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 Name: Shivani Goel Title: QC Signature/Date: Shivani Goel 04 Jan 2022