

Product Number:	101635
Lot Number:	503165
Expiration:	2-Oct-2022
Platform(s):	HD-1/ HD-X



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

Component	Part Number	Lot Number
Bead Reagent	101637	127504
Detector Reagent	101636	127510
SBG Reagent	101638	134402
Sample Diluent	101639	127602
Calibrator Diluent	101640	127608
Calibrator Concentrate	100661	127604
RGP Reagent	103159	N/A ¹

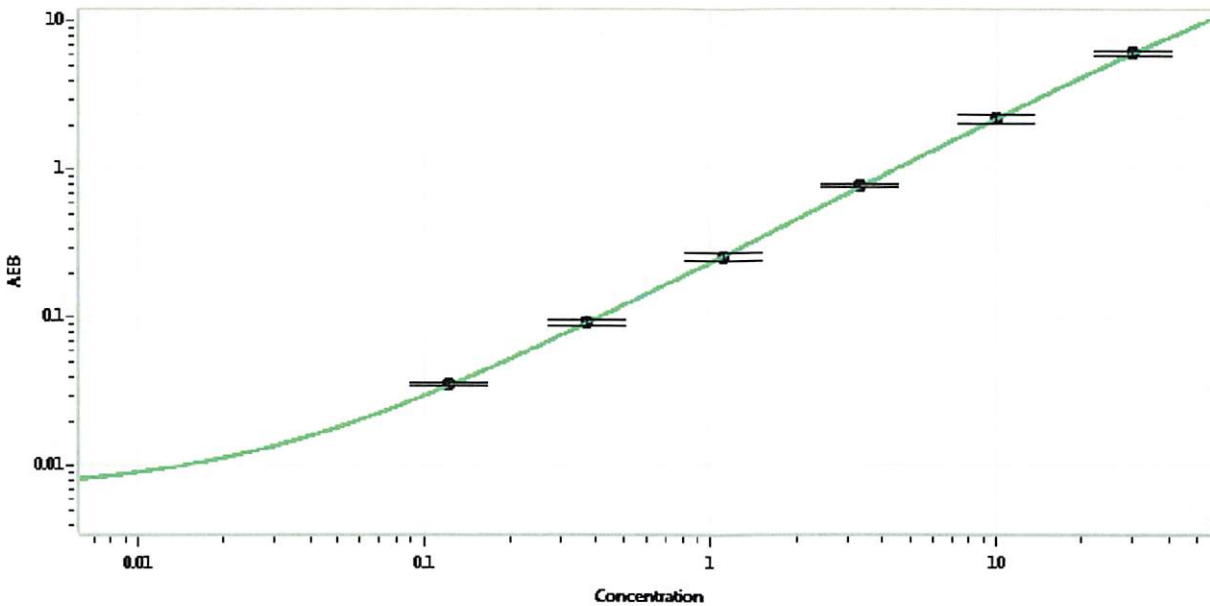
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	129010	0.543	0.384 - 0.631
Control 2	129010	58.3	37.8 - 62.4
Panel 1	035301	0.425	0.262 - 0.456
Panel 2	035301	4.62	3.25 - 4.88
Panel 3	035301	48.0	35.5 - 53.2

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

Calibrator Concentrate Value Assignment (pg/mL)	2119
---	------

Calibration Curve



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

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel
Signature/Date

05Jan2022
05Jan2022