

Product Number:	102329
Lot Number:	502956
Expiration:	21-Apr-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	102322	111702
Detector Reagent	102324	111818
SBG Reagent	102325	111717
Sample Diluent	102326	111117
Calibrator Diluent	102327	111722
Calibrator Concentrate	100849	108120
RGP Reagent	103159	N/A ¹

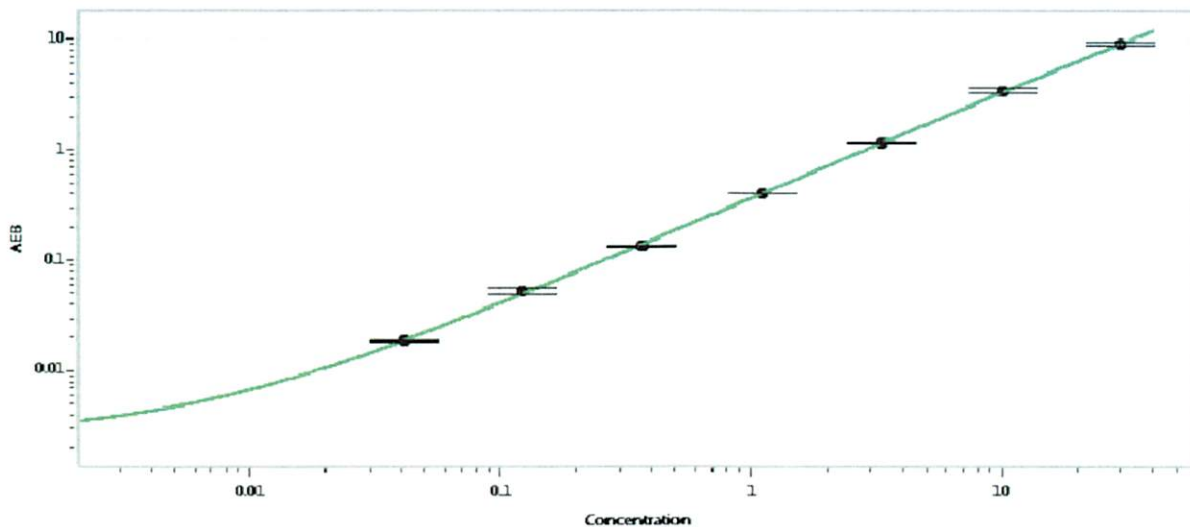
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	023702	0.211	0.114-0.249
Control 2	023702	27.6	17.6-33.7
Panel 1	027607	0.156	0.050-0.176
Panel 2	027607	0.171	0.071-0.213
Panel 3	027607	14.7	7.45-18.2

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

Calibrator Concentrate Value Assignment (pg/mL)	5585
---	------

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

Junpil Park
Name

QC Supervisor
Title

30 Jul 2021

Signature/Date