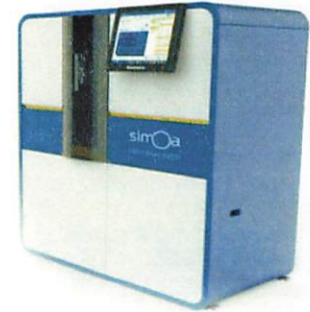


Product Number:	100196
Lot Number:	503378
Expiration:	5-Dec-2022
Platform(s):	HD-1/HD-X



Data below represents results generated on the Simoa® HD-1/ HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	100391	207206
Detector Reagent	100315	207305
SBG Reagent	100316	207306
Sample Diluent	100317	133939
Calibrator Diluent	100665	204515
Calibrator Concentrate	100664	121803
RGP Reagent	103159	N/A ¹

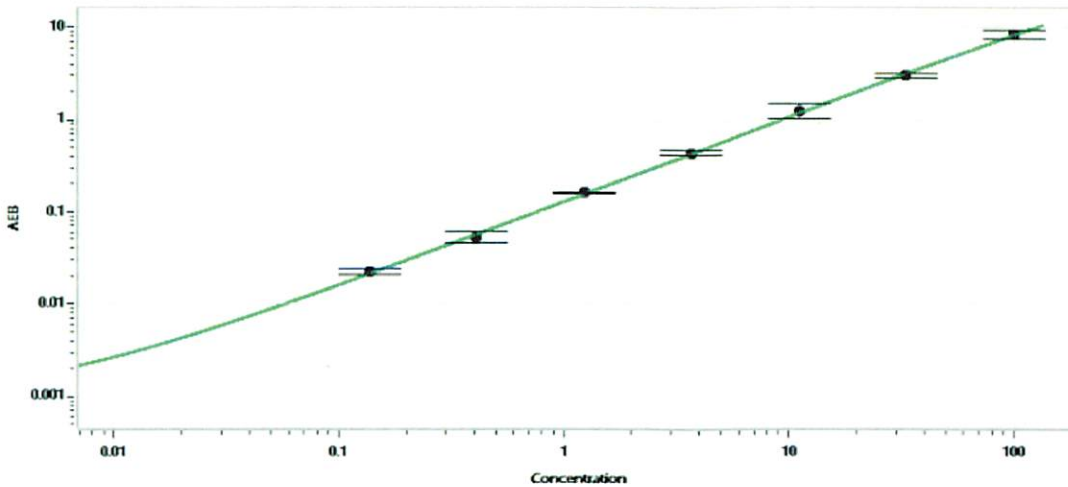
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	123410	10.2	6.13-11.5
Control 2	123410	127	48.0-142
Panel 1	029002	1.80	1.22-2.58
Panel 2	029002	10.0	5.51-12.5
Panel 3	029002	117	42.5-138

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

Calibrator Concentrate Value Assignment (pg/mL)	2110
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.24
E	3.70
F	11.1
G	33.3
H	100

Review/Approval	Shivani Goel	QC Manager	Shivani Goel	12 Apr 2022
	Name	Title	Signature/Date	