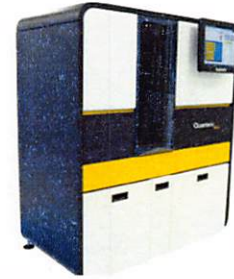


Product Number:	101580
Lot Number:	503364
Expiration:	15-Sep-2022
Platform(s):	HD-1/HD-x & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101582	125807
Detector Reagent	101581	126001
SBG Reagent	101583	126002
Sample Diluent	103735	126008
Calibrator Diluent	101585	126004
Calibrator Concentrate	100655	126005
RGP Reagent	103159	N/A ¹

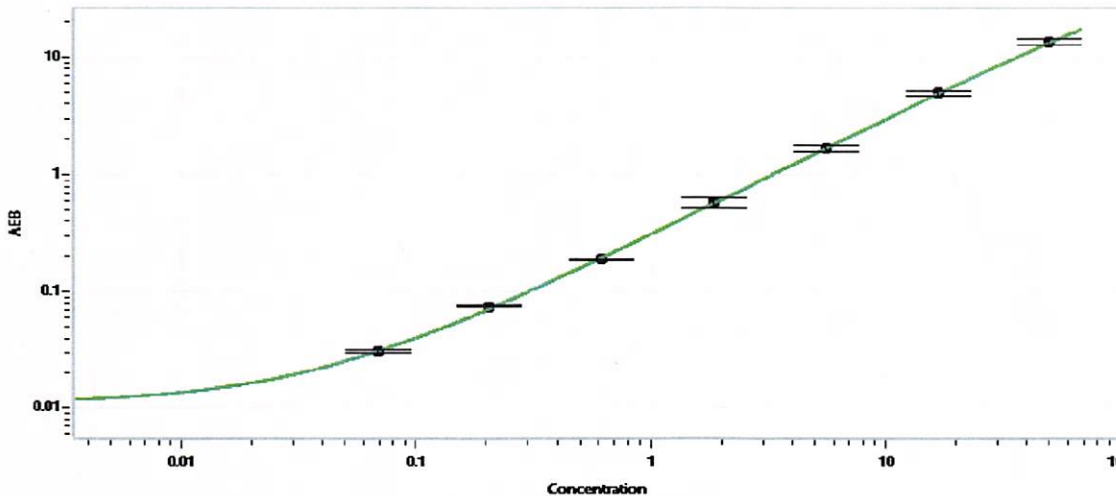
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	206502	3.16	2.16-3.40
Control 2	206502	52.0	35.3-57.4
Panel 2	115317	2.70	1.51-3.72
Panel 3	116301	30.1	15.5-39.3

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

Calibrator Concentrate Value Assignment (pg/mL)	1120
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.069
C	0.206
D	0.617
E	1.85
F	5.56
G	16.7
H	50.0

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 05 May 2022
Signature/Date