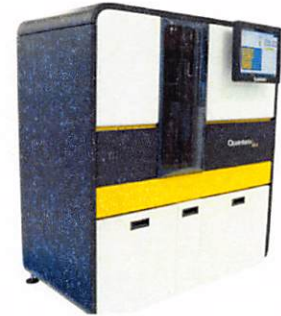


Product Number:	101664
Lot Number:	503408
Expiration:	27-Feb-2023
Platform(s):	HD-1/HD-X & SR-X



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

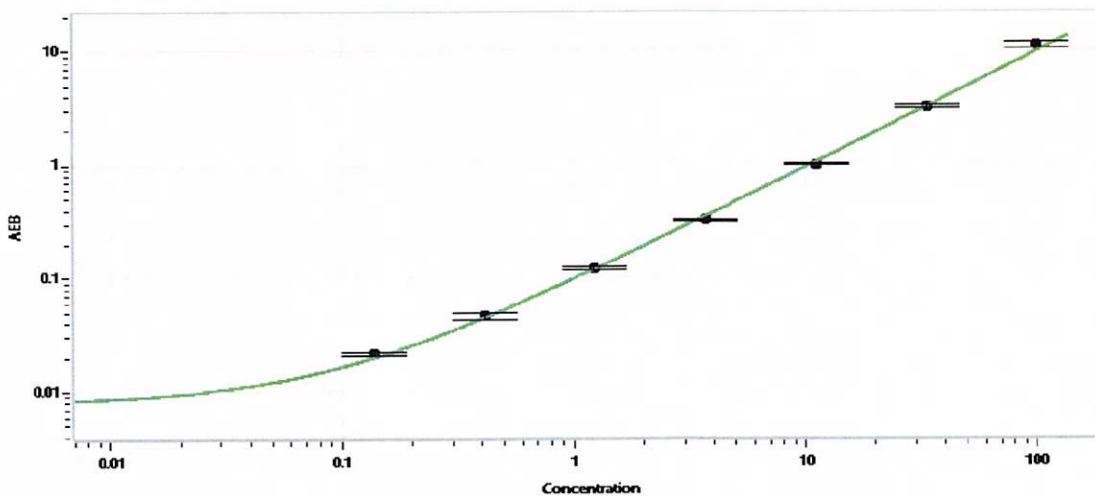
Component	Part Number	Lot Number
Bead Reagent	101666	207701
Detector Reagent	101665	207801
SBG Reagent	101667	207802
Sample Diluent	101668	207803
Calibrator Diluent	101669	207804
Calibrator Concentrate	103217	202410
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	201802	1.26	0.681-1.45	
Control 2	201802	70.8	45.8-78.9	
Panel 1	202604	8.82	5.87-9.46	
Panel 2	202604	50.6	31.3-58.5	
Panel 3	202604	148	81.7-158	

Calibrator Concentrate Value Assignment (pg/mL)	1417
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval: Shivani Goel QC Manager Shivani Goel 06 Apr 2022

Name Title Signature/Date