

Product Number:	101478
Lot Number:	503261
Expiration:	5-Dec-2022
Platform(s):	HD-1/HD-X & SR-X



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

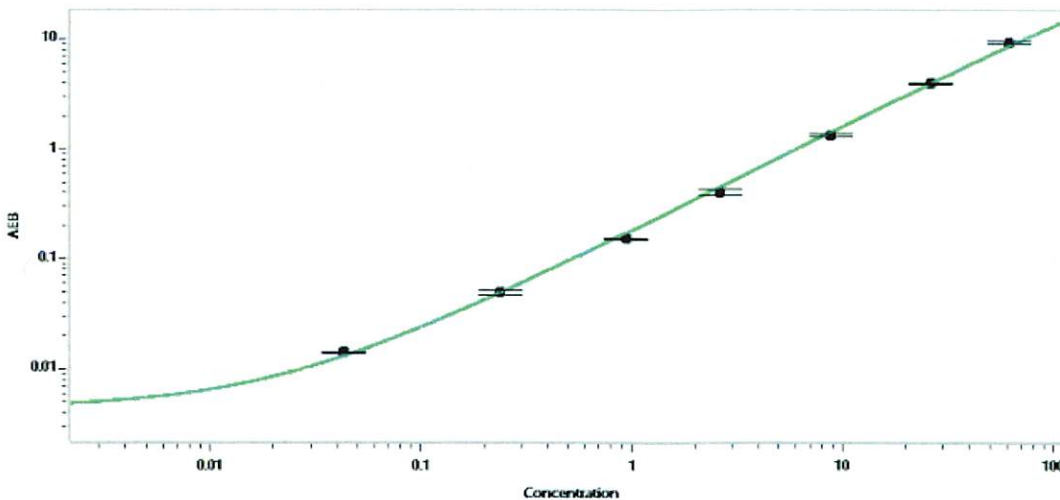
Component	Part Number	Lot Number
Bead Reagent	101480	133907
Detector Reagent	101481	134011
SBG Reagent	101482	134012
Sample Diluent	101483	134013
Calibrators	100758-100765	121603
Control 1	100768	121604
Control 2	100769	121604
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	121604	2.41	1.70-3.55
Control 2	121604	55.4	43.1-64.6
Panel 1	122912	0.228	0.155-0.315
Panel 2	122912	4.06	3.49-5.97
Panel 3	122912	51.2	40.2-70.5

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

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.043
C	0.237
D	0.937
E	2.62
F	8.76
G	26.2
H	61.6

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 12 Apr 2022  
Signature/Date