

Product Number:	103464
Lot Number:	503066
Expiration:	19-Jul-2022
Platform(s):	HD-1 / HD-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

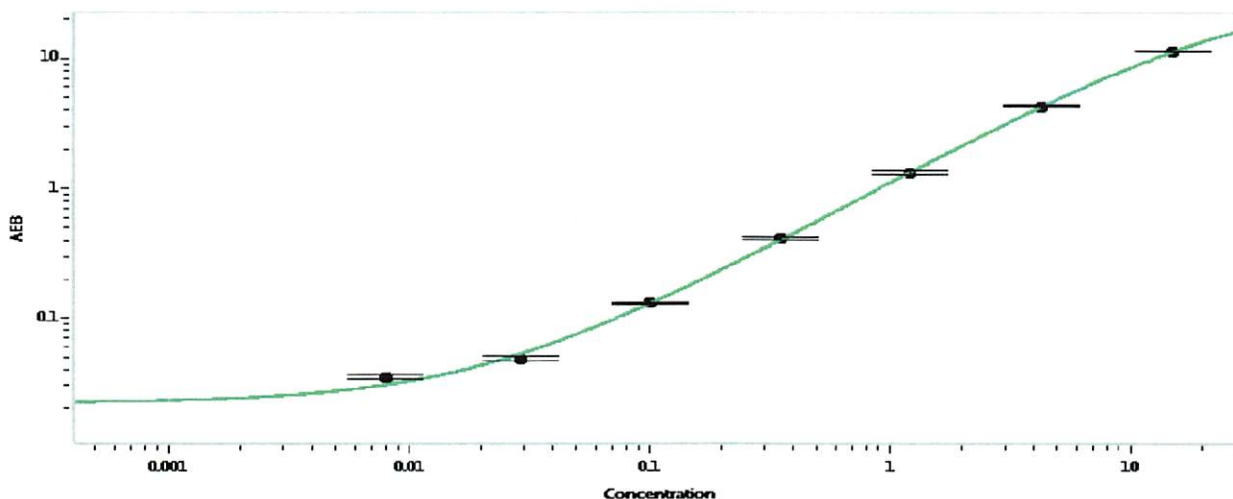
Component	Part Number	Lot Number
Bead Reagent	103452	120203
Detector Reagent	103454	120402
SBG Reagent	103456	120403
Sample Diluent	103458	120027
Calibrator Diluent	103460	120028
Calibrator Concentrate	103459	120904
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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	027502	0.220	0.130 - 0.313	
Control 2	027502	11.8	6.59 - 13.2	
Panel 1	106413	0.107	0.085 - 0.128	
Panel 2	106413	0.039	0.032 - 0.072	
Panel 3	106413	5.59	3.00 - 7.86	
Panel 4	106413	21.1	10.3 - 29.9	

Calibrator Concentrate Value Assignment (pg/mL)	259
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.008
C	0.029
D	0.100
E	0.350
F	1.22
G	4.29
H	15.0

Review/Approval

*Shivani Goel*  
Name

QC  
Title

*Shivani Goel*  
Signature/Date

*23 Aug 2021*  
*SG*  
23 Aug 2021