

Product Number:	103186
Lot Number:	502496
Expiration:	27-May-2021
Platform(s):	HD-1/ HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	102246	022302
Detector Reagent	102248	022404
SBG Reagent	102250	022403
Sample Diluent	102252	017008
Calibrators	102255	022401
Control 1	102256	022402
Control 2	102257	022402
RGP Reagent	103159	N/A <sup>1</sup>

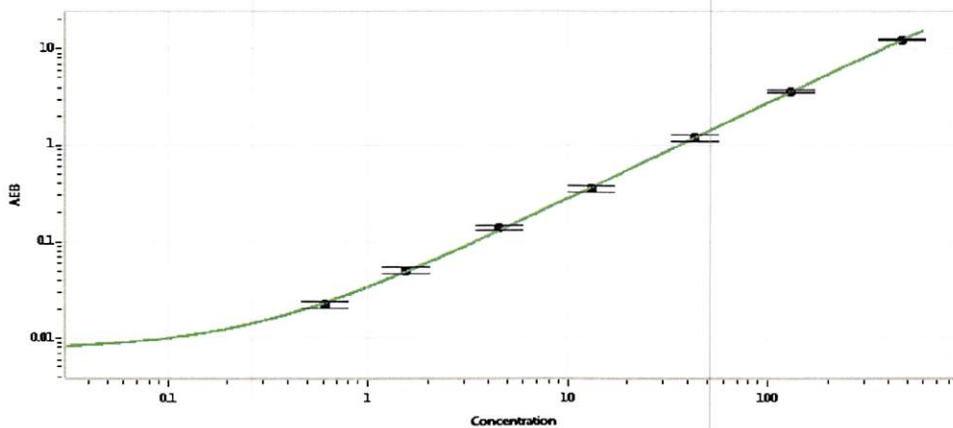
<sup>1</sup> RGP is not Kit Lot Specific



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

	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	022402	4.04	3.33-5.18	
Control 2	022402	188	134-212	
Panel 1	926601	18.7	14.1-21.1	
Panel 2	922602	6.56	4.94-7.41	
Panel 3	922603	42.3	28.4-42.6	
Panel 4	922604	160	129-204	

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.609
C	1.56
D	4.56
E	13.1
F	43.1
G	131
H	468

Review/Approval: Shivani Goel QC Manager  
Name Title

Shivani Goel 24 Aug 2020  
Signature/Date