

Product Number:	100860
Lot Number:	502576
Expiration:	9-Jun-2021
Platform(s):	HD-1/ HD-X



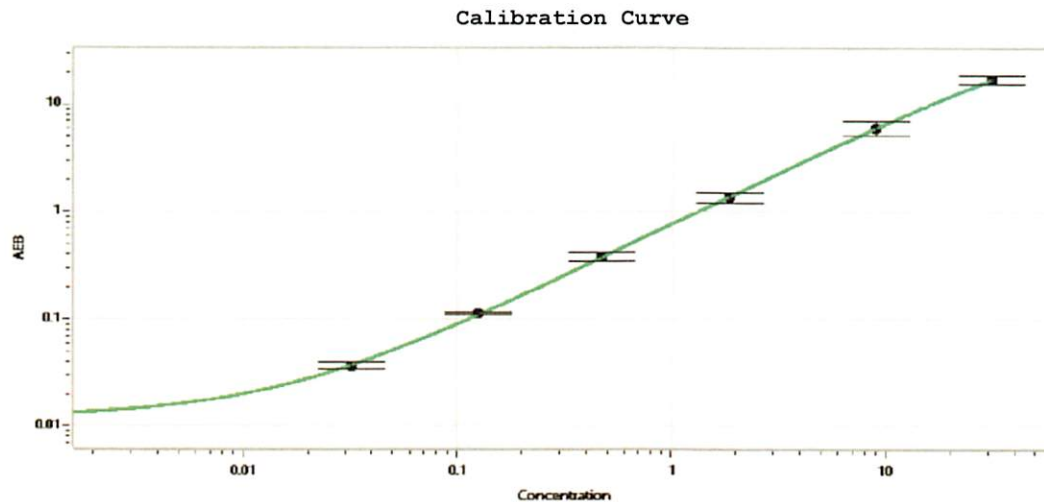
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	100865	028809
Detector Reagent	100862	029307
SBG Reagent	100868	016108
Sample Diluent	100870	028703
Calibrators	101111	016901
Control 1	100875	016902
Control 2	100876	016902
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	016902	6.73	4.99-7.48
Control 2	016902	0.112	0.076-0.118
Panel 1	908501	1.01	0.631-2.10
Panel 2	908502	6.27	4.38-12.5
Panel 3	908503	21.6	12.1-47.2

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



Calibrator Levels (pg/mL)	
A	0.000
B	0.032
C	0.126
D	0.470
E	1.85
F	8.88
G	30.9

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	25 Nov 2020
	Name	Title	Signature/Date	