

Product Number:	100860
Lot Number:	502385
Expiration:	29-Apr-2021
Platform(s):	HD-1/ HD-X



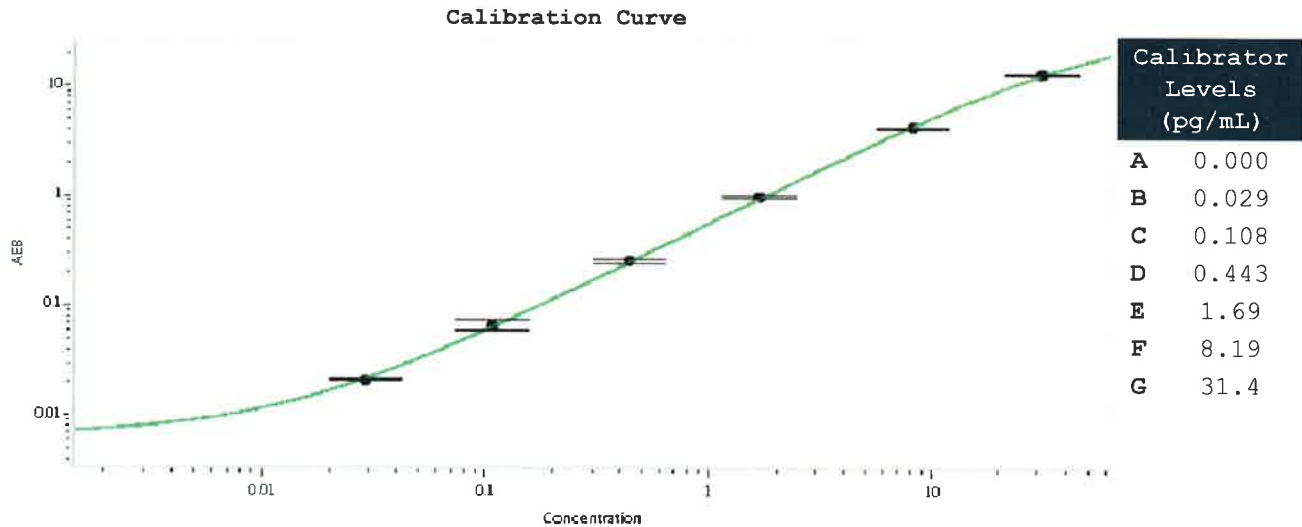
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	100865	013702
Detector Reagent	100862	013206
SBG Reagent	100868	013801
Sample Diluent	100870	012911
Calibrators	101111	013501
Controls	100875 & 100876	013502
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	013502	5.92	4.01-7.68
Control 2	013502	0.099	0.072-0.108
Panel 1	908501	1.30	0.631-2.10
Panel 2	908502	8.60	4.38-12.5
Panel 3	908503	26.6	12.1-47.2

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



Review/Approval

Shivani Goel

QC Manager

*Shivani Goel 22 May 2020*

Name

Title

Signature/Date