

Product Number:	100860
Lot Number:	502384
Expiration:	30-Jan-2021
Platform(s):	HD-1/ HD-X



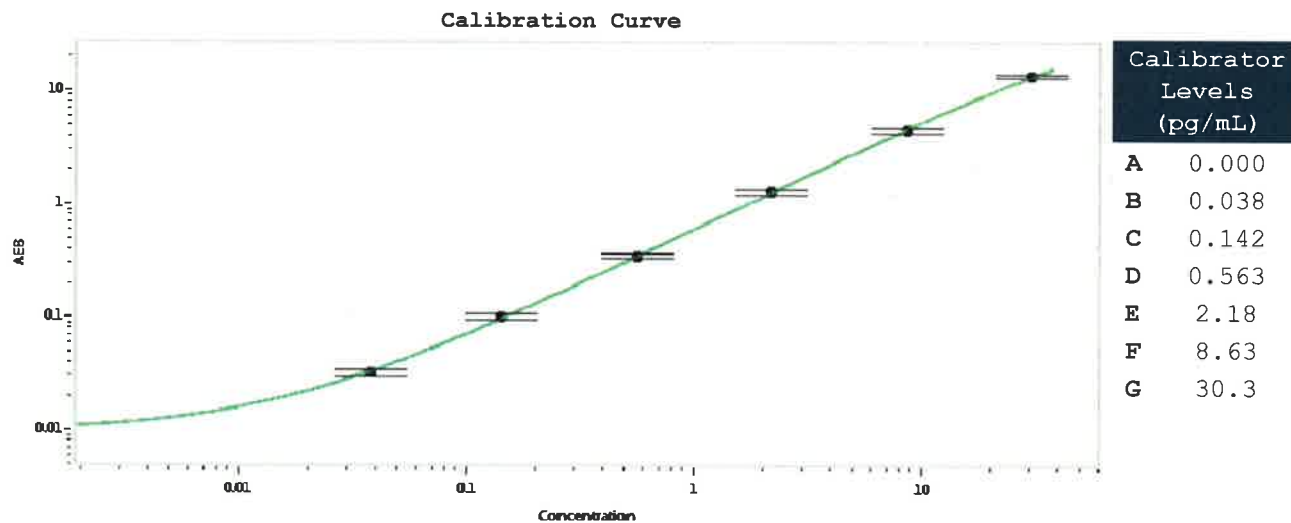
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	100865	013401
Detector Reagent	100862	007206
SBG Reagent	100868	013801
Sample Diluent	100870	003004
Calibrators	101111	923904
Controls	100875 & 100876	923905
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	923905	5.91	2.97-8.92
Control 2	923905	0.102	0.053-0.158
Panel 1	908501	1.14	0.631-2.10
Panel 2	908502	7.64	4.38-12.5
Panel 3	908503	25.3	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