

Product Number:	100860
Lot Number:	502025
Expiration:	3-Sep-2020
Platform(s):	HD-1/ HD-X



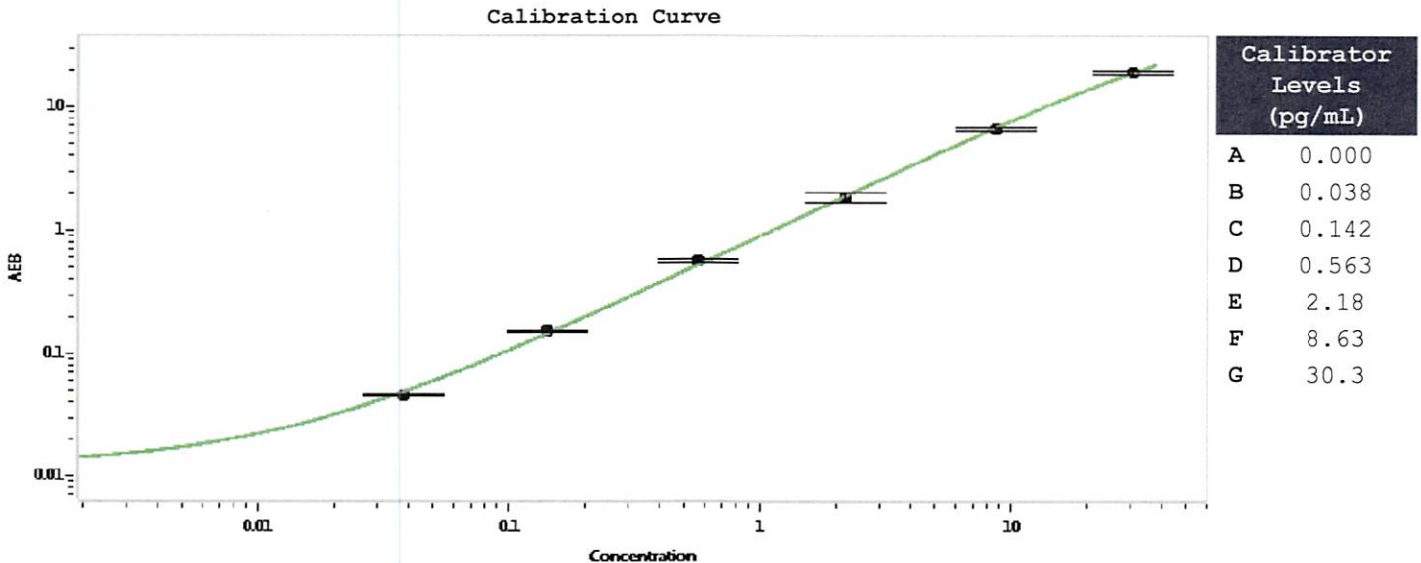
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	100865	924608
Detector Reagent	100862	924605
SBG Reagent	100868	924607
Sample Diluent	100870	924610
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	6.04	2.97-8.92
Control 2	923905	0.110	0.053-0.158
Panel 1	908501	1.18	1.04-1.95
Panel 2	908502	8.12	6.92-12.1
Panel 3	908503	26.1	21.5-45.4

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



Review/Approval	Patrizia Stadler Name	Assoc. Director, QC Title	<i>P. Stadler</i> Signature/Date	17 Sep 2019 Date
-----------------	--------------------------	------------------------------	-------------------------------------	---------------------