

Product Number:	103186
Lot Number:	502342
Expiration:	24-Mar-2021
Platform(s):	HD-1/ HD-X



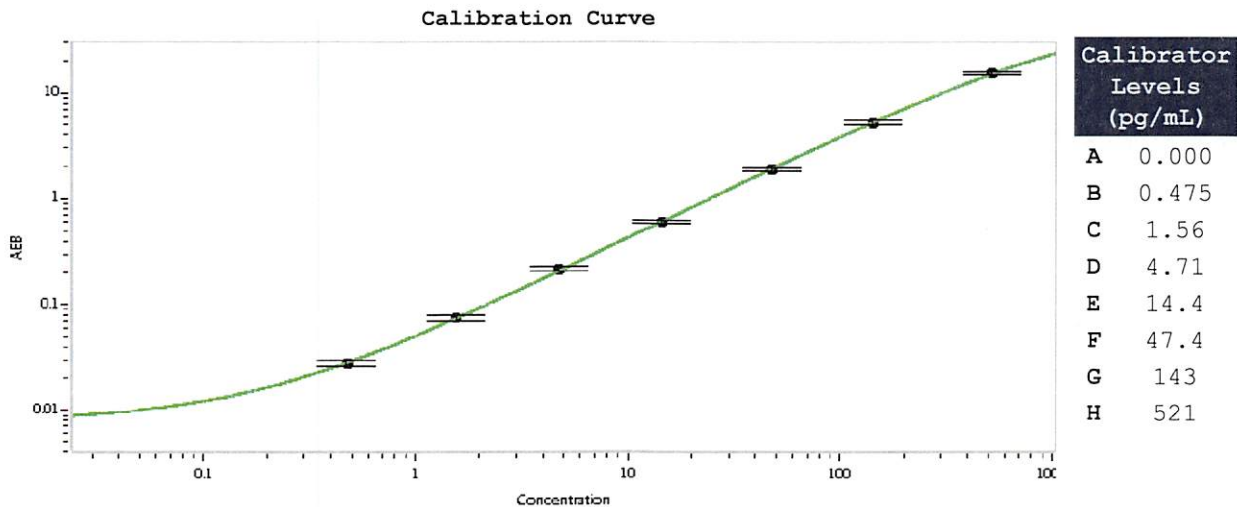
Data below represents results generated on the Simoa™ HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102246	010802
Detector Reagent	102248	010803
SBG Reagent	102250	010902
Sample Diluent	102252	010903
Calibrators	102255	011201
Controls	102256 - 102257	011301
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	011301	4.65	3.73-5.68
Control 2	011301	182	149 - 223
Panel 1	922601	18.0	14.1 - 21.1
Panel 2	922602	6.55	4.94 - 7.41
Panel 3	922603	40.9	28.4 - 42.6
Panel 4	922604	171	129 - 204

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



Review/Approval: Shivani Goel @CManager Shivani Goel 06 May 2020  
 Name Title Signature/Date