

Product Number:	103180
Lot Number:	502024
Expiration:	9-Sep-2020
Platform(s):	HD-1/HD-x & SR-X

Component	Part Number	Lot Number
Bead Reagent	102246	925415
Detector Reagent	102248	925205
SBG Reagent	102250	925204
Sample Diluent	102252	925202
Calibrators	102255	924701
Controls	102256 - 102257	924803
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

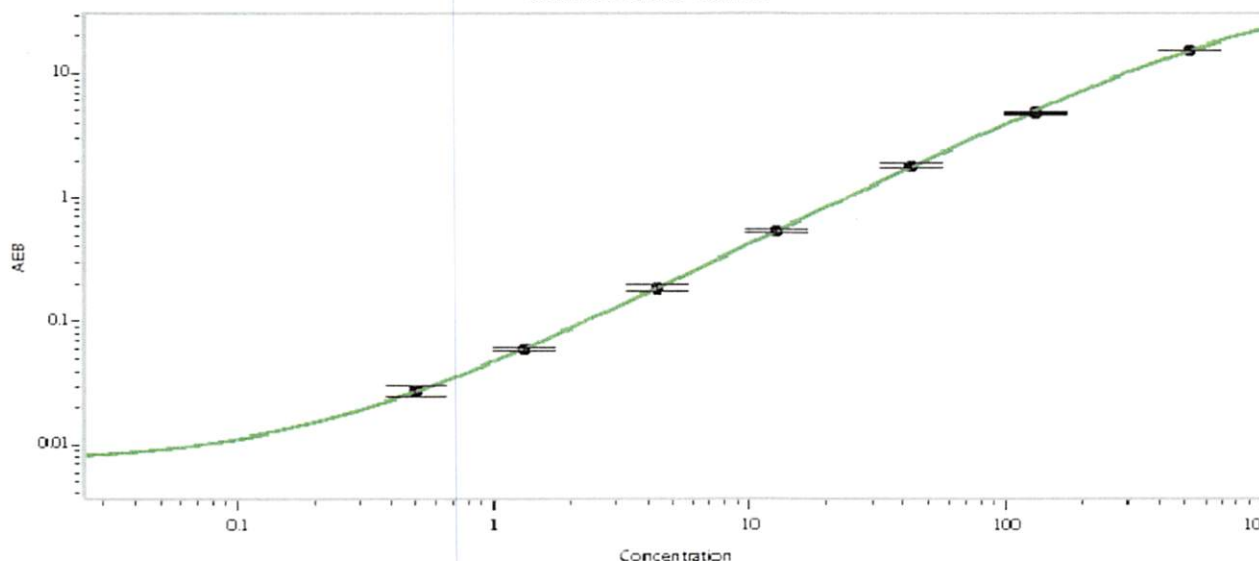


Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	924803	3.27	2.50-4.49
Control 2	924803	144	116-173
Panel 1a	823301	4.70	4.08-6.68
Panel 1b	823302	5.93	4.90-7.61
Panel 2	823601	80.1	59.3-101
Panel 3	823602	303	236-402

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.503
C	1.33
D	4.32
E	12.8
F	42.3
G	130
H	521

Review/Approval Patrizia Stadler Assoc. Director, QC P. Stoller 16 Sep 2019
 Name Title Signature/Date