

Product Number:	103186
Lot Number:	502057
Expiration:	23-Sep-2020
Platform(s):	HD-1/HD-x

Component	Part Number	Lot Number
Bead Reagent	102246	927001
Detector Reagent	102248	926801
SBG Reagent	102250	926608
Sample Diluent	102252	926609
Calibrators	102255	927405
Controls	102256 - 102257	927502
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

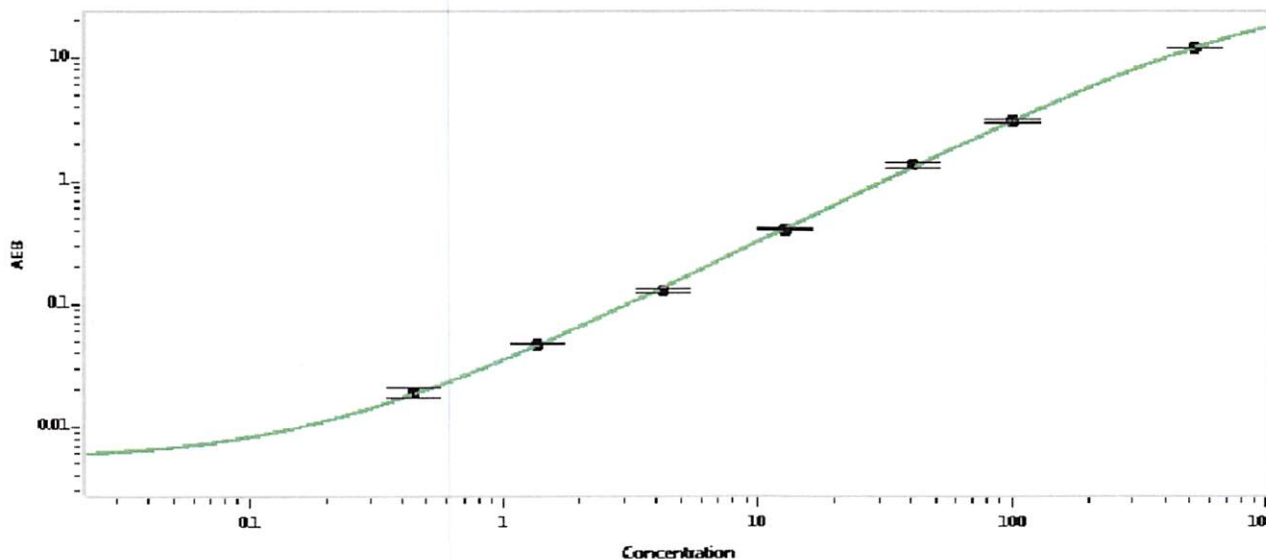


Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	927502	4.13	3.20-4.80
Control 2	927502	147	130-195
Panel 1a	922601	16.7	14.1-21.1
Panel 1b	922602	5.64	4.94-7.41
Panel 2	922603	34.7	28.4-42.6
Panel 3	922604	147	129-204

* 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	0.445
C	1.36
D	4.22
E	12.8
F	40.7
G	100
H	527

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

15 Oct 2019

Name

Title

Signature/Date