

Product Number:	103186
Lot Number:	502036
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	925601
Controls	102256 - 102257	925901
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

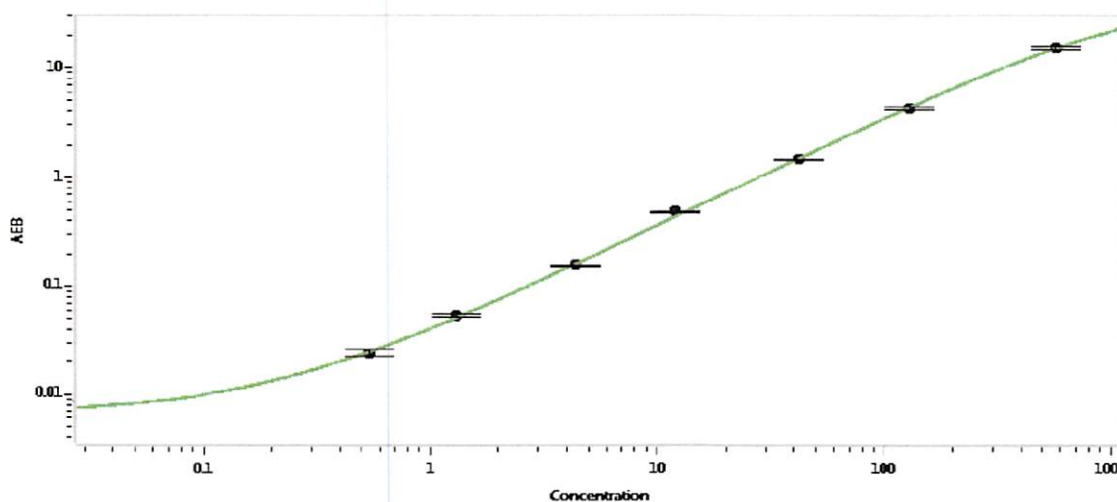


Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	925901	3.89	3.07-4.61
Control 2	925901	152	111-170
Panel 1a	922601	18.8	14.0-21.0
Panel 1b	922602	6.55	4.95-7.42
Panel 2	922603	38.0	28.1-42.1
Panel 3	922604	175	134-202

* 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	0.542
C	1.31
D	4.34
E	12.0
F	41.8
G	129
H	567

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	26 Sep 2019
	Name	Title	Signature/Date	