

Product Number:	103186
Lot Number:	502255
Expiration:	29-Jan-2021
Platform(s):	HD-1/HD-x, SRx

Component	Part Number	Lot Number
Bead Reagent	102246	005504
Detector Reagent	102248	005005
SBG Reagent	102250	005001
Sample Diluent	102252	005002
Calibrators	102255	003501
Controls	102256 - 102257	003601
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

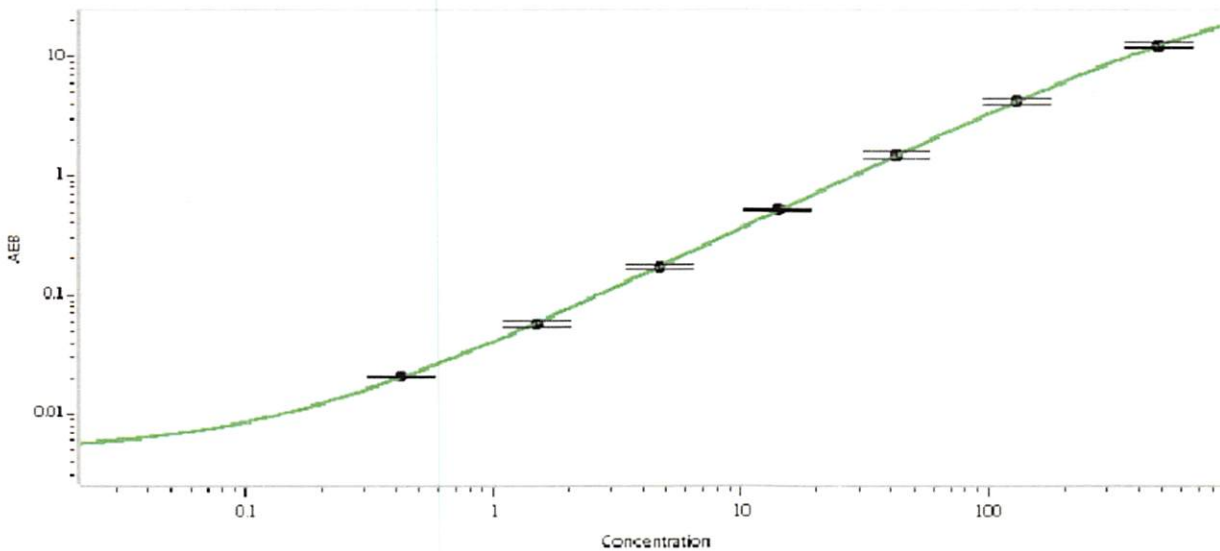


Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	003601	4.29	3.33-4.99
Control 2	003601	166	114-234
Panel 1	922601	15.9	14.1-21.1
Panel 2	922602	5.81	4.94-7.41
Panel 3	922603	33.8	28.4-42.6
Panel 4	922604	156	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	0.419
C	1.50
D	4.66
E	14.0
F	42.5
G	130
H	484

Review/Approval Patrizia Stadler Assoc. Director, QC *P. Stadler* 27 Feb 2020
Name Title Signature/Date