

Product Number:	103186
Lot Number:	502234
Expiration:	23-Sep-2020
Platform(s):	HD-1/HD-x, SRx



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

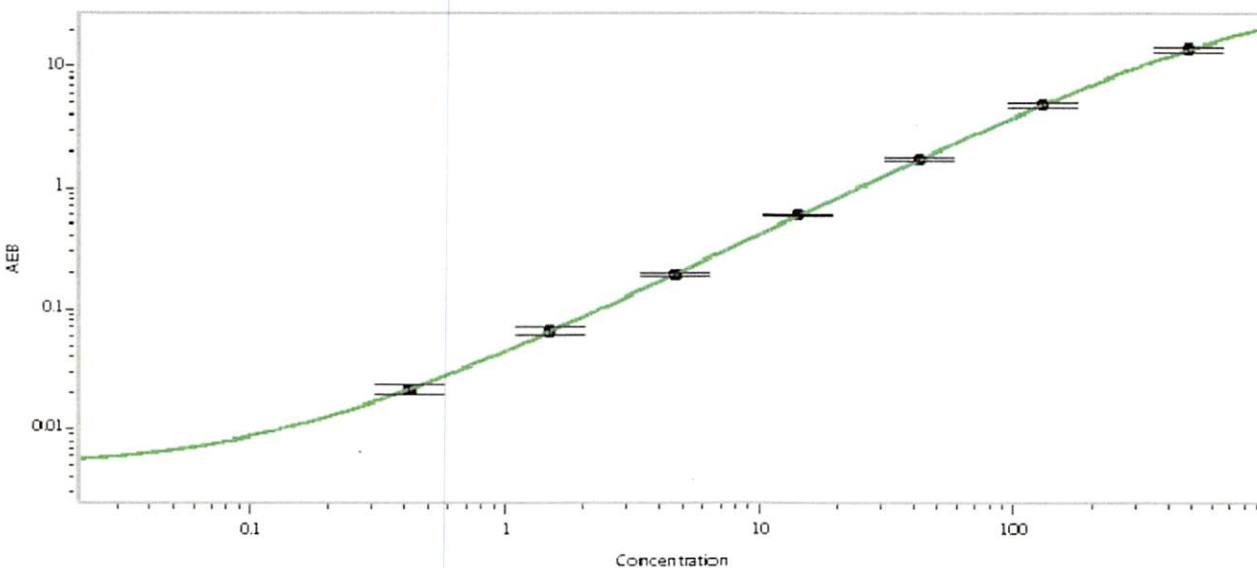
Component	Part Number	Lot Number
Bead Reagent	102246	002705
Detector Reagent	102248	002701
SBG Reagent	102250	002711
Sample Diluent	102252	926609
Calibrators	102255	003501
Controls	102256 - 102257	003601
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	003601	4.70	3.29-4.93
Control 2	003601	179	108-244
Panel 1	922601	17.9	14.1-21.1
Panel 2	922602	6.50	4.94-7.41
Panel 3	922603	36.6	28.4-42.6
Panel 4	922604	164	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* 11 Feb 2020  
 Name Title Signature/Date