

<b>Product Number:</b>	102958
<b>Lot Number:</b>	502303
<b>Expiration:</b>	18-Nov-2020
<b>Platform(s):</b>	HD-X and SR-X

Component	Part Number	Lot Number
<b>Bead Reagent</b>	102949	008405
<b>Detector Reagent</b>	102951	932210
<b>SBG Reagent</b>	102953	005503
<b>Sample Diluent</b>	102954	005501
<b>Calibrators</b>	102960	929405
<b>Controls</b>	102961 & 102962	929406
<b>RGP Reagent</b>	103159	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific



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

<b>Review/Approval</b>	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	<i>03 April 2020</i>
	<b>Name</b>	<b>Title</b>	<b>Signature/Date</b>	

**Release Materials**

**IFN $\gamma$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	929406	2.40	1.01-2.58
Control 2	929406	58.7	28.8-62.2
Panel 1	927401	0.306	0.154-0.440
Panel 2	927703	26.4	13.8-28.6
Panel 3	927704	84.2	46.9-90.3

**IL-6**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	929406	1.38	1.10-1.72
Control 2	929406	38.4	29.2-43.8
Panel 1	927401	5.45	3.87-5.80
Panel 2	927703	13.4	10.0-14.9
Panel 3	927704	53.1	32.3-56.5

**IL-10**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	929406	1.11	0.700-1.27
Control 2	929406	18.9	16.8-26.7
Panel 1	927401	1.13	0.207-1.29
Panel 2	927703	12.6	6.20-14.0
Panel 3	927704	41.9	17.8-45.3

**IL-12p70**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	929406	1.03	0.813-1.22
Control 2	929406	23.2	17.8-26.7
Panel 1	927401	0.216	0.161-0.287
Panel 2	927703	6.92	5.00-7.62
Panel 3	927704	24.4	17.6-26.4

**IL-17A**

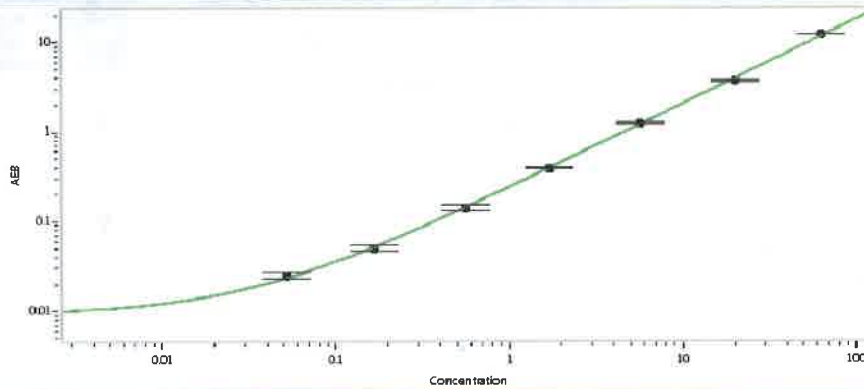
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	929406	0.361	0.305-0.516
Control 2	929406	13.6	10.3-15.4
Panel 1	927401	0.117	0.085-0.196
Panel 2	927703	7.35	5.47-8.45
Panel 3	927704	23.6	16.9-25.4

**TNF $\alpha$**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	929406	1.03	0.809-1.32
Control 2	929406	47.4	34.8-52.2
Panel 1	927401	2.38	1.65-2.65
Panel 2	927703	10.9	7.22-11.6
Panel 3	927704	51.6	29.4-55.2

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

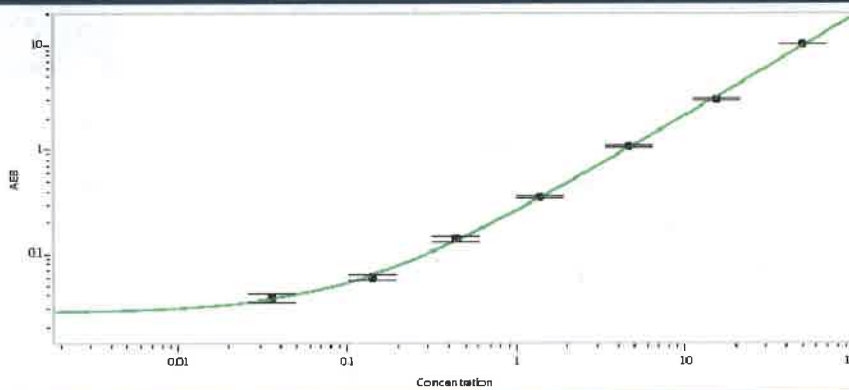
**IFN $\gamma$  Calibration Curve**



**Calibrator Levels (pg/mL)**

<b>A</b>	0.000
<b>B</b>	0.052
<b>C</b>	0.165
<b>D</b>	0.566
<b>E</b>	1.70
<b>F</b>	5.75
<b>G</b>	20.1
<b>H</b>	63.4

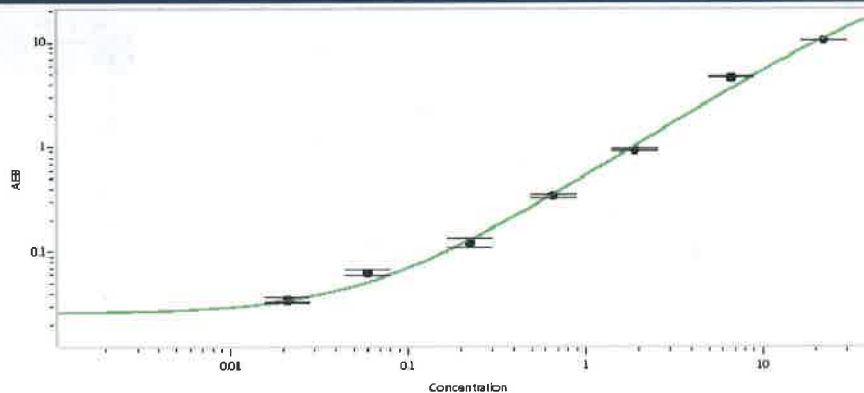
**IL-6 Calibration Curve**



**Calibrator Levels (pg/mL)**

<b>A</b>	0.000
<b>B</b>	0.036
<b>C</b>	0.141
<b>D</b>	0.440
<b>E</b>	1.37
<b>F</b>	4.68
<b>G</b>	15.4
<b>H</b>	50.4

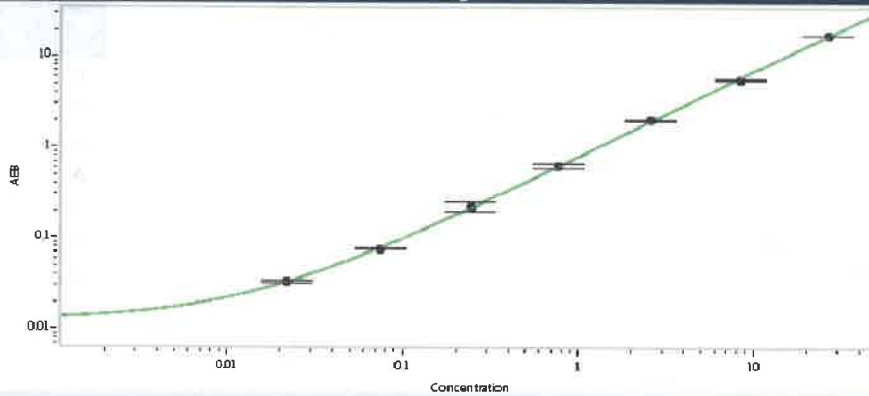
**IL-10 Calibration Curve**



**Calibrator Levels (pg/mL)**

<b>A</b>	0.000
<b>B</b>	0.021
<b>C</b>	0.060
<b>D</b>	0.224
<b>E</b>	0.661
<b>F</b>	1.89
<b>G</b>	6.71
<b>H</b>	22.4

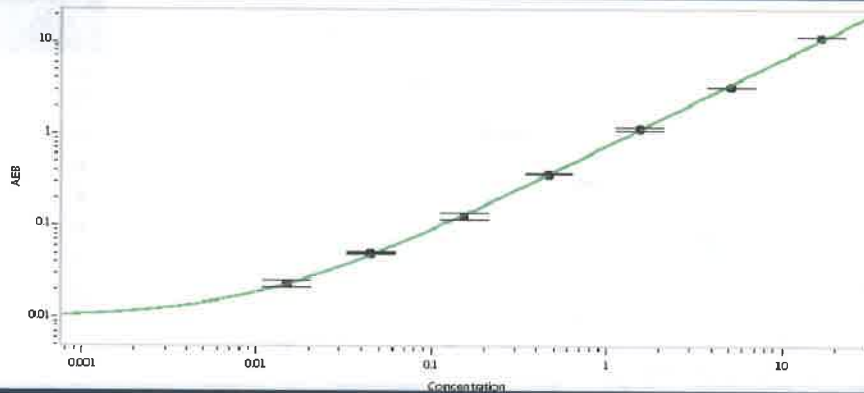
**IL-12p70 Calibration Curve**



**Calibrator Levels  
(pg/mL)**

<b>A</b>	0.000
<b>B</b>	0.022
<b>C</b>	0.075
<b>D</b>	0.243
<b>E</b>	0.766
<b>F</b>	2.59
<b>G</b>	8.39
<b>H</b>	26.4

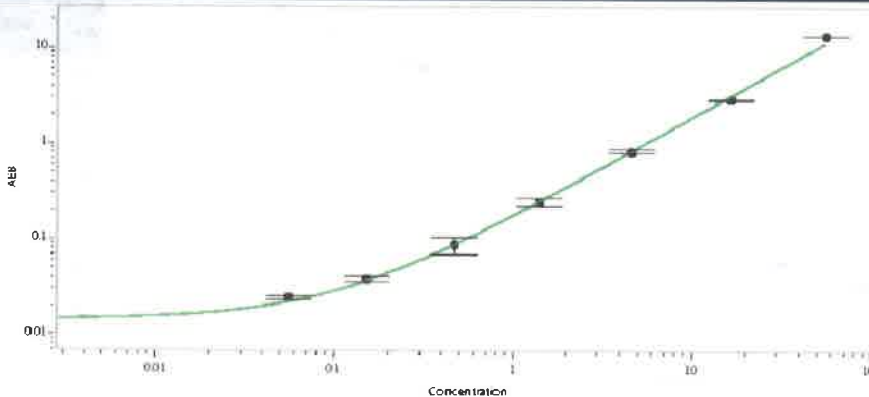
**IL-17A Calibration Curve**



**Calibrator Levels  
(pg/mL)**

<b>A</b>	0.000
<b>B</b>	0.015
<b>C</b>	0.045
<b>D</b>	0.153
<b>E</b>	0.468
<b>F</b>	1.55
<b>G</b>	5.16
<b>H</b>	16.8

**TNFα Calibration Curve**



**Calibrator Levels  
(pg/mL)**

<b>A</b>	0.000
<b>B</b>	0.056
<b>C</b>	0.153
<b>D</b>	0.468
<b>E</b>	1.42
<b>F</b>	4.64
<b>G</b>	16.7
<b>H</b>	56.0