

<b>Product Number:</b>	101643
<b>Lot Number:</b>	504169
<b>Expiration:</b>	18-Jan-2025
<b>Platform(s):</b>	SR-X and HD-1/HD-X

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	101645	327605	N/A
Detector Reagent	101644	327104	N/A
SBG Reagent	101646	327105	327105A
Sample Diluent	102200	319920	319920A
Calibrator Diluent	101648	324512	N/A
Calibrator Concentrate	100670	315305	315305A
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>	N/A <sup>1</sup>



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

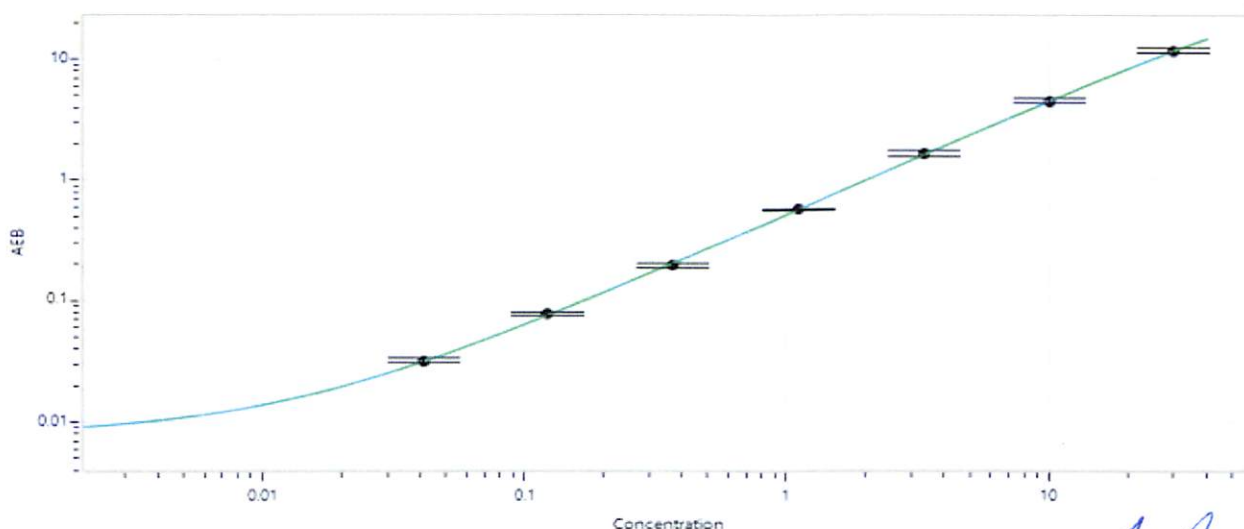
<sup>1</sup> Reagents are not Kit Lot Specific

† Specific to use on the SRX platform.  
Refer to the IL-10 Kit Instructions - SR-X (KI-0036).

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	315305A	1.49	1.20 - 2.00
Control 2	315305A	22.6	18.0 - 30.0
Panel 1	215202	0.395	0.213 - 0.561
Panel 2	215202	0.625	0.425 - 0.738
Panel 3	306507	68.7	47.1 - 71.5

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

**Calibrator Concentrate Value Assignment (pg/mL)** 603



**Calibrator Levels (pg/mL)**

A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Cody Carlson QC Manager  
Name Title

*Cody Carlson* 24 APR 2024  
Signature/Date