

|                 |               |
|-----------------|---------------|
| Product Number: | 101643        |
| Lot Number:     | 501690        |
| Expiration:     | 21-Dec-2019   |
| Platform(s):    | HD-1 and SR-X |

| Component              | Part Number | Lot Number       |
|------------------------|-------------|------------------|
| Bead Reagent           | 101645      | 905812           |
| Detector Reagent       | 101644      | 905808           |
| SBG Reagent (HD-1)     | 101646      | 905809           |
| Sample Diluent         | 102200      | 902931           |
| Calibrator Diluent     | 101648      | 835505           |
| Calibrator Concentrate | 100670      | 833805           |
| SBG Diluent (SRX)†     | 103158      | N/A <sup>1</sup> |
| RGP Reagent            | 103159      | N/A <sup>1</sup> |

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

† Specific for use on the SR-X platform - refer to the Simoa™ Human IL-10 Advantage SR-X Kit Instructions (KI-0036).



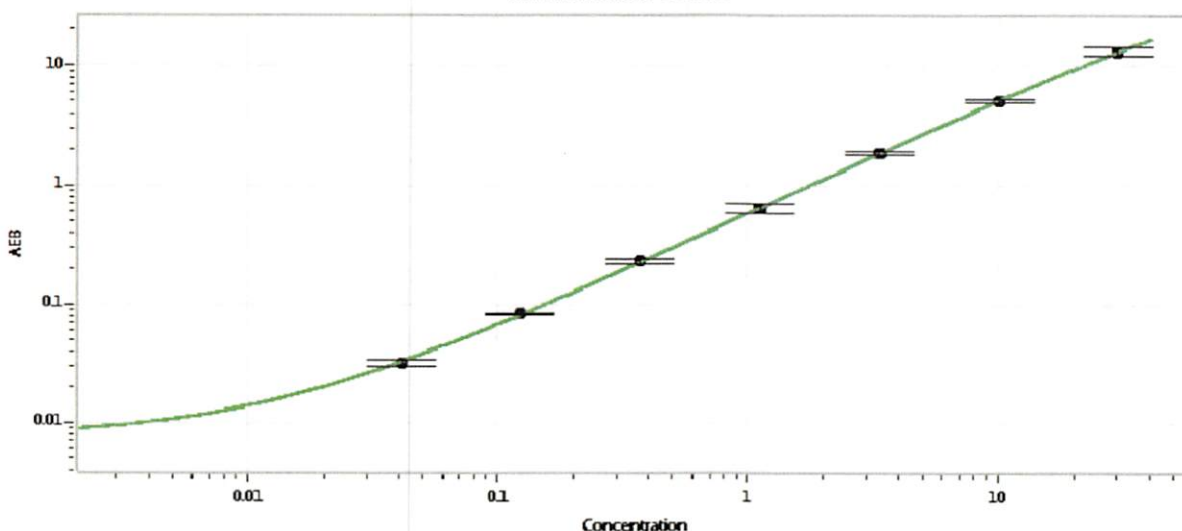
Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Release Materials |              |                |                    | Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges. |
|-------------------|--------------|----------------|--------------------|--------------------------------------------------------------------------------------------------------------------|
|                   | Lot          | Result (pg/mL) | Mean Range (pg/mL) |                                                                                                                    |
| Control 1         | MF-060518-03 | 1.37           | 0.740-1.52         |                                                                                                                    |
| Control 2         | MF-060518-03 | 28.4           | 11.1-32.3          |                                                                                                                    |
| Panel 1           | MF-060718-01 | 0.519          | 0.188-0.563        |                                                                                                                    |
| Panel 2           | MF-060718-02 | 1.17           | 0.536-1.32         |                                                                                                                    |
| Panel 3           | MF-060718-03 | 83.2           | 49.8-89.0          |                                                                                                                    |

#### Calibrator Concentrate Value Assignment

209 pg/mL

Calibration Curve



#### 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

Patrizia Stadler Assoc. Director, QC  
Name Title

P. Stadler 14 Nov 2019  
Signature/Date