

|                 |                  |
|-----------------|------------------|
| Product Number: | 100198           |
| Lot Number:     | 502427           |
| Expiration:     | 15-Jan-2021      |
| Platform(s):    | HD-1/HD-x & SR-X |



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

| Component              | Part Number | Lot Number       |
|------------------------|-------------|------------------|
| Bead Reagent           | 100815      | 016810           |
| Detector Reagent       | 100816      | 016805           |
| SBG Reagent            | 100817      | 016804           |
| Sample Diluent         | 103204      | 016807           |
| Calibrator Diluent     | 100829      | 016803           |
| Calibrator Concentrate | 100828      | 016905           |
| SBG Diluent (SR-X) †   | 103158      | N/A <sup>1</sup> |
| RGP Reagent            | 103159      | N/A <sup>1</sup> |

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

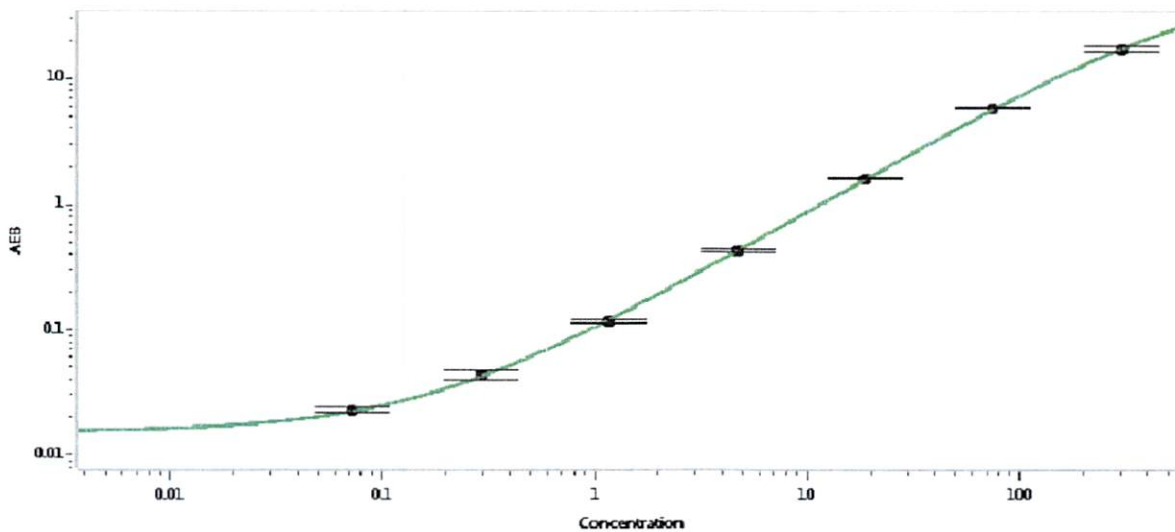
† Specific for use on the SRX platform - refer to the Simoa™ IL-8 SR-X Kit Instructions (KI-0092).

| Release Materials* |              |                |                    |
|--------------------|--------------|----------------|--------------------|
|                    | Lot          | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1          | MF-091217-02 | 9.97           | 7.70-11.6          |
| Control 2          | MF-091217-02 | 186            | 143-220            |
| Panel 1            | MF-061617-01 | 5.05           | 3.17-6.91          |
| Panel 2            | MF-061617-02 | 54.5           | 42.6-63.9          |
| Panel 3            | MF-061617-03 | 311            | 267-459            |

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

|   |       |
|---|-------|
| Calibrator Concentrate Value Assignment (pg/mL) | 5,050 |
|---|-------|

Calibration Curve



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 0.073 |
| C                         | 0.293 |
| D                         | 1.17  |
| E                         | 4.69  |
| F                         | 18.8  |
| G                         | 75.0  |
| H                         | 300   |

Review/Approval: Shivani Goel QC Manager Shivani Goel 02 Jul 2020

Name: Shivani Goel Title: QC Manager Signature/Date: Shivani Goel 02 Jul 2020