

|                 |             |
|-----------------|-------------|
| Product Number: | 103400      |
| Lot Number:     | 502995      |
| Expiration:     | 16-Apr-2022 |
| Platform(s):    | SR-X        |

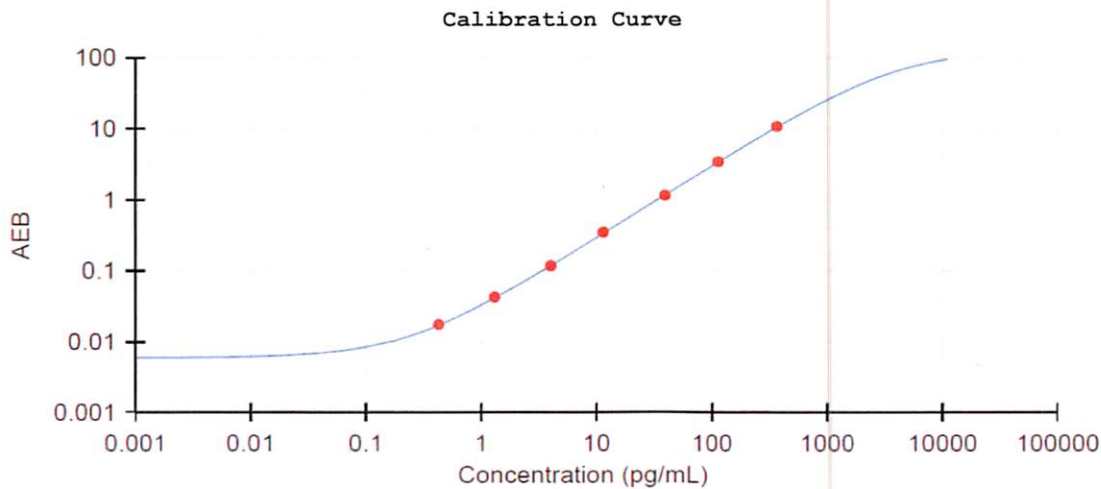


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

| Component        | Part Number | Lot Number       |
|------------------|-------------|------------------|
| Bead Reagent     | 102246      | 115708           |
| Detector Reagent | 102248      | 115809           |
| SBG Reagent      | 102250      | 115810           |
| Sample Diluent   | 103399      | 110609           |
| Calibrators      | 102255      | 115906           |
| Control 1        | 102256      | 115907           |
| Control 2        | 102257      | 115907           |
| RGP Reagent      | 103159      | N/A <sup>1</sup> |

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

| 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         | 115907 | 4.36           | 3.06-5.03          |  |
| Control 2         | 115907 | 252            | 140-314            |  |
| Panel 1           | 008501 | 11.5           | 8.59-20.5          |  |
| Panel 2           | 008501 | 9.00           | 7.39-15.1          |  |
| Panel 3           | 008501 | 66.7           | 43.8-91.8          |  |
| Panel 4           | 009401 | 174            | 128-302            |  |



| Calibrator Levels (pg/mL) |       |
|---------------------------|-------|
| A                         | 0.000 |
| B                         | 0.425 |
| C                         | 1.30  |
| D                         | 3.97  |
| E                         | 11.4  |
| F                         | 38.9  |
| G                         | 113   |
| H                         | 364   |

Review/Approval: Junpil Park QC Supervisor

*[Signature]* 02 Jul 2021