

|                 |             |
|-----------------|-------------|
| Product Number: | 103400      |
| Lot Number:     | 502343      |
| Expiration:     | 24-Mar-2021 |
| Platform(s):    | SR-X        |

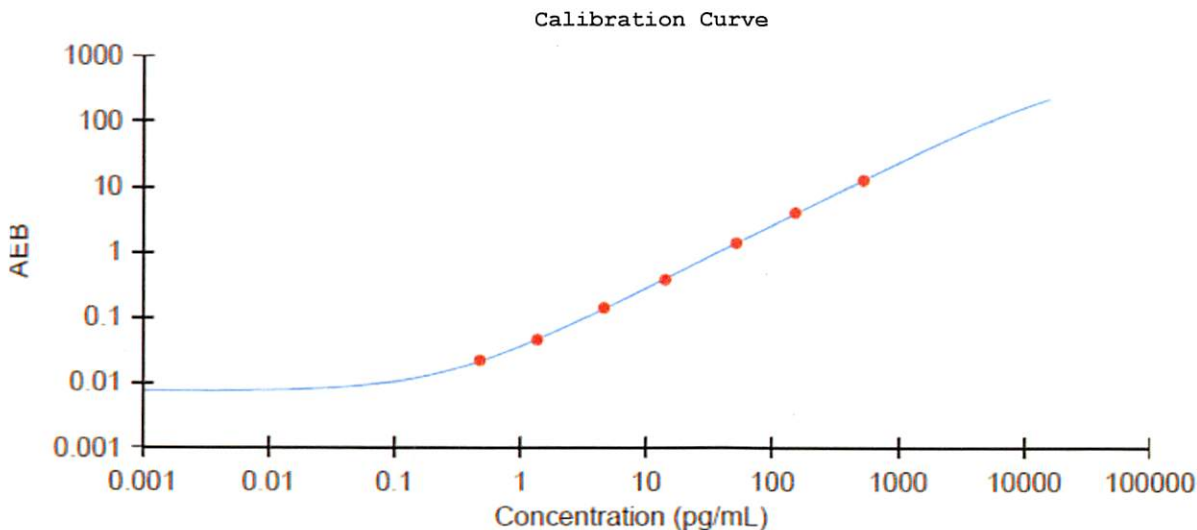


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive Biomarker Detection System

| Component        | Part Number   | Lot Number       |
|------------------|---------------|------------------|
| Bead Reagent     | 102246        | 010001           |
| Detector Reagent | 102248        | 009701           |
| SBG Reagent      | 102250        | 009707           |
| Sample Diluent   | 103399        | 011101           |
| Calibrators      | 102255        | 007001           |
| Controls         | 102256/102257 | 007002           |
| 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         | 007002 | 4.09           | 2.52-5.10          |  |
| Control 2         | 007002 | 169            | 124-186            |  |
| Panel 1           | 922601 | 19.6           | 14.1-21.1          |  |
| Panel 2           | 922602 | 6.41           | 4.94-7.41          |  |
| Panel 3           | 922603 | 40.8           | 28.4-42.6          |  |
| Panel 4           | 922604 | 167            | 129-204            |  |



Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 30 Apr 2020  
Signature/Date