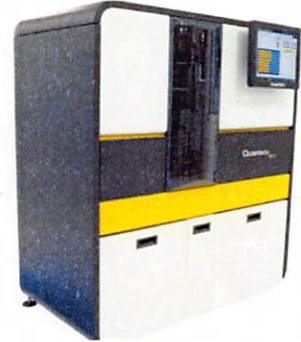


|                 |                  |
|-----------------|------------------|
| Product Number: | 101552           |
| Lot Number:     | 503287           |
| Expiration:     | 14-Oct-2022      |
| Platform(s):    | HD-1/HD-x & SR-X |



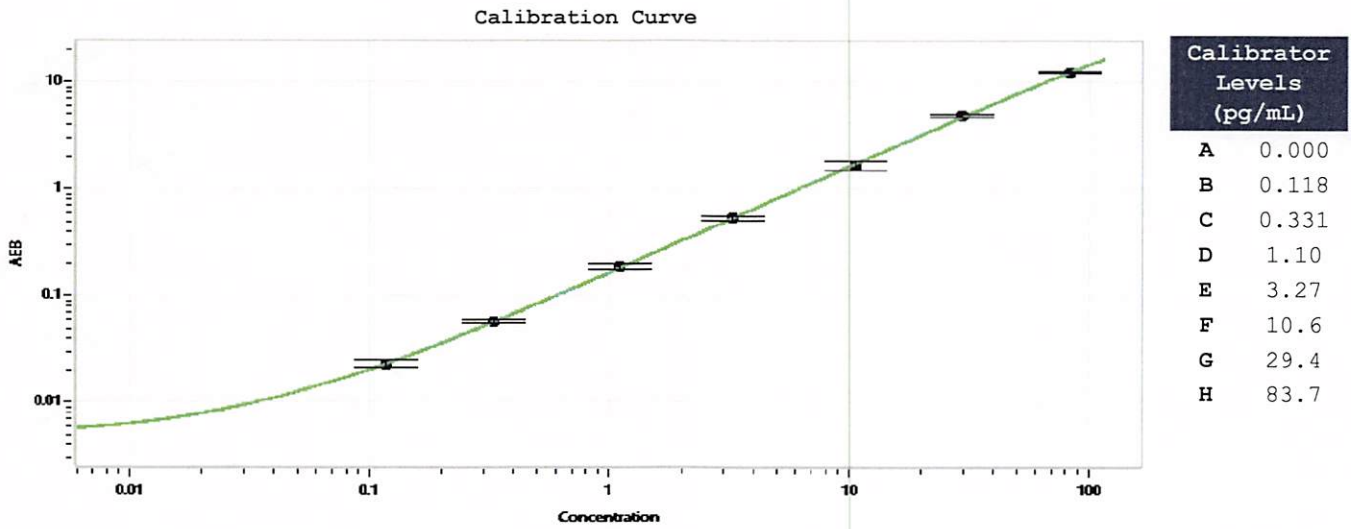
Data below represents results generated on the Simoa™ HD-X Analyzer.

| Component            | Part Number | Lot Number       |
|----------------------|-------------|------------------|
| Bead Reagent         | 101554      | 128704           |
| Detector Reagent     | 101553      | 135505           |
| SBG Reagent          | 101555      | 128804           |
| Sample Diluent       | 101556      | 128805           |
| Calibrators          | 101559      | 127104           |
| Control 1            | 101560      | 127105           |
| Control 2            | 101561      | 127105           |
| 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

| Release Materials* |        |                |                    |
|--------------------|--------|----------------|--------------------|
|                    | Lot    | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1          | 127105 | 1.27           | 0.824-1.53         |
| Control 2          | 127105 | 66.0           | 48.8-73.2          |
| Panel 1            | 117205 | 5.06           | 3.46 - 5.48        |
| Panel 2            | 117205 | 3.11           | 2.03 - 4.76        |
| Panel 3            | 117205 | 22.6           | 17.9 - 26.9        |
| Panel 4            | 119612 | 38.1           | 22.5 - 43.7        |

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



|                 |                            |                           |  |
|-----------------|----------------------------|---------------------------|--|
| Review/Approval | Shivani Goel<br><hr/> Name | QC Manager<br><hr/> Title | Shivani Goel 20 Jan 2022<br><hr/> Signature/Date |
|-----------------|----------------------------|---------------------------|--|