

| | |
|-----------------|--------------------|
| Product Number: | 103714 |
| Lot Number: | 503006 |
| Expiration: | 9-Mar-2022 |
| Platform(s): | HD-1 / HD-X & SR-X |



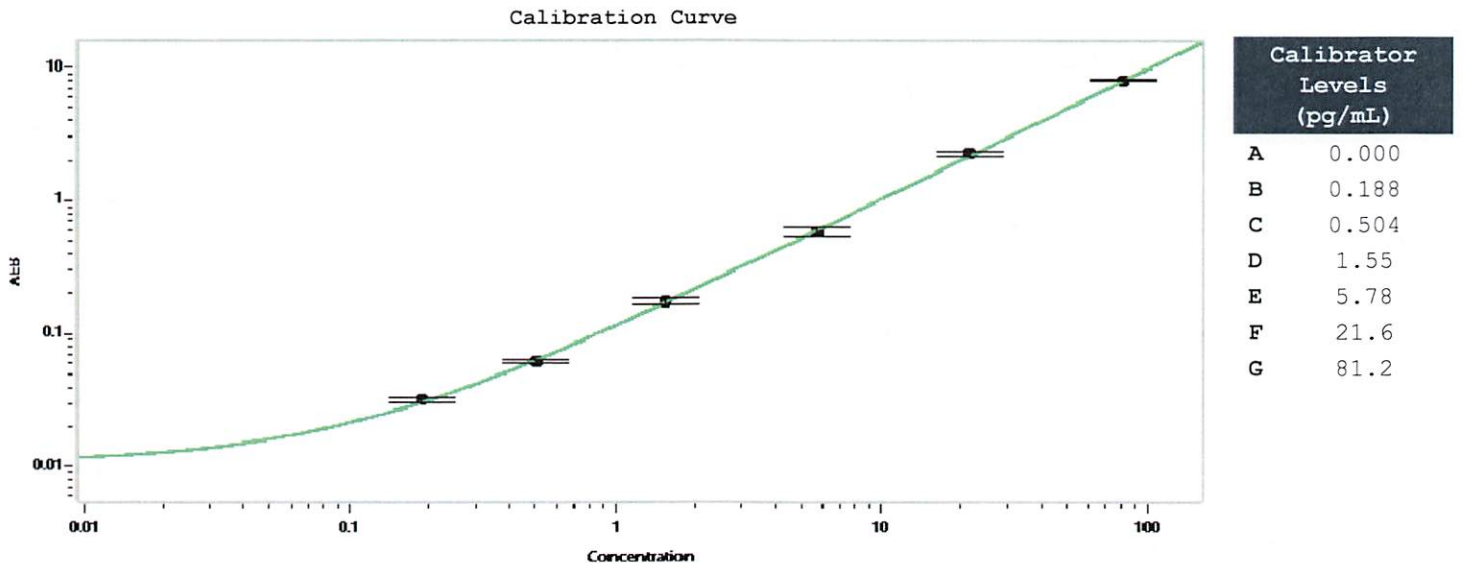
Data below represents results generated on the Simoa™ HD-X Analyzer.

| Component | Part Number | Lot Number |
|----------------------|-------------|------------------|
| Bead Reagent | 103698 | 112903 |
| Detector Reagent | 103700 | 112904 |
| SBG Reagent | 103702 | 112905 |
| Sample Diluent | 103704 | 113013 |
| Reference Calibrator | 103711 | 116604 |
| Control 1 | 103712 | 116605 |
| Control 2 | 103713 | 116605 |
| RGP Reagent | 103159 | N/A ¹ |

¹ RGP is not Kit Lot Specific

| Release Materials* | | | |
|--------------------|--------|----------------|--------------------|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 116605 | 3.27 | 2.55-3.82 |
| Control 2 | 116605 | 112 | 89.3-134 |
| Panel 1 | 013903 | 0.652 | 0.258-1.11 |
| Panel 2 | 013903 | 1.43 | 0.841-2.44 |
| Panel 3 | 013903 | 2.41 | 1.79-4.31 |

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



Review/Approval

Junpil Park
Name

QC Supervisor
Title

25 Jun 2021
Signature/Date