

| | |
|-----------------|------------|
| Product Number: | 103806 |
| Lot Number: | 503175 |
| Expiration: | 6-Oct-2022 |
| Bulk Kit | |
| Product Number: | 103872 |
| Lot Number: | 503176 |
| Platform(s): | HD-X |



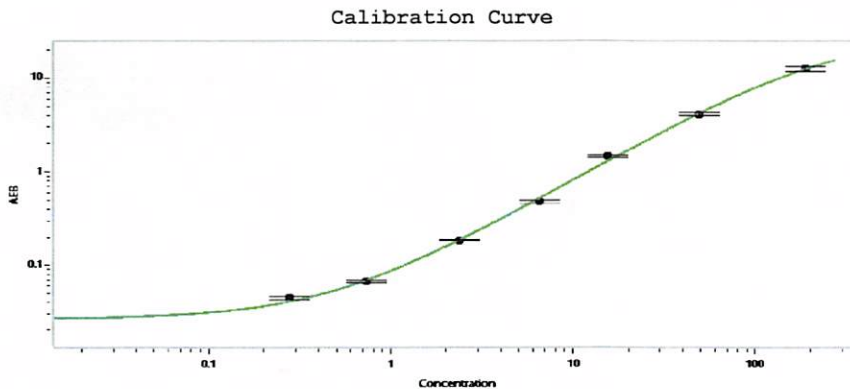
Data below represents results generated on the Simoa® HD-X Analyzer.

| Component | Part Number | Lot Number |
|------------------|-------------|------------------|
| Bead Reagent | 103794 | 127906 |
| Detector Reagent | 103796 | 201807 |
| SBG Reagent | 103798 | 128007 |
| Sample Diluent | 103800 | 128011 |
| Calibrators | 103803 | 207302 |
| Control 1 | 103804 | 207303 |
| Control 2 | 103805 | 207303 |
| RGP Reagent | 103159 | N/A ¹ |

¹ RGP is not Kit Lot Specific

| Release Materials* | | | |
|--------------------|--------|----------------|--------------------|
| Sample | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | 207303 | 7.29 | 5.51-8.26 |
| Control 2 | 207303 | 124 | 92.4-163 |
| Panel 1 | 021305 | 3.39 | 2.55-4.27 |
| Panel 2 | 021111 | 63.8 | 42.2-90.4 |
| Panel 3 | 021111 | 131 | 89.1-202 |

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



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.278 |
| C | 0.729 |
| D | 2.35 |
| E | 6.49 |
| F | 15.4 |
| G | 48.8 |
| H | 189 |

Review/Approval Shivani Goel QC Manager Shivani Goel 14 Apr 2022
 Name Title Signature/Date