

| | |
|-----------------|------------------|
| Product Number: | 103337 |
| Lot Number: | 504001 |
| Expiration: | 4-Oct-2024 |
| Platform(s): | HD-1/HD-X & SR-X |



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

| Component | Item Number | Lot Number |
|------------------------|-------------|------------------|
| Bead Reagent | 103334 | 328411 |
| Detector Reagent | 101091 | 328503 |
| SBG Reagent | 101101 | 328504 |
| Sample Diluent | 103336 | 328505 |
| Calibrator Diluent | 101095 | 327712 |
| Calibrator Concentrate | 101093 | 328513 |
| RGP Reagent | 103159 | N/A ¹ |

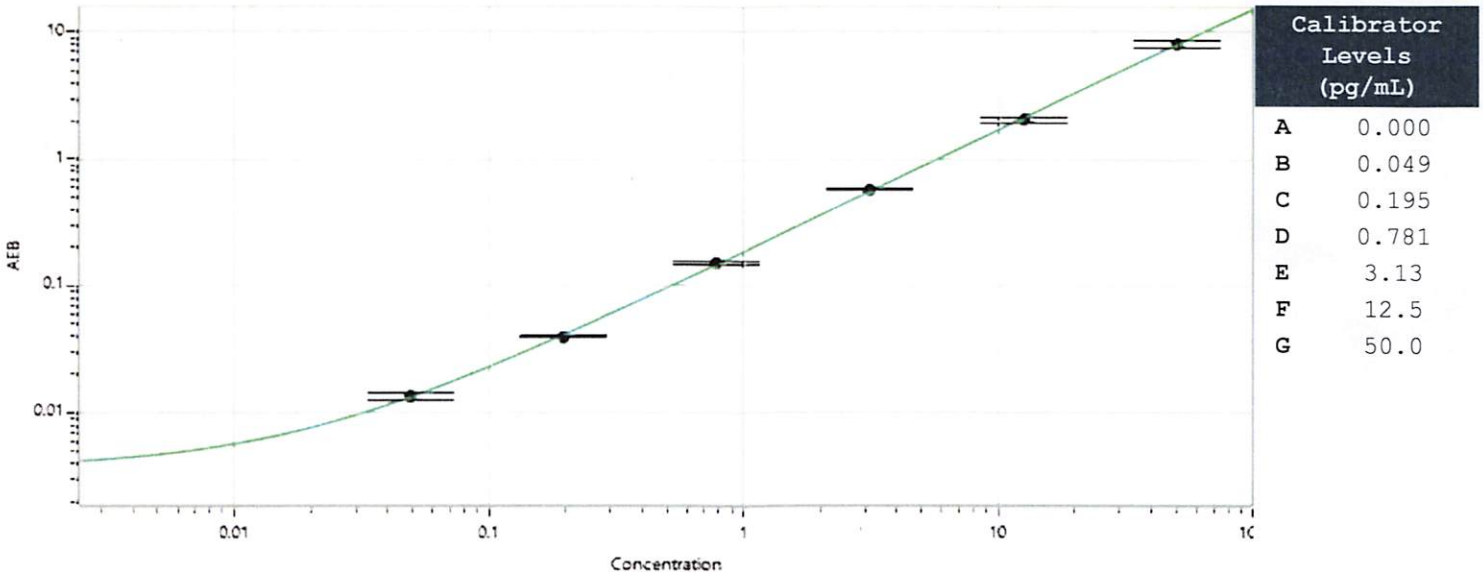
¹ RGP is not Kit Lot Specific

| Release Materials* | | | |
|--------------------|--------|----------------|---------------------|
| | Lot | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1 | 328513 | 0.786 | 0.750 - 1.25 |
| Control 2 | 328513 | 22.6 | 15.0 - 25.0 |
| Panel 1 | 211906 | 1.94 | 1.26 - 2.41 |
| Panel 2 | 211906 | 3.40 | 2.37 - 4.51 |
| Panel 3 | 211906 | 33.6 | 21.0 - 45.2 |

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

| | |
|---|-------|
| Calibrator Concentrate Value Assignment (pg/mL) | 2,499 |
|---|-------|

Calibration Curve



Review/Approval

Sonia Khillan
Name

QC
Title

Sonia Khillan 03NOV2023
Signature/Date