

| | |
|-----------------|------------------|
| Product Number: | 103293 |
| Lot Number: | 503401 |
| Expiration: | 24-Mar-2023 |
| 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 | 103281 | 208906 |
| Detector Reagent | 103283 | 208909 |
| SBG Reagent | 103285 | 208910 |
| Sample Diluent | 103286 | 217221 |
| Calibrators Diluent | 103288 | 208311 |
| Calibrator Concentrate | 103287 | 233612 |
| SBG Diluent (SR-X) † | 103158 | N/A ¹ |
| RGP Reagent | 103159 | N/A ¹ |

¹ RGP is not Kit Lot Specific

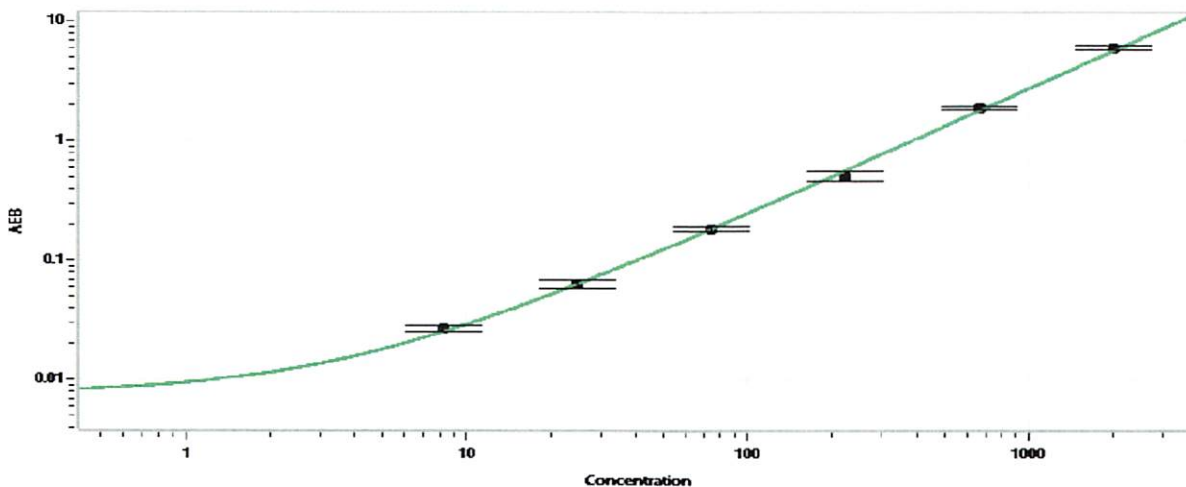
† Specific for use on the SR-X platform - refer to the Simoa™ TDP-43 SR-X Kit Instructions (KI-0132).

| Release Materials* | | | |
|--------------------|--------|----------------|---------------------|
| | Lot | Result (pg/mL) | Spec. Range (pg/mL) |
| Control 1 | 106304 | 83.2 | 70.8 - 127 |
| Control 2 | 106304 | 929 | 510 - 1320 |
| Panel 2 | 231407 | 41.3 | 18.9 - 71.0 |
| Panel 3 | 232608 | 175 | 122 - 239 |

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

| | |
|---|--------|
| Calibrator Concentrate Value Assignment (pg/mL) | 22,287 |
|---|--------|

Calibration Curve



Calibrator Levels (pg/mL)

| | |
|---|-------|
| A | 0.000 |
| B | 8.23 |
| C | 24.7 |
| D | 74.1 |
| E | 222 |
| F | 667 |
| G | 2000 |

Review/Approval

Sonia Khillan
Name

QC
Title

Sonia Della / 04 JAN 2023
Signature/Date