

| | |
|-----------------|------------------|
| Product Number: | 100199 |
| Lot Number: | 502080 |
| Expiration: | 30-Mar-2020 |
| Platform(s): | HD-1/HD-X & SR-X |

| Component | Part Number | Lot Number |
|------------------------|-------------|------------------|
| Bead Reagent | 101045 | 929003 |
| Detector Reagent | 101047 | 928812 |
| SBG Reagent | 101057 | 928811 |
| Sample Diluent | 101055 | 908925 |
| Calibrator Diluent | 101051 | 928012 |
| Calibrator Concentrate | 101049 | 81302 |
| SBG Diluent (SR-X)† | 103158 | N/A ¹ |
| RGP Reagent | 103159 | N/A ¹ |

¹ Reagents are not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ C-Peptide Advantage SR-X Kit Instructions (KI-0109).



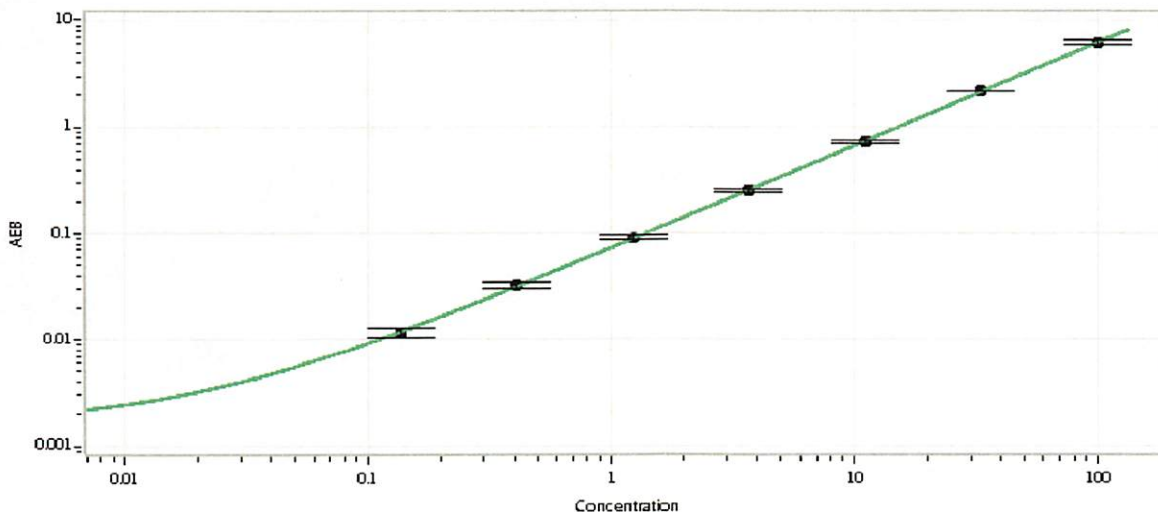
Data below represents results generated on the Simoa™ HD-1 Analyzer.

| Release Materials* | | | |
|--------------------|--------------|----------------|--------------------|
| | Lot | Result (pg/mL) | Mean Range (pg/mL) |
| Control 1 | MF-032017-04 | 1.65 | 1.32-1.97 |
| Control 2 | MF-032017-04 | 50.9 | 42.4-63.5 |
| Panel 1 | 818704 | 129 | 91.6-137 |
| Panel 2 | MF-060118-02 | 1885 | 1458-2290 |
| Panel 3 | MF-060118-03 | 2828 | 2193-3290 |

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

| | |
|---|------|
| Calibrator Concentrate Value Assignment (pg/mL) | 1832 |
|---|------|

Calibration Curve



| Calibrator Levels (pg/mL) | |
|---------------------------|-------|
| A | 0.000 |
| B | 0.137 |
| C | 0.412 |
| D | 1.23 |
| E | 3.70 |
| F | 11.1 |
| G | 33.3 |
| H | 100 |

Review/Approval

Patrizia Stadler Assoc. Director, QC

P. Stoollev

13 NOV 2019

Name

Title

Signature/Date