



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 5

Catalog Number: C0065-2

Lot Number: A00U0009

Expiration Date: 31 MAY 2020

Storage Condition: 2-8°C

| Analyte Description | Amount of Analyte in Calibrator (pg) | Stock Concentration of Reconstituted Calibrator (Vol = 250 µL) (pg/mL) | Concentration of Standard 1 (after 5-fold Dilution) (pg/mL) |
|---------------------|--------------------------------------|--|---|
| ms IFN- γ | 3 600 | 14 400 | 2 880 |
| ms IL-1 β | 16 250 | 65 000 | 13 000 |
| ms IL-2 | 13 625 | 54 500 | 10 900 |
| ms IL-4 | 12 450 | 49 800 | 9 960 |
| ms IL-5 | 3 325 | 13 300 | 2 660 |
| ms IL-6 | 20 000 | 80 000 | 16 000 |
| ms KC | 2 800 | 11 200 | 2 240 |
| ms IL-10 | 28 500 | 114 000 | 22 800 |
| ms IL-12p70 | 106 750 | 427 000 | 85 400 |
| ms TNF- α | 7 688 | 30 750 | 6 150 |
| ms GM-CSF | 1 263 | 5 050 | 1 010 |
| ms IL-13 | 27 250 | 109 000 | 21 800 |
| ms EPO | 15 500 | 62 000 | 12 400 |
| ms VEGF-A | 14 500 | 58 000 | 11 600 |

Statement: The above calibrator lot is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|------------------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 19 JUL 2017 |