



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 1

**Catalog Number:** C0060-2

**Lot Number:** A00U0027

**Expiration Date:** 30 SEP 2021

**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) |
|---------------------|--------------------------------------|--|---|
| Human IFN- $\gamma$ | 22 250                               | 89 000   | 17 800  |
| Human IL-1 $\beta$  | 4 950                                | 19 800   | 3 960   |
| Human IL-2          | 2 488                                | 9 950  | 1 990   |
| Human IL-4          | 2 738                                | 10 950   | 2 190   |
| Human IL-6          | 2 638                                | 10 550   | 2 110   |
| Human IL-8          | 2 850                                | 11 400   | 2 280   |
| Human IL-10         | 4 825                                | 19 300   | 3 860   |
| Human IL-12p70      | 7 088                                | 28 350   | 5 670   |
| Human IL-13         | 2 400                                | 9 600  | 1 920   |
| Human TNF- $\alpha$ | 4 425                                | 17 700   | 3 540   |
| Human IL-5          | 5 163                                | 20 650   | 4 130   |
| Human GM-CSF        | 12 250                               | 49 000   | 9 800   |
| Human IL-17A        | 30 500                               | 122 000  | 24 400  |
| Human VEGF-A        | 6 125                                | 24 500   | 4 900   |

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

|                        | Name         | Function          | Signature           | Date        |
|------------------------|--------------|-------------------|---------------------|-------------|
| <b>Review/Approval</b> | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 17 JUL 2017 |