



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0096

Expiration Date: 29 FEB 2024

Storage Condition: 2-8°C

| Analyte Description | Amount of Analyte in Calibrator | Units | Stock Concentration of Reconstituted Calibrator (Vol = 250 µL) | Units |
|---------------------|---------------------------------|-------|--|-------|
| hu IFN- γ | 31 375 | pg | 125 500 | pg/mL |
| hu IL-1 β | 4 575 | pg | 18 300 | pg/mL |
| hu IL-2 | 2 563 | pg | 10 250 | pg/mL |
| hu IL-4 | 2 425 | pg | 9 700 | pg/mL |
| hu IL-6 | 2 563 | pg | 10 250 | pg/mL |
| hu IL-8 | 2 538 | pg | 10 150 | pg/mL |
| hu IL-10 | 4 613 | pg | 18 450 | pg/mL |
| hu IL-12p70 | 8 088 | pg | 32 350 | pg/mL |
| hu IL-13 | 2 888 | pg | 11 550 | pg/mL |
| hu TNF- α | 4 300 | pg | 17 200 | pg/mL |
| hu IL-5 | 4 950 | pg | 19 800 | pg/mL |
| hu GM-CSF | 12 188 | pg | 48 750 | pg/mL |
| hu IL-17A | 49 250 | pg | 197 000 | pg/mL |
| hu VEGF-A | 6 813 | pg | 27 250 | pg/mL |

To calculate the **Calibrator Standard 1 (top of curve) concentration**, please refer to the dilution-factor used in the **Prepare Calibrator Standards** section of the respective product insert. Divide the stock concentration of the reconstituted calibrator by this dilution factor to generate the working calibrator standard 1 concentration.

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 | | 03 APR 2019 |