



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0030

Expiration Date: 28 FEB 2022

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 | 25 750 | pg | 103 000 | pg/mL |
| hu IL-1 β | 4 925 | pg | 19 700 | pg/mL |
| hu IL-2 | 2 425 | pg | 9 700 | pg/mL |
| hu IL-4 | 2 213 | pg | 8 850 | pg/mL |
| hu IL-6 | 2 700 | pg | 10 800 | pg/mL |
| hu IL-8 | 2 450 | pg | 9 800 | pg/mL |
| hu IL-10 | 4 925 | pg | 19 700 | pg/mL |
| hu IL-12p70 | 7 150 | pg | 28 600 | pg/mL |
| hu IL-13 | 2 425 | pg | 9 700 | pg/mL |
| hu TNF- α | 4 538 | pg | 18 150 | pg/mL |
| hu IL-5 | 5 213 | pg | 20 850 | pg/mL |
| hu GM-CSF | 12 625 | pg | 50 500 | pg/mL |
| hu IL-17A | 22 750 | pg | 91 000 | pg/mL |
| hu VEGF-A | 7 113 | pg | 28 450 | 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 | <i>Dawn Stollar</i> | 29 AUG 2018 |