



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 3

Catalog Number: C0062-2

Lot Number: A00U0154

Expiration Date: 30 APR 2025

Storage Temperature: 2-8°C

Intended Assay Species: Human

Concentration Assignment:

| Analyte | Amount of Analyte in Calibrator | Unit of Measure | Stock Concentration of Reconstituted Calibrator | Unit of Measure |
|------------------|---------------------------------|-----------------|---|-----------------|
| IL-1 α | 7800 | pg | 31200 | pg/mL |
| IL-7 | 7738 | pg | 30950 | pg/mL |
| IL-12/IL-23p40 | 26500 | pg | 106000 | pg/mL |
| IL-15 | 3663 | pg | 14650 | pg/mL |
| IL-16 | 31625 | pg | 126500 | pg/mL |
| TNF- β | 4713 | pg | 18850 | pg/mL |
| G-CSF | 23375 | pg | 93500 | pg/mL |
| IFN- α 2a | 42625 | pg | 170500 | pg/mL |
| IL-18 | 48625 | pg | 194500 | pg/mL |
| TPO | 38625 | pg | 154500 | pg/mL |

Reconstitution Volume: 250 μ L

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.

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|------------------------|---------------|----------|-----------|-------------|
| Review/Approval | Karen Hamilla | Quality | | 19 MAY 2020 |