



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 1 (human) Calibrator Blend

Catalog Number: C0050-2

Lot Number: A0080318

Expiration Date: 31 JAN 2027

Storage Temperature: 2-8°C

Concentration Assignment:

| Analyte | Stock Concentration of Reconstituted Calibrator | Unit of Measure |
|----------------|---|-----------------|
| GM-CSF | 1180 | pg/mL |
| IL-1 α | 418 | pg/mL |
| IL-5 | 853 | pg/mL |
| IL-7 | 782 | pg/mL |
| IL-12/IL-23p40 | 3050 | pg/mL |
| IL-15 | 763 | pg/mL |
| IL-16 | 2410 | pg/mL |
| IL-17A | 5660 | pg/mL |
| TNF- β | 745 | pg/mL |
| VEGF | 1040 | pg/mL |

Reconstitution Volume: 1000 μ L

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

| | Name | Function | Signature | Date |
|------------------------|---------------|----------|-----------|-------------|
| Review/Approval | Karen Hamilla | Quality | | 02 JUN 2022 |