



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 3

**Catalog Number:** C0062-2

**Lot Number:** A00U0147

**Expiration Date:** 31 MAR 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 $\alpha$	6988	pg	27950	pg/mL
IL-7	7713	pg	30850	pg/mL
IL-12/IL-23p40	28750	pg	115000	pg/mL
IL-15	4088	pg	16350	pg/mL
IL-16	33250	pg	133000	pg/mL
TNF- $\beta$	4813	pg	19250	pg/mL
G-CSF	22375	pg	89500	pg/mL
IFN- $\alpha$ 2a	44375	pg	177500	pg/mL
IL-18	52500	pg	210000	pg/mL
TPO	36000	pg	144000	pg/mL

**Reconstitution Volume:** 250  $\mu$ 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		20 MAR 2020