



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 3

**Catalog Number:** C0062-2

**Lot Number:** A00U0179

**Expiration Date:** 30 NOV 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 $\alpha$	7363	pg	29450	pg/mL
IL-7	8313	pg	33250	pg/mL
IL-12/IL-23p40	28750	pg	115000	pg/mL
IL-15	3950	pg	15800	pg/mL
IL-16	24875	pg	99500	pg/mL
TNF- $\beta$	5063	pg	20250	pg/mL
G-CSF	24125	pg	96500	pg/mL
IFN- $\alpha$ 2a	44625	pg	178500	pg/mL
IL-18	46250	pg	185000	pg/mL
TPO	39750	pg	159000	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
<b>Review/Approval</b>	Karen Hamilla	Quality		25 JAN 2021