



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 3

**Catalog Number:** C0062-2

**Lot Number:** A00U0115

**Expiration Date:** 30 JUN 2024

**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 $\alpha$	6 538	pg	26 150	pg/mL
IL-7	7 363	pg	29 450	pg/mL
IL-12/IL-23p40	31 875	pg	127 500	pg/mL
IL-15	4 163	pg	16 650	pg/mL
IL-16	34 125	pg	136 500	pg/mL
TNF- $\beta$	4 288	pg	17 150	pg/mL
G-CSF	23 000	pg	92 000	pg/mL
IFN- $\alpha$ 2a	44 500	pg	178 000	pg/mL
IL-18	50 625	pg	202 500	pg/mL
TPO	43 250	pg	173 000	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		16 JUL 2019