

**CERTIFICATE OF ANALYSIS**

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0395
Expiration Date: 30 SEP 2029
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
IFN- γ	40375	pg	161500	pg/mL
IL-1 β	5575	pg	22300	pg/mL
IL-2	2387.5	pg	9550	pg/mL
IL-4	2337.5	pg	9350	pg/mL
IL-6	2512.5	pg	10050	pg/mL
IL-8	2812.5	pg	11250	pg/mL
IL-10	5062.5	pg	20250	pg/mL
IL-12p70	7350	pg	29400	pg/mL
IL-13	2325	pg	9300	pg/mL
TNF- α	4762.5	pg	19050	pg/mL
IL-5	5687.5	pg	22750	pg/mL
GM-CSF	11250	pg	45000	pg/mL
IL-17A	24750	pg	99000	pg/mL
VEGF-A	3375	pg	13500	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	Michelle Likiam	Quality		11 NOV 2024