



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0300
Expiration Date: 31 OCT 2027
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	38500	pg	154000	pg/mL
IL-1 β	6300	pg	25200	pg/mL
IL-2	2600	pg	10400	pg/mL
IL-4	2587.5	pg	10350	pg/mL
IL-6	2462.5	pg	9850	pg/mL
IL-8	3175	pg	12700	pg/mL
IL-10	5075	pg	20300	pg/mL
IL-12p70	7437.5	pg	29750	pg/mL
IL-13	2637.5	pg	10550	pg/mL
TNF- α	3800	pg	15200	pg/mL
IL-5	5775	pg	23100	pg/mL
GM-CSF	12875	pg	51500	pg/mL
IL-17A	38500	pg	154000	pg/mL
VEGF-A	3225	pg	12900	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 was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		12 DEC 2022