



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0322
Expiration Date: 31 MAR 2028
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	36875	pg	147500	pg/mL
IL-1 β	6437.5	pg	25750	pg/mL
IL-2	2787.5	pg	11150	pg/mL
IL-4	2912.5	pg	11650	pg/mL
IL-6	2800	pg	11200	pg/mL
IL-8	3100	pg	12400	pg/mL
IL-10	5662.5	pg	22650	pg/mL
IL-12p70	8437.5	pg	33750	pg/mL
IL-13	2887.5	pg	11550	pg/mL
TNF- α	4850	pg	19400	pg/mL
IL-5	5862.5	pg	23450	pg/mL
GM-CSF	13750	pg	55000	pg/mL
IL-17A	37000	pg	148000	pg/mL
VEGF-A	6225	pg	24900	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		02 MAY 2023