



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0328
Expiration Date: 31 MAY 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	37500	pg	150000	pg/mL
IL-1 β	6062.5	pg	24250	pg/mL
IL-2	2587.5	pg	10350	pg/mL
IL-4	2025	pg	8100	pg/mL
IL-6	2437.5	pg	9750	pg/mL
IL-8	2712.5	pg	10850	pg/mL
IL-10	4912.5	pg	19650	pg/mL
IL-12p70	7137.5	pg	28550	pg/mL
IL-13	2250	pg	9000	pg/mL
TNF- α	3650	pg	14600	pg/mL
IL-5	5325	pg	21300	pg/mL
GM-CSF	12875	pg	51500	pg/mL
IL-17A	34875	pg	139500	pg/mL
VEGF-A	3737.5	pg	14950	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		13 JUN 2023