



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 1
Catalog Number: C0060-2
Lot Number: A00U0317
Expiration Date: 29 FEB 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	40625	pg	162500	pg/mL
IL-1 β	6625	pg	26500	pg/mL
IL-2	2712.5	pg	10850	pg/mL
IL-4	2575	pg	10300	pg/mL
IL-6	2400	pg	9600	pg/mL
IL-8	2487.5	pg	9950	pg/mL
IL-10	4975	pg	19900	pg/mL
IL-12p70	7212.5	pg	28850	pg/mL
IL-13	2662.5	pg	10650	pg/mL
TNF- α	3600	pg	14400	pg/mL
IL-5	5575	pg	22300	pg/mL
GM-CSF	12875	pg	51500	pg/mL
IL-17A	40375	pg	161500	pg/mL
VEGF-A	5087.5	pg	20350	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		23 MAR 2023