



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 1  
**Catalog Number:** C0060-2  
**Lot Number:** A00U0335  
**Expiration Date:** 31 JUL 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- $\gamma$	35875	pg	143500	pg/mL
IL-1 $\beta$	6625	pg	26500	pg/mL
IL-2	2700	pg	10800	pg/mL
IL-4	2450	pg	9800	pg/mL
IL-6	2550	pg	10200	pg/mL
IL-8	3012.5	pg	12050	pg/mL
IL-10	5087.5	pg	20350	pg/mL
IL-12p70	7537.5	pg	30150	pg/mL
IL-13	2500	pg	10000	pg/mL
TNF- $\alpha$	3575	pg	14300	pg/mL
IL-5	5400	pg	21600	pg/mL
GM-CSF	13250	pg	53000	pg/mL
IL-17A	41500	pg	166000	pg/mL
VEGF-A	3962.5	pg	15850	pg/mL

**Reconstitution Volume:** 250  $\mu$ 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
<b>Review/Approval</b>	Karen Hamilla	Quality		23 AUG 2023