



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 1  
**Catalog Number:** C0060-2  
**Lot Number:** A00U0327  
**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- $\gamma$	42375	pg	169500	pg/mL
IL-1 $\beta$	6362.5	pg	25450	pg/mL
IL-2	2837.5	pg	11350	pg/mL
IL-4	2475	pg	9900	pg/mL
IL-6	2625	pg	10500	pg/mL
IL-8	2950	pg	11800	pg/mL
IL-10	5337.5	pg	21350	pg/mL
IL-12p70	8162.5	pg	32650	pg/mL
IL-13	2687.5	pg	10750	pg/mL
TNF- $\alpha$	4662.5	pg	18650	pg/mL
IL-5	5737.5	pg	22950	pg/mL
GM-CSF	13500	pg	54000	pg/mL
IL-17A	37250	pg	149000	pg/mL
VEGF-A	4800	pg	19200	pg/mL

**Reconstitution Volume:** 250 $\mu$ L  $\mu$ 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		20 JUN 2023