



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 5  
**Catalog Number:** C0065-2  
**Lot Number:** A00U0309  
**Expiration Date:** 29 FEB 2028  
**Storage Temperature:** 2-8°C  
**Intended Assay Species:** Mouse

**Concentration Assignment:**

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	3512.5	pg	14050	pg/mL
IL-1 $\beta$	16875	pg	67500	pg/mL
IL-2	11300	pg	45200	pg/mL
IL-4	12175	pg	48700	pg/mL
IL-5	3712.5	pg	14850	pg/mL
IL-6	20000	pg	80000	pg/mL
KC/GRO	2537.5	pg	10150	pg/mL
IL-10	27000	pg	108000	pg/mL
IL-12p70	124125	pg	496500	pg/mL
TNF- $\alpha$	7475	pg	29900	pg/mL
GM-CSF	1325	pg	5300	pg/mL
IL-13	21625	pg	86500	pg/mL
EPO	16000	pg	64000	pg/mL
VEGF-A	12475	pg	49900	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		03 MAY 2023