



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 5  
**Catalog Number:** C0065-2  
**Lot Number:** A00U0325  
**Expiration Date:** 30 APR 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- $\gamma$	3925	pg	15700	pg/mL
IL-1 $\beta$	18500	pg	74000	pg/mL
IL-2	12625	pg	50500	pg/mL
IL-4	13125	pg	52500	pg/mL
IL-5	4025	pg	16100	pg/mL
IL-6	22375	pg	89500	pg/mL
KC/GRO	2712.5	pg	10850	pg/mL
IL-10	32500	pg	130000	pg/mL
IL-12p70	137500	pg	550000	pg/mL
TNF- $\alpha$	8562.5	pg	34250	pg/mL
GM-CSF	1375	pg	5500	pg/mL
IL-13	26500	pg	106000	pg/mL
EPO	19000	pg	76000	pg/mL
VEGF-A	14125	pg	56500	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>	Xiomara Talavera	Quality		25 MAY 2023