



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 5  
**Catalog Number:** C0065-2  
**Lot Number:** A00U0285  
**Expiration Date:** 30 NOV 2027  
**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- $\gamma$	3762.5	pg	15050	pg/mL
IL-1 $\beta$	16750	pg	67000	pg/mL
IL-2	11375	pg	45500	pg/mL
IL-4	12400	pg	49600	pg/mL
IL-5	3812.5	pg	15250	pg/mL
IL-6	21000	pg	84000	pg/mL
KC/GRO	2600	pg	10400	pg/mL
IL-10	29000	pg	116000	pg/mL
IL-12p70	133750	pg	535000	pg/mL
TNF- $\alpha$	8150	pg	32600	pg/mL
GM-CSF	1337.5	pg	5350	pg/mL
IL-13	24625	pg	98500	pg/mL
EPO	16625	pg	66500	pg/mL
VEGF-A	13125	pg	52500	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		29 DEC 2022



Page 2 of 2

**CERTIFICATE OF ANALYSIS**

COA-17612-A

QA-FM-184-F