



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 5  
**Catalog Number:** C0065-2  
**Lot Number:** A00U0384  
**Expiration Date:** 31 DEC 2029  
**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- $\gamma$	3200	pg	12800	pg/mL
IL-1 $\beta$	15750	pg	63000	pg/mL
IL-2	11325	pg	45300	pg/mL
IL-4	10962.5	pg	43850	pg/mL
IL-5	3437.5	pg	13750	pg/mL
IL-6	21000	pg	84000	pg/mL
KC/GRO	2437.5	pg	9750	pg/mL
IL-10	27750	pg	111000	pg/mL
IL-12p70	109000	pg	436000	pg/mL
TNF- $\alpha$	7537.5	pg	30150	pg/mL
GM-CSF	1312.5	pg	5250	pg/mL
IL-13	24125	pg	96500	pg/mL
EPO	15000	pg	60000	pg/mL
VEGF-A	12450	pg	49800	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 is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Michelle Likiam	Quality		26 FEB 2025