



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 5

**Catalog Number:** C0065-2

**Lot Number:** A00U0273

**Expiration Date:** 30 JUN 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- $\gamma$	3988	pg	15950	pg/mL
IL-1 $\beta$	20000	pg	80000	pg/mL
IL-2	15500	pg	62000	pg/mL
IL-4	13000	pg	52000	pg/mL
IL-5	3988	pg	15950	pg/mL
IL-6	26125	pg	104500	pg/mL
KC-GRO	2713	pg	10850	pg/mL
IL-10	33000	pg	132000	pg/mL
IL12p70	123625	pg	494500	pg/mL
TNF- $\alpha$	8475	pg	33900	pg/mL
GM-CSF	1488	pg	5950	pg/mL
IL-13	23250	pg	93000	pg/mL
EPO	17000	pg	68000	pg/mL
VEGF-A	13250	pg	53000	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>	Karen Hamilla	Quality		15 JUL 2022