



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 5

Catalog Number: C0065-2

Lot Number: A00U0229

Expiration Date: 31 AUG 2026

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- γ	3950	pg	15800	pg/mL
IL-1 β	19250	pg	77000	pg/mL
IL-2	14875	pg	59500	pg/mL
IL-4	12750	pg	51000	pg/mL
IL-5	5213	pg	20850	pg/mL
IL-6	25625	pg	102500	pg/mL
KC-GRO	2663	pg	10650	pg/mL
IL-10	32500	pg	130000	pg/mL
IL12p70	126250	pg	505000	pg/mL
TNF- α	8875	pg	35500	pg/mL
GM-CSF	1450	pg	5800	pg/mL
IL-13	26625	pg	106500	pg/mL
EPO	16750	pg	67000	pg/mL
VEGF-A	13500	pg	54000	pg/mL

Reconstitution Volume: 250 μ 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
Review/Approval	Karen Hamilla	Quality		09 NOV 2021