



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 5

Catalog Number: C0065-2

Lot Number: A00U0244

Expiration Date: 30 APR 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- γ	3938	pg	15750	pg/mL
IL-1 β	17625	pg	70500	pg/mL
IL-2	15500	pg	62000	pg/mL
IL-4	13125	pg	52500	pg/mL
IL-5	5175	pg	20700	pg/mL
IL-6	25625	pg	102500	pg/mL
KC-GRO	2713	pg	10850	pg/mL
IL-10	31875	pg	127500	pg/mL
IL12p70	123000	pg	492000	pg/mL
TNF- α	8913	pg	35650	pg/mL
GM-CSF	1475	pg	5900	pg/mL
IL-13	22750	pg	91000	pg/mL
EPO	17125	pg	68500	pg/mL
VEGF-A	13125	pg	52500	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		16 JUN 2022