



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 5

Catalog Number: C0065-2

Lot Number: A00U0217

Expiration Date: 31 JUL 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	3988	pg	15950	pg/mL
IL-1 β	19750	pg	79000	pg/mL
IL-2	15250	pg	61000	pg/mL
IL-4	13500	pg	54000	pg/mL
IL-5	5238	pg	20950	pg/mL
IL-6	26500	pg	106000	pg/mL
KC-GRO	2825	pg	11300	pg/mL
IL-10	32625	pg	130500	pg/mL
IL12p70	126250	pg	505000	pg/mL
TNF- α	9213	pg	36850	pg/mL
GM-CSF	1463	pg	5850	pg/mL
IL-13	28250	pg	113000	pg/mL
EPO	17625	pg	70500	pg/mL
VEGF-A	13875	pg	55500	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 SEP 2021