



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 5

Catalog Number: C0065-2

Lot Number: A00U0199

Expiration Date: 31 JAN 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	4675	pg	18700	pg/mL
IL-1 β	18875	pg	75500	pg/mL
IL-2	14750	pg	59000	pg/mL
IL-4	13750	pg	55000	pg/mL
IL-5	3925	pg	15700	pg/mL
IL-6	24750	pg	99000	pg/mL
KC-GRO	2675	pg	10700	pg/mL
IL-10	32375	pg	129500	pg/mL
IL12p70	119250	pg	477000	pg/mL
TNF- α	9075	pg	36300	pg/mL
GM-CSF	1400	pg	5600	pg/mL
IL-13	28875	pg	115500	pg/mL
EPO	17625	pg	70500	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		12 MAY 2021