



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 5

Catalog Number: C0065-2

Lot Number: A00U0159

Expiration Date: 31 MAY 2025

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	4438	pg	17750	pg/mL
IL-1 β	17625	pg	70500	pg/mL
IL-2	13875	pg	55500	pg/mL
IL-4	13750	pg	55000	pg/mL
IL-5	3875	pg	15500	pg/mL
IL-6	22000	pg	88000	pg/mL
KC/GRO	2588	pg	10350	pg/mL
IL-10	29500	pg	118000	pg/mL
IL-12p70	121750	pg	487000	pg/mL
TNF- α	9200	pg	36800	pg/mL
GM-CSF	1338	pg	5350	pg/mL
IL-13	31500	pg	126000	pg/mL
EPO	17125	pg	68500	pg/mL
VEGF-A	13000	pg	52000	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		07 JUL 2020