



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 5
Catalog Number: C0065-2
Lot Number: A00U0442
Expiration Date: 31 MAY 2030
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	3300	pg	13200	pg/mL
IL-1 β	16625	pg	66500	pg/mL
IL-2	12175	pg	48700	pg/mL
IL-4	12750	pg	51000	pg/mL
IL-5	3787.5	pg	15150	pg/mL
IL-6	21750	pg	87000	pg/mL
KC/GRO	2625	pg	10500	pg/mL
IL-10	28375	pg	113500	pg/mL
IL-12p70	130000	pg	520000	pg/mL
TNF- α	8850	pg	35400	pg/mL
GM-CSF	1275	pg	5100	pg/mL
IL-13	23250	pg	93000	pg/mL
EPO	15125	pg	60500	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		24 JUL 2025