



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 5
Catalog Number: C0065-2
Lot Number: A00U0417
Expiration Date: 30 APR 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	3125	pg	12500	pg/mL
IL-1 β	15750	pg	63000	pg/mL
IL-2	11475	pg	45900	pg/mL
IL-4	10775	pg	43100	pg/mL
IL-5	3437.5	pg	13750	pg/mL
IL-6	20750	pg	83000	pg/mL
KC/GRO	2400	pg	9600	pg/mL
IL-10	26250	pg	105000	pg/mL
IL-12p70	103875	pg	415500	pg/mL
TNF- α	8587.5	pg	34350	pg/mL
GM-CSF	1287.5	pg	5150	pg/mL
IL-13	26375	pg	105500	pg/mL
EPO	15875	pg	63500	pg/mL
VEGF-A	12175	pg	48700	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		23 MAY 2025