



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 5
Catalog Number: C0065-2
Lot Number: A00U0377
Expiration Date: 31 MAY 2029
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	4062.5	pg	16250	pg/mL
IL-1 β	19875	pg	79500	pg/mL
IL-2	13375	pg	53500	pg/mL
IL-4	12500	pg	50000	pg/mL
IL-5	3762.5	pg	15050	pg/mL
IL-6	22625	pg	90500	pg/mL
KC/GRO	2712.5	pg	10850	pg/mL
IL-10	31750	pg	127000	pg/mL
IL-12p70	130000	pg	520000	pg/mL
TNF- α	8925	pg	35700	pg/mL
GM-CSF	1400	pg	5600	pg/mL
IL-13	27750	pg	111000	pg/mL
EPO	16875	pg	67500	pg/mL
VEGF-A	12750	pg	51000	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		08 JUL 2024