



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 5
Catalog Number: C0065-2
Lot Number: A00U0349
Expiration Date: 31 OCT 2028
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	4025	pg	16100	pg/mL
IL-1 β	18750	pg	75000	pg/mL
IL-2	12750	pg	51000	pg/mL
IL-4	12013	pg	48050	pg/mL
IL-5	3800	pg	15200	pg/mL
IL-6	20875	pg	83500	pg/mL
KC/GRO	3038	pg	12150	pg/mL
IL-10	30250	pg	121000	pg/mL
IL-12p70	12750	pg	510000	pg/mL
TNF- α	8775	pg	35100	pg/mL
GM-CSF	1350	pg	5400	pg/mL
IL-13	26875	pg	107500	pg/mL
EPO	16250	pg	65000	pg/mL
VEGF-A	12125	pg	48500	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		10 APR 2024