



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 5

Catalog Number: C0065-2

Lot Number: A00U0171

Expiration Date: 30 NOV 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	4563	pg	18250	pg/mL
IL-1 β	14500	pg	58000	pg/mL
IL-2	14375	pg	57500	pg/mL
IL-4	13750	pg	55000	pg/mL
IL-5	3950	pg	15800	pg/mL
IL-6	23375	pg	93500	pg/mL
KC/GRO	2700	pg	10800	pg/mL
IL-10	31250	pg	125000	pg/mL
IL-12p70	121125	pg	484500	pg/mL
TNF- α	9188	pg	36750	pg/mL
GM-CSF	1388	pg	5550	pg/mL
IL-13	25000	pg	100000	pg/mL
EPO	17750	pg	71000	pg/mL
VEGF-A	12875	pg	51500	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 JAN 2021