



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 5

Catalog Number: C0065-2

Lot Number: A00U0135

Expiration Date: 30 NOV 2024

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	4263	pg	17050	pg/mL
IL-1 β	13125	pg	52500	pg/mL
IL-2	12875	pg	51500	pg/mL
IL-4	13000	pg	52000	pg/mL
IL-5	3413	pg	13650	pg/mL
IL-6	21875	pg	87500	pg/mL
KC/GRO	2713	pg	10850	pg/mL
IL-10	28625	pg	114500	pg/mL
IL-12p70	119250	pg	477000	pg/mL
TNF- α	8413	pg	33650	pg/mL
GM-CSF	1400	pg	5600	pg/mL
IL-13	25125	pg	100500	pg/mL
EPO	16625	pg	66500	pg/mL
VEGF-A	13625	pg	54500	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		07 JAN 2019