



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 2
Catalog Number: C0061-2
Lot Number: A00U0437
Expiration Date: 31 MAY 2030
Storage Temperature: 2-8°C
Intended Assay Species: Human, NHP


Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	7200	pg	28800	pg/mL
Eotaxin-3	28125	pg	112500	pg/mL
IP-10	14000	pg	56000	pg/mL
MCP-1	9075	pg	36300	pg/mL
MCP-4	5300	pg	21200	pg/mL
MDC	30250	pg	121000	pg/mL
MIP-1 α	9075	pg	36300	pg/mL
MIP-1 β	2800	pg	11200	pg/mL
TARC/CCL17	4275	pg	17100	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		03 JUL 2025