



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 2
Catalog Number: C0061-2
Lot Number: A00U0468
Expiration Date: 28 FEB 2031
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	6825	pg	27300	pg/mL
Eotaxin-3	20700	pg	82800	pg/mL
IP-10	13575	pg	54300	pg/mL
MCP-1	6825	pg	27300	pg/mL
MCP-4	5150	pg	20600	pg/mL
MDC	24675	pg	98700	pg/mL
MIP-1 α	7000	pg	28000	pg/mL
MIP-1 β	2225	pg	8900	pg/mL
TARC/CCL17	3400	pg	13600	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		13 APR 2026