



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 2
Catalog Number: C0061-2
Lot Number: A00U0451
Expiration Date: 31 OCT 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	6900	pg	27600	pg/mL
Eotaxin-3	23425	pg	93700	pg/mL
IP-10	11100	pg	44400	pg/mL
MCP-1	7675	pg	30700	pg/mL
MCP-4	4850	pg	19400	pg/mL
MDC	24450	pg	97800	pg/mL
MIP-1 α	7400	pg	29600	pg/mL
MIP-1 β	2230	pg	8920	pg/mL
TARC/CCL17	3600	pg	14400	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		20 NOV 2025