



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 2
Catalog Number: C0061-2
Lot Number: A00U0484
Expiration Date: 30 JUN 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	7275	pg	29100	pg/mL
Eotaxin-3	23775	pg	95100	pg/mL
IP-10	12250	pg	49000	pg/mL
MCP-1	7650	pg	30600	pg/mL
MCP-4	5175	pg	20700	pg/mL
MDC	27750	pg	111000	pg/mL
MIP-1 α	7675	pg	30700	pg/mL
MIP-1 β	2357.5	pg	9430	pg/mL
TARC/CCL17	3550	pg	14200	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		24 JUL 2026