



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 2  
**Catalog Number:** C0061-2  
**Lot Number:** A00U0403  
**Expiration Date:** 31 MAR 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	6437.5	pg	25750	pg/mL
Eotaxin-3	23750	pg	95000	pg/mL
IP-10	12275	pg	49100	pg/mL
MCP-1	7875	pg	31500	pg/mL
MCP-4	5075	pg	20300	pg/mL
MDC	25625	pg	102500	pg/mL
MIP-1 $\alpha$	7850	pg	31400	pg/mL
MIP-1 $\beta$	2387.5	pg	9550	pg/mL
TARC/CCL17	3712.5	pg	14850	pg/mL

**Reconstitution Volume:** 250  $\mu$ 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
<b>Review/Approval</b>	Karen Hamila	Quality		14 APR 2025