



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 2  
**Catalog Number:** C0061-2  
**Lot Number:** A00U0380  
**Expiration Date:** 31 JUL 2029  
**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	5662.5	pg	22650	pg/mL
Eotaxin-3	23625	pg	94500	pg/mL
IP-10	14750	pg	59000	pg/mL
MCP-1	7625	pg	30500	pg/mL
MCP-4	4925	pg	19700	pg/mL
MDC	24750	pg	99000	pg/mL
MIP-1 $\alpha$	7575	pg	30300	pg/mL
MIP-1 $\beta$	2137.5	pg	8550	pg/mL
TARC/CCL17	3337.5	pg	13350	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 Hamilla	Quality		21 AUG 2024