



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 2

**Catalog Number:** C0061-2

**Lot Number:** A00U0214

**Expiration Date:** 30 JUN 2026

**Storage Temperature:** 2-8°C

**Intended Assay Species:** Human

### Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin	7763	pg	31050	pg/mL
Eotaxin-3	25500	pg	102000	pg/mL
IP-10	12625	pg	50500	pg/mL
MCP-1	7863	pg	31450	pg/mL
MCP-4	4838	pg	19350	pg/mL
MDC	22500	pg	90000	pg/mL
MIP-1 $\alpha$	7175	pg	28700	pg/mL
MIP-1 $\beta$	2300	pg	9200	pg/mL
TARC	3763	pg	15050	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>	Carissa B. Hubbard	Quality		19 AUG 2021