



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 2  
**Catalog Number:** C0061-2  
**Lot Number:** A00U0320  
**Expiration Date:** 31 MAY 2028  
**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	7562.5	pg	30250	pg/mL
Eotaxin-3	26750	pg	107000	pg/mL
IP-10	13250	pg	53000	pg/mL
MCP-1	7075	pg	28300	pg/mL
MCP-4	4600	pg	18400	pg/mL
MDC	27250	pg	109000	pg/mL
MIP-1 $\alpha$	6925	pg	27700	pg/mL
MIP-1 $\beta$	2775	pg	11100	pg/mL
TARC/CCL17	3412.5	pg	13650	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 was manufactured and tested according to MSD documents. This lot meets MSD's specifications. 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		09 AUG 2023