



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 2  
**Catalog Number:** C0061-2  
**Lot Number:** A00U0345  
**Expiration Date:** 30 SEP 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	5262.5	pg	21050	pg/mL
Eotaxin-3	21875	pg	87500	pg/mL
IP-10	10787.5	pg	43150	pg/mL
MCP-1	6525	pg	26100	pg/mL
MCP-4	4275	pg	17100	pg/mL
MDC	22625	pg	90500	pg/mL
MIP-1 $\alpha$	6525	pg	26100	pg/mL
MIP-1 $\beta$	2062.5	pg	8250	pg/mL
TARC/CCL17	2962.5	pg	11850	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		20 NOV 2023