



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 10  
**Catalog Number:** C0091-2  
**Lot Number:** A00U0365  
**Expiration Date:** 30 APR 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-2	6075	pg	24300	pg/mL
GRO- $\alpha$	3425	pg	13700	pg/mL
I-309	4175	pg	16700	pg/mL
MCP-2	2262.5	pg	9050	pg/mL
MCP-3	6550	pg	26200	pg/mL
M-CSF	2287.5	pg	9150	pg/mL
MIF	26625	pg	106500	pg/mL
MIP-5	40875	pg	163500	pg/mL
TRAIL	11437.5	pg	45750	pg/mL
YKL-40	6887.5	pg	27550	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		11 JUN 2024