



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 10

**Catalog Number:** C0091-2

**Lot Number:** A00U0200

**Expiration Date:** 31 JAN 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-2	7250	pg	29000	pg/mL
GRO- $\alpha$	3125	pg	12500	pg/mL
I-309	3088	pg	12350	pg/mL
MCP-2	2263	pg	9050	pg/mL
MCP-3	5213	pg	20850	pg/mL
M-CSF	2200	pg	8800	pg/mL
MIF	28500	pg	114000	pg/mL
MIP-5	33750	pg	135000	pg/mL
TRAIL	11800	pg	47200	pg/mL
YKL-40	6450	pg	25800	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		19 MAR 2021