



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 10

**Catalog Number:** C0091-2

**Lot Number:** A00U0269

**Expiration Date:** 30 APR 2027

**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	6913	pg	27650	pg/mL
GRO- $\alpha$	2700	pg	10800	pg/mL
I-309	2663	pg	10650	pg/mL
MCP-2	2075	pg	8300	pg/mL
MCP-3	4313	pg	17250	pg/mL
M-CSF	2200	pg	8800	pg/mL
MIF	26125	pg	104500	pg/mL
MIP-5	24625	pg	98500	pg/mL
TRAIL	10613	pg	42450	pg/mL
YKL-40	5675	pg	22700	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		05 JUL 2022