



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 10
Catalog Number: C0091-2
Lot Number: A00U0316
Expiration Date: 31 MAR 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-2	6800	pg	27200	pg/mL
GRO- α	3125	pg	12500	pg/mL
I-309	3950	pg	15800	pg/mL
MCP-2	2500	pg	10000	pg/mL
MCP-3	6462.5	pg	25850	pg/mL
M-CSF	2200	pg	8800	pg/mL
MIF	27875	pg	111500	pg/mL
MIP-5	47250	pg	189000	pg/mL
TRAIL	11500	pg	46000	pg/mL
YKL-40	5862.5	pg	23450	pg/mL

Reconstitution Volume: 250 μ 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
Review/Approval	Karen Hamilla	Quality		24 APR 2023