



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 10
Catalog Number: C0091-2
Lot Number: A00U0311
Expiration Date: 31 DEC 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	5137.5	pg	20550	pg/mL
GRO- α	2537.5	pg	10150	pg/mL
I-309	2612.5	pg	10450	pg/mL
MCP-2	2000	pg	8000	pg/mL
MCP-3	4662.5	pg	18650	pg/mL
M-CSF	2087.5	pg	8350	pg/mL
MIF	25875	pg	103500	pg/mL
MIP-5	24875	pg	99500	pg/mL
TRAIL	10312.5	pg	41250	pg/mL
YKL-40	5200	pg	20800	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		28 FEB 2023