



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 12
Catalog Number: C0092-2
Lot Number: A00U0341
Expiration Date: 31 AUG 2028
Storage Temperature: 2-8°C
Intended Assay Species: Mouse

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-9	9650	pg	38600	pg/mL
IL-17A/F	13875	pg	55500	pg/mL
IP-10	4437.5	pg	17750	pg/mL
MCP-1	1262.5	pg	5050	pg/mL
MIP-1 α	2537.5	pg	10150	pg/mL
MIP-1 β	36000	pg	144000	pg/mL
MIP-2	2912.5	pg	11650	pg/mL
MIP-3 α	3300	pg	13200	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		27 SEP 2023