



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 12
Catalog Number: C0092-2
Lot Number: A00U0297
Expiration Date: 31 DEC 2027
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	9412.5	pg	37650	pg/mL
IL-17A/F	14125	pg	56500	pg/mL
IP-10	6025	pg	24100	pg/mL
MCP-1	1687.5	pg	6750	pg/mL
MIP-1 α	2587.5	pg	10350	pg/mL
MIP-1 β	34625	pg	138500	pg/mL
MIP-2	3050	pg	12200	pg/mL
MIP-3 α	2937.5	pg	11750	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		31 JAN 2023