



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 16
Catalog Number: C0295-2
Lot Number: A00U0331
Expiration Date: 31 DEC 2025
Storage Temperature: $\leq -70^{\circ}\text{C}$
Intended Assay Species: Mouse

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Calibrator	Unit of Measure
BAFF	4387.5	pg	17550	pg/mL
6CKine/CCL21	1737.5	pg	6950	pg/mL
CD40/TNFRSF5	9137.5	pg	36550	pg/mL
BCA-1/BLC	22625	pg	90500	pg/mL
IFN- β	6875	pg	27500	pg/mL
MDC	4600	pg	18400	pg/mL
MCP-5/CCL12	1152.5	pg	4610	pg/mL

Volume: 250 μL

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 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		25 JUL 2023