



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 16  
**Catalog Number:** C0295-2  
**Lot Number:** A00U0390  
**Expiration Date:** 28 FEB 2027  
**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	4112.5	pg	16450	pg/mL
6CKine/CCL21	2000	pg	8000	pg/mL
CD40/TNFRSF5	9012.5	pg	36050	pg/mL
BCA-1/BLC	23125	pg	92500	pg/mL
IFN- $\beta$	6712.5	pg	26850	pg/mL
MDC	4975	pg	19900	pg/mL
MCP-5/CCL12	1136.25	pg	4545	pg/mL

**Volume:** 250  $\mu\text{L}$

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 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
<b>Review/Approval</b>	Karen Hamilla	Quality		10 OCT 2024