



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 16  
**Catalog Number:** C0295-2  
**Lot Number:** A00U0433  
**Expiration Date:** 30 NOV 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	4000	pg	16000	pg/mL
6CKine/CCL21	1975	pg	7900	pg/mL
CD40/TNFRSF5	8925	pg	35700	pg/mL
BCA-1/BLC	24250	pg	97000	pg/mL
IFN- $\beta$	7237.5	pg	28950	pg/mL
MDC	5462.5	pg	21850	pg/mL
MCP-5/CCL12	1130	pg	4520	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		02 JUL 2025