



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 16  
**Catalog Number:** C0295-2  
**Lot Number:** A00U0298  
**Expiration Date:** 30 JUN 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	3800	pg	15200	pg/mL
6CKine/CCL21	1800	pg	7200	pg/mL
CD40/TNFRSF5	8137.5	pg	32550	pg/mL
BCA-1/BLC	22375	pg	89500	pg/mL
IFN- $\beta$	6450	pg	25800	pg/mL
MDC	5450	pg	21800	pg/mL
MCP-5/CCL12	1061.25	pg	4245	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		01 FEB 2023