



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 17  
**Catalog Number:** C0296-2  
**Lot Number:** A00U0439  
**Expiration Date:** 30 APR 2030  
**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
MMP-9 (total)	85250	pg	341000	pg/mL
TNF-RI	2575	pg	10300	pg/mL
TARC/CCL17	1387.5	pg	5550	pg/mL
NGAL/LCN2	54000	pg	216000	pg/mL
SDF-1 $\alpha$	45500	pg	182000	pg/mL
RANTES	2475	pg	9900	pg/mL
Eotaxin	15750	pg	63000	pg/mL

**Reconstitution Volume:** 250  $\mu$ 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 is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		18 JUN 2025