



CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 17  
**Catalog Number:** C0296-2  
**Lot Number:** A00U0400  
**Expiration Date:** 30 NOV 2029  
**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)	77500	pg	310000	pg/mL
TNF-RI	2875	pg	11500	pg/mL
TARC/CCL17	1625	pg	6500	pg/mL
NGAL/LCN2	52750	pg	211000	pg/mL
SDF-1 $\alpha$	67000	pg	268000	pg/mL
RANTES	2325	pg	9300	pg/mL
Eotaxin	16750	pg	67000	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		17 JAN 2025