



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 17

Catalog Number: C0296-2

Lot Number: A00U0223

Expiration Date: 30 JUN 2026

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)	103750	pg	415000	pg/mL
TNF-RI	2788	pg	11150	pg/mL
TARC	1538	pg	6150	pg/mL
NGAL/LCN2	66875	pg	267500	pg/mL
SDF-1 α	61125	pg	244500	pg/mL
RANTES	2300	pg	9200	pg/mL
Eotaxin	13500	pg	54000	pg/mL

Reconstitution Volume: 250 μ 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
Review/Approval	Karen Hamilla	Quality		10 SEP 2021