



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 17
Catalog Number: C0296-2
Lot Number: A00U0470
Expiration Date: 30 JUN 2031
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)	87750	pg	351000	pg/mL
TNF-RI	2525	pg	10100	pg/mL
TARC/CCL17	1395	pg	5580	pg/mL
NGAL/LCN2	65500	pg	262000	pg/mL
SDF-1 α	57000	pg	228000	pg/mL
RANTES	2550	pg	10200	pg/mL
Eotaxin	16950	pg	67800	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	Michelle Likiam	Quality		04 AUG 2026