



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 17
Catalog Number: C0296-2
Lot Number: A00U0355
Expiration Date: 31 MAR 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)	82375	pg	329500	pg/mL
TNF-RI	2850	pg	11400	pg/mL
TARC/CCL17	1387.5	pg	5550	pg/mL
NGAL/LCN2	54125	pg	216500	pg/mL
SDF-1 α	65625	pg	262500	pg/mL
RANTES	2387.5	pg	9550	pg/mL
Eotaxin	15625	pg	62500	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		11 APR 2024