



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 17

Catalog Number: C0296-2

Lot Number: A00U0105

Expiration Date: 30 APR 2024

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)	109 500	pg	438 000	pg/mL
TNF-RI	2 588	pg	10 350	pg/mL
TARC	1 450	pg	5 800	pg/mL
NGAL/LCN2	68 875	pg	275 500	pg/mL
SDF-1 α	60 500	pg	242 000	pg/mL
RANTES	2 813	pg	11 250	pg/mL
Eotaxin	15 250	pg	61 000	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	Dawn Stollar	Quality		30 MAY 2019