



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 17

**Catalog Number:** C0296-2

**Lot Number:** A00U0114

**Expiration Date:** 30 JUN 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)	106 375	pg	425 500	pg/mL
TNF-RI	2 788	pg	11 150	pg/mL
TARC	1 575	pg	6 300	pg/mL
NGAL/LCN2	67 375	pg	269 500	pg/mL
SDF-1 $\alpha$	65 375	pg	261 500	pg/mL
RANTES	2 375	pg	9 500	pg/mL
Eotaxin	14 500	pg	58 000	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		07 AUG 2019