



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 29
Catalog Number: C0412-2
Lot Number: A00U0412
Expiration Date: 31 JAN 2030
Storage Temperature: 2-8°C
Intended Assay Species: Human

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
ICOS-L/B7-H2	27250	pg	109000	pg/mL
MMP-2	109500	pg	438000	pg/mL
proMMP-9	35500	pg	142000	pg/mL
P-Selectin	108000	pg	432000	pg/mL
RANTES	3975	pg	15900	pg/mL
S100A12	3650	pg	14600	pg/mL
TNF-RI	3025	pg	12100	pg/mL
TNF-RII	17475	pg	69900	pg/mL
MMP-9 (total)	35500	pg	142000	pg/mL

Reconstitution Volume: 250 µL

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		04 MAR 2025