



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 29  
**Catalog Number:** C0412-2  
**Lot Number:** A00U0362  
**Expiration Date:** 28 FEB 2029  
**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	41500	pg	166000	pg/mL
MMP-2	108750	pg	435000	pg/mL
proMMP-9	34500	pg	138000	pg/mL
MMP-9 (total)	34500	pg	138000	pg/mL
P-Selectin	187750	pg	751000	pg/mL
RANTES	3675	pg	14700	pg/mL
S100A12	3725	pg	14900	pg/mL
TNF-RI	4500	pg	18000	pg/mL
TNF-RII	18975	pg	75900	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
<b>Review/Approval</b>	Karen Hamilla	Quality		17 APR 2024