



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 29  
**Catalog Number:** C0412-2  
**Lot Number:** A00U0266  
**Expiration Date:** 30 APR 2027  
**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	36750	pg	147000	pg/mL
MMP-2	97000	pg	388000	pg/mL
proMMP-9	60750	pg	243000	pg/mL
MMP-9 (total)	60750	pg	243000	pg/mL
P-Selectin	175000	pg	700000	pg/mL
RANTES	4500	pg	18000	pg/mL
S100A12	3375	pg	13500	pg/mL
TNF-RI	4125	pg	16500	pg/mL
TNF-RII	18025	pg	72100	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 was manufactured and tested according to MSD documents. This lot meets MSD's specifications. 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		08 SEP 2023