



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 28
Catalog Number: C0411-2
Lot Number: A00U0455
Expiration Date: 31 AUG 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	38500	pg	154000	pg/mL
LIGHT/TNFSF14	11100	pg	44400	pg/mL
MMP-1	23075	pg	92300	pg/mL
MMP-7	20675	pg	82700	pg/mL
Nectin-4	24375	pg	97500	pg/mL
Perforin	46250	pg	185000	pg/mL
VEGFR-1/Flt-1	36000	pg	144000	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		13 OCT 2025