



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 28
Catalog Number: C0411-2
Lot Number: A00U0438
Expiration Date: 31 JUL 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	35750	pg	143000	pg/mL
LIGHT/TNFSF14	11350	pg	45400	pg/mL
MMP-1	23150	pg	92600	pg/mL
MMP-7	20575	pg	82300	pg/mL
Nectin-4	24525	pg	98100	pg/mL
Perforin	49000	pg	196000	pg/mL
VEGFR-1/Flt-1	36500	pg	146000	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		02 SEP 2025