



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 28
Catalog Number: C0411-2
Lot Number: A00U0482
Expiration Date: 31 MAY 2031
Storage Temperature: 2-8°C
Intended Assay Species: Human

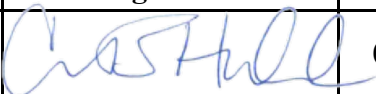
Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
ICOS	33500	pg	134000	pg/mL
LIGHT/TNFSF14	12025	pg	48100	pg/mL
MMP-1	24125	pg	96500	pg/mL
MMP-7	18150	pg	72600	pg/mL
Nectin-4	24525	pg	98100	pg/mL
Perforin	47250	pg	189000	pg/mL
VEGFR-1/Flt-1	32250	pg	129000	pg/mL

Reconstitution Volume:

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	Carissa B. Hubbard	Quality		09 JUL 2026