



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 28

Catalog Number: C0411-2

Lot Number: A00U0264

Expiration Date: 31 MAR 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	36500	pg	146000	pg/mL
LIGHT/TNFSF14	12275	pg	49100	pg/mL
MMP-1	23825	pg	95300	pg/mL
MMP-7	20675	pg	82700	pg/mL
Nectin-4	23900	pg	95600	pg/mL
Perforin	47000	pg	188000	pg/mL
VEGFR-1/Flt-1	30250	pg	121000	pg/mL

Reconstitution Volume: 250 μ L

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		03 JUN 2022