



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 28
Catalog Number: C0411-2
Lot Number: A00U0330
Expiration Date: 31 MAY 2028
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	36750	pg	147000	pg/mL
LIGHT/TNFSF14	12850	pg	51400	pg/mL
MMP-1	25500	pg	102000	pg/mL
MMP-7	21300	pg	85200	pg/mL
Nectin-4	24875	pg	99500	pg/mL
Perforin	47750	pg	191000	pg/mL
VEGFR-1/Flt-1	30250	pg	121000	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
Review/Approval	Karen Hamilla	Quality		15 AUG 2023