



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 28  
**Catalog Number:** C0411-2  
**Lot Number:** A00U0370  
**Expiration Date:** 30 JUN 2029  
**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	40250	pg	161000	pg/mL
LIGHT/TNFSF14	15575	pg	62300	pg/mL
MMP-1	26750	pg	107000	pg/mL
MMP-7	22100	pg	88400	pg/mL
Nectin-4	26000	pg	104000	pg/mL
Perforin	49750	pg	199000	pg/mL
VEGFR-1/Flt-1	42500	pg	170000	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
<b>Review/Approval</b>	Karen Hamilla	Quality		24 JUL 2024