



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 30
Catalog Number: C0717-2
Lot Number: A00U0461
Expiration Date: 31 DEC 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
E-Cadherin	53750	pg	215000	pg/mL
Fas (soluble)	4025	pg	16100	pg/mL
proMMP-8	18250	pg	73000	pg/mL
MMP-8 (total)	18250	pg	73000	pg/mL
PDGF-B	2500	pg	10000	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	Michelle Likiam	Quality		24 FEB 2026