



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 30
Catalog Number: C0717-2
Lot Number: A00U0346
Expiration Date: 31 OCT 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
E-Cadherin	56500	pg	226000	pg/mL
Fas (soluble)	4000	pg	16000	pg/mL
proMMP-8	18275	pg	73100	pg/mL
MMP-8 (total)	18275	pg	73100	pg/mL
PDGF-B	2625	pg	10500	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		02 FEB 2024