



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 11
Catalog Number: C0244-2
Lot Number: A00U0233
Expiration Date: 30 SEP 2026
Storage Temperature: 2-8°C
Intended Assay Species: Human, Mouse, NHP

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
TGF- β 1	49500	pg	198000	pg/mL
TGF- β 2	46500	pg	186000	pg/mL
TGF- β 3	44125	pg	176500	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		07 DEC 2021