



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 11  
**Catalog Number:** C0244-2  
**Lot Number:** A00U0207  
**Expiration Date:** 31 MAY 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- $\beta$ 1	47000	pg	188000	pg/mL
TGF- $\beta$ 2	48000	pg	192000	pg/mL
TGF- $\beta$ 3	57875	pg	231500	pg/mL

**Reconstitution Volume:** 250  $\mu$ 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
<b>Review/Approval</b>	Karen Hamilla	Quality		04 AUG 2021