



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 11

**Catalog Number:** C0244-2

**Lot Number:** A00U0034

**Expiration Date:** 31 DEC 2021

**Storage Condition:** 2-8°C

Analyte Description	Amount of Analyte in Calibrator (pg)	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL) (pg/mL)	Concentration of Standard 1 (after 5-fold Dilution) (pg/mL)
TGF-β1	70 750	283 000	56 600
TGF-β2	56 750	227 000	45 400
TGF-β3	57 250	229 000	45 800

**Statement:** The above calibrator lot was tested using MSD procedures and conforms to MSD's specifications. It is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	15 FEB 2017