



## CERTIFICATE OF ANALYSIS

**Product Description:** Calibrator 23  
**Catalog Number:** C0331-2  
**Lot Number:** A00U0449  
**Expiration Date:** 30 NOV 2027  
**Storage Temperature:**  $\leq -70^{\circ}\text{C}$   
**Intended Assay Species:** Human


**Concentration Assignment:**

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Calibrator	Unit of Measure
Granzyme A	8900	pg	35600	pg/mL
Granzyme B	1812.5	pg	7250	pg/mL
LAG3	89500	pg	358000	pg/mL

**Volume:** 250  $\mu\text{L}$

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 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
<b>Review/Approval</b>	Karen Hamilla	Quality		07 JUL 2025