



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 23
Catalog Number: C0331-2
Lot Number: A00U0456
Expiration Date: 31 MAR 2028
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	8875	pg	35500	pg/mL
Granzyme B	1780	pg	7120	pg/mL
LAG3	92750	pg	371000	pg/mL

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 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		14 OCT 2025