



CERTIFICATE OF ANALYSIS

Product Description: Calibrator 23
Catalog Number: C0331-2
Lot Number: A00U0319
Expiration Date: 30 SEP 2025
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	8525	pg	34100	pg/mL
Granzyme B	1855	pg	7420	pg/mL
LAG3	71250	pg	285000	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		17 APR 2023