



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 23  
**Catalog Number:** C0331-2  
**Lot Number:** A00U0208  
**Expiration Date:** 30 SEP 2023  
**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	9250	pg	37000	pg/mL
Granzyme B	1968	pg	7870	pg/mL
LAG3	109000	pg	436000	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 is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Carissa B. Hubbard	Quality		07 JUL 2021