



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 23

**Catalog Number:** C0331-2

**Lot Number:** A00U0251

**Expiration Date:** 31 JUL 2024

**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	8800	pg	35200	pg/mL
Granzyme B	1900	pg	7600	pg/mL
LAG3	94500	pg	378000	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
Review/Approval	Karen Hamilla	Quality		06 MAY 2022