



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Calibrator 23

**Catalog Number:** C0331-2

**Lot Number:** A00U0241

**Expiration Date:** 30 APR 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	9525	pg	38100	pg/mL
Granzyme B	2023	pg	8090	pg/mL
LAG3	102250	pg	409000	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>	Karen Hamilla	Quality		19 JAN 2022