



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 23

Catalog Number: C0331-2

Lot Number: A00U0188

Expiration Date: 31 JUL 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	8725	pg	34900	pg/mL
Granzyme B	1883	pg	7530	pg/mL
LAG3	101750	pg	407000	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 is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Carissa B. Hubbard	Quality		19 FEB 2021