



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Human IFN- $\gamma$  Calibrator

**Catalog Number:** C01AE-2

**Lot Number:** A0032472

**Expiration Date:** 30 JUN 2023

**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

### Concentration Assignments:

	Nominal Concentration		Measured Concentration	
	Stock Concentration ( $\mu\text{g/mL}$ )	Top of Curve Concentration ( $\text{pg/mL}$ )	Stock Concentration ( $\mu\text{g/mL}$ )	Top of Curve Concentration ( $\text{pg/mL}$ )
IFNg	1.00	10,000	1.02	10200

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		16 DEC 2020