



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Human IFN- γ Calibrator

Catalog Number: C01AE-2

Lot Number: A0032308

Expiration Date: 30SEP2019

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

Concentration Assignments:

	Nominal Concentration		Measured Concentration	
	Stock Concentration ($\mu\text{g/mL}$)	Top of Curve Concentration (pg/mL)	Stock Concentration ($\mu\text{g/mL}$)	Top of Curve Concentration (pg/mL)
IFN- γ	1.00	10,000	0.965	9650

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	16 MAY 2017