



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA - 07325

Product Description: Human IFN- γ Calibrator

Catalog Number: C01XX-2

Lot Number: A0032266

Expiration Date: 28Feb2019

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

Concentration Assignments:

	Nominal Concentration		Measured Concentration	
	Stock Concentration ($\mu\text{g}/\text{mL}$)	Top of Curve Concentration (pg/mL)	Stock Concentration ($\mu\text{g}/\text{mL}$)	Top of Curve Concentration (pg/mL)
IFN- γ	1.00	10,000	0.98	9,790

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	28 SEP 2016