



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA - 07555

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

**Catalog Number:** C01XX-2

**Lot Number:** A0032280

**Expiration Date:** 30 APR 2019

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

### Concentration Assignments:

	Nominal Concentration		Measured Concentration	
	Stock Concentration ( $\mu\text{g}/\text{mL}$ )	Top of Curve Concentration ( $\text{pg}/\text{mL}$ )	Stock Concentration ( $\mu\text{g}/\text{mL}$ )	Top of Curve Concentration ( $\text{pg}/\text{mL}$ )
IFN- $\gamma$	1.00	10,000	0.98	9,840

**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	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	01 DEC 2016