



## CERTIFICATE OF ANALYSIS


**Product Description:** Human IFN- $\gamma$  Calibrator  
**Catalog Number:** C01X9-2  
**Lot Number:** A00S0117  
**Expiration Date:** 30 NOV 2029  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\gamma$	33400	fg/mL

**Reconstitution Volume:** 1000  $\mu$ L

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		17 JAN 2025