



CERTIFICATE OF ANALYSIS

Product Description: Human IFN- γ Calibrator
Catalog Number: C01X9-2
Lot Number: A00S0158
Expiration Date: 31 MAY 2031
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	33500	fg/mL

Reconstitution Volume: 1000 μ 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
Review/Approval	Karen Hamilla	Quality		16 JUL 2026