



CERTIFICATE OF ANALYSIS

Product Description: Human IFN- α 2a Calibrator
Catalog Number: C01P3-2
Lot Number: A00S0109
Expiration Date: 31 AUG 2029
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- α 2a	50500000	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		06 NOV 2024