



## CERTIFICATE OF ANALYSIS

**Product Description:** Human IFN- $\alpha$ 2a Calibrator  
**Catalog Number:** C01P3-2  
**Lot Number:** A00S0146  
**Expiration Date:** 31 JAN 2031  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- $\alpha$ 2a	52000000	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
Review/Approval	Karen Hamilla	Quality		26 MAR 2026