



CERTIFICATE OF ANALYSIS

Product Description: Human IFN- ω Calibrator
Catalog Number: C01AXD-2
Lot Number: A00S0149
Expiration Date: 31 AUG 2028
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
IFN- ω	1158000	fg/mL

Dilution to Top of Curve: 20 X

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		17 APR 2026