



CERTIFICATE OF ANALYSIS

Product Description: Mouse IFN- α Calibrator
Catalog Number: C02W1-2
Lot Number: A00U0334
Expiration Date: 31 DEC 2026
Storage Temperature: 2-8°C
Intended Assay Species: Mouse

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- α	117375	pg	469500	pg/mL

Reconstitution Volume: 250 μ L μ L

To calculate the **Calibrator Standard 1 (top of curve) concentration**, please refer to the dilution-factor used in the **Prepare Calibrator Standards** section of the respective product insert. Divide the stock concentration of the reconstituted calibrator by this dilution factor to generate the working calibrator standard 1 concentration.

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		31 JUL 2023