



CERTIFICATE OF ANALYSIS

Product Description: Mouse IFN- α Calibrator
Catalog Number: C02W1-2
Lot Number: A00U0394
Expiration Date: 30 SEP 2028
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	123500	pg	494000	pg/mL

Reconstitution Volume: 250 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 is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		15 APR 2025