



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Mouse IFN- $\alpha$  Calibrator

**Catalog Number:** C02W1-2

**Lot Number:** A00U0174

**Expiration Date:** 30 SEP 2025

**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- $\alpha$	128750	pg	515000	pg/mL

**Reconstitution Volume:** 250  $\mu$ 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
<b>Review/Approval</b>	Karen Hamilla	Quality		17 NOV 2020