

**CERTIFICATE OF ANALYSIS**


Product Description: Human TNF- β Calibrator
Catalog Number: C01ADP-2
Lot Number: A00S0122
Expiration Date: 31 JAN 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
TNF- β	35800	fg/mL

Reconstitution Volume: 1000 μ 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		21 JUL 2025