

**CERTIFICATE OF ANALYSIS**


Product Description: Human IL-1 β Calibrator
Catalog Number: C01ADS-2
Lot Number: A00S0129
Expiration Date: 31 MAR 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-1 β	115000	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		15 MAY 2025