



CERTIFICATE OF ANALYSIS

Product Description: Human IL-6 Calibrator
Catalog Number: C012M-2
Lot Number: A00S0143
Expiration Date: 30 SEP 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IL-6	744000	fg/mL

Reconstitution Volume: 1000 µL

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		14 OCT 2025