

**CERTIFICATE OF ANALYSIS**


Product Description: Human Eotaxin-3 Calibrator
Catalog Number: C01Q3-2
Lot Number: A00S0132
Expiration Date: 31 MAY 2030
Storage Temperature: 2-8°C

Concentration Assignment:

Analyte	Stock Concentration of Reconstituted Calibrator	Unit of Measure
Eotaxin-3	655000	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		27 JUN 2025