

**CERTIFICATE OF ANALYSIS**


Product Description: Human IL-5 Calibrator
Catalog Number: C01J3-2
Lot Number: A00S0131
Expiration Date: 31 MAR 2030
Storage Temperature: 2-8°C

Concentration Assignment:

| Analyte | Stock Concentration of Reconstituted Calibrator | Unit of Measure |
|---------|---|-----------------|
| IL-5 | 964000 | 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 |  | 06 MAY 2025 |