



## CERTIFICATE OF ANALYSIS

**Product Description:** Human IL-1 $\beta$  Calibrator  
**Catalog Number:** C01ADS-2  
**Lot Number:** 590649  
**Expiration Date:** 30 JUN 2031  
**Storage Temperature:** 2-8°C

**Concentration Assignment:**

| Analyte      | Stock Concentration of Reconstituted Calibrator | Unit of Measure |
|--------------|-------------------------------------------------|-----------------|
| IL-1 $\beta$ | 107000                                          | fg/mL           |

**Reconstitution Volume:** 1000  $\mu$ 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  |  | 14 AUG 2026 |