




### CERTIFICATE OF CONFORMANCE

**Product Description:** Diluent 3  
**Catalog Number:** R51BA-Series; R50AP-Series  
**Lot Number:** M0030127  
**Expiration Date** 30 SEP 2026  
**Storage Temperature:** ≤ -10°C

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.

|                        | <b>Name</b>   | <b>Function</b> | <b>Signature</b>   | <b>Date</b> |
|------------------------|---------------|-----------------|--|-------------|
| <b>Review/Approval</b> | Karen Hamilla | Quality         |  | 23 APR 2024 |