



CERTIFICATE OF CONFORMANCE

Product Description: U-PLEX Linker 10
Catalog Number: E2235-Series
Lot Number: L0100025
Expiration Date: 31 AUG 2025
Storage Temperature: 2-8°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.

| | Name | Function | Signature | Date |
|------------------------|---------------|-----------------|----------------------|-------------|
| Review/Approval | Karen Hamilla | Quality | <i>Karen Hamilla</i> | 03 AUG 2023 |