




CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 3
Catalog Number: R51BA-Series; R50AP-Series
Lot Number: M0030107
Expiration Date: 30 SEP 2023
Storage Temperature: $\leq -10^{\circ}\text{C}$

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 |  | 30 SEP 2021 |