

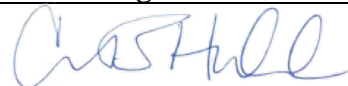


**CERTIFICATE OF CONFORMANCE**

Page 1 of 1

**Product Description:** Diluent 3  
**Catalog Number:** R50AP- Series; R51BA-Series  
**Lot Number:** M0030088  
**Expiration Date:** 30 SEP 2021  
**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> | Carissa B. Hubbard | Quality         |  | 06 APR 2021 |