




CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 43
Catalog Number: R50AG - Series
Lot Number: M0430033
Expiration Date: 31 JAN 2022
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.

| | Name | Function | Signature | Date |
|------------------------|---------------|-----------------|--|-------------|
| Review/Approval | Karen Hamilla | Quality |  | 06 SEP 2019 |