




### CERTIFICATE OF CONFORMANCE

Page 1 of 1

**Product Description:** Diluent 43  
**Catalog Number:** R50AG-Series  
**Lot Number:** M0430045  
**Expiration Date:** 30 NOV 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 | Xiomara Talavera | Quality  |  | 17 AUG 2021 |