



### CERTIFICATE OF CONFORMANCE

Page 1 of 1

**Product Description:** Diluent 7  
**Catalog Number:** R54BB-Series  
**Lot Number:** M0070066  
**Expiration Date:** 30 JUN 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.

c

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality	<i>Karen Hamilla</i>	22 FEB 2021