



### CERTIFICATE OF CONFORMANCE

Page 1 of 1

**Product Description:** Diluent 7  
**Catalog Number:** R54BB-Series  
**Lot Number:** M0070063  
**Expiration Date:** 31 OCT 2022  
**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	<i>Karen Hamilla</i>	08 JUN 2020