



## CERTIFICATE OF CONFORMANCE

Page 1 of 1

**Product Description:** Diluent 11  
**Catalog Number:** R55BA - Series  
**Lot Number:** M0110044  
**Expiration Date:** 30 JUN 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.

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Karen Hamilla	Quality	<i>Karen Hamilla</i>	21 JAN 2020