



## CERTIFICATE OF CONFORMANCE

Page 1 of 1

**Product Description:** Diluent 100


**Catalog Number:** R50AA-Series

**Lot Number:** Y0270284

**Expiration Date:** 31 MAR 2023

**Storage Temperature:** 2-8°C

This product was manufactured according to MSD documents. 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		09 NOV 2020