



## CERTIFICATE OF CONFORMANCE

Page 1 of 1

**Product Description:** Diluent 100


**Catalog Number:** R50AA-Series

**Lot Number:** Y0270283

**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.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		15 OCT 2020