



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 100


Catalog Number: R50AA-Series

Lot Number: Y0270299

Expiration Date: 31 JAN 2024

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		24 AUG 2021