



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 11
Catalog Number: R55BA-Series
Lot Number: M0110050
Expiration Date: 30 SEP 2023
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>	12 MAY 2021