



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 13
Catalog Number: R56BB - Series
Lot Number: M0130050
Expiration Date: 31 AUG 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.

c

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality	<i>Karen Hamilla</i>	12 APR 2021