



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 12

Catalog Number: R50JA-Series

Lot Number: M0120069

Expiration Date: 31 MAR 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>	09 NOV 2020