




CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 3
Catalog Number: R50AP-Series, R51BA-Series
Lot Number: M0030111
Expiration Date: 31 MAY 2024
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		10 JAN 2022