




CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 7
Catalog Number: R54BB-Series
Lot Number: M0070064
Expiration Date: 30 NOV 2022
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		29 JUN 2020