




CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 37
Catalog Number: R50AF - Series
Lot Number: M0370024
Expiration Date: 31 OCT 2021
Storage Temperature: $\leq -10^{\circ}\text{C}$

This product was manufactured according to MSD documents. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		25 JUL 2019