



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 37

Catalog Number: R50AF-Series

Lot Number: M0370032

Expiration Date: 30 SEP 2023

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	Xiomara Talavera	Quality	<i>X Talavera</i>	16 AUG 2021