



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 40
Catalog Number: R50AJ-Series
Lot Number: M0400016
Expiration Date: 30 NOV 2022
Storage Temperature: ≤ -10°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	Xiomara Talavera	Quality	<i>Talavera</i>	11 OCT 2021