



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 43

Catalog #/Size: R50AG - Series


Lot Number: M0430030

Expiration Date: 30 JUN 2021

Storage Condition: $\leq -10^{\circ}\text{C}$

Statement: This product was manufactured according to approved MSD documents.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance		20 MAR 2019