



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: Diluent 2

Catalog #/Size: R51BB - Series

Lot Number: M0020096

Expiration Date: 30 NOV 2019

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	<i>Dawn Stollar</i>	31 OCT 2017