



CERTIFICATE OF CONFORMANCE

Product Description: Diluent 101
Catalog Number: R51AD-Series
Lot Number: M1010052
Expiration Date: 31 MAY 2026
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	Michelle Likiam	Quality		02 JAN 2024