




CERTIFICATE OF CONFORMANCE

Product Description: Diluent 7
Catalog Number: R54BB-Series
Lot Number: M0070076
Expiration Date 30 NOV 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	Karen Hamilla	Quality		17 JUN 2024