



CERTIFICATE OF CONFORMANCE

Product Description: Diluent 58
Catalog Number: R50CA-Series
Lot Number: M0580010
Expiration Date 30 APR 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	<i>Karen Hamilla</i>	20 NOV 2023