



CERTIFICATE OF CONFORMANCE

Product Description: Diluent 58
Catalog Number: R50CA-Series
Lot Number: M0580006
Expiration Date: 30 APR 2025
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	Xiomara Talavera	Quality	<i>X Talavera</i>	21 FEB 2023