



CERTIFICATE OF CONFORMANCE

Product Description: Diluent 11
Catalog Number: R55BA-Series
Lot Number: M0110058
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>Talavera</i>	05 JAN 2023