



**CERTIFICATE OF CONFORMANCE**

**Product Description:** Diluent 12  
**Catalog Number:** R50JA-Series  
**Lot Number:** M0120072  
**Expiration Date:** 31 DEC 2023  
**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.

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Xiomara Talavera-Herrera	Quality	<i>Talavera</i>	08 JUL 2022