




**CERTIFICATE OF CONFORMANCE**

Page 1 of 1

**Product Description:** Diluent 3  
**Catalog Number:** R50AP - Series  
**Lot Number:** M0030094  
**Expiration Date:** 31 MAY 2022  
**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>	Carissa Hubbard	Quality		23 DEC 2019