



**CERTIFICATE OF CONFORMANCE**

Page 1 of 1

**Product Description:** U-PLEX® Linker 8


**Catalog Number:** E2233-Series

**Lot Number:** L0080014

**Expiration Date:** 31 DEC 2022

**Storage Temperature:** 2-8°C

This product was manufactured according to MSD documents. 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 B. Hubbard	Quality		22 DEC 2020