



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: U-PLEX Linker 8


Catalog Number: E2233-Series

Lot Number: L0080015

Expiration Date: 30 JUN 2023

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.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		26 MAY 2021