



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: U-PLEX® Linker 3

Catalog Number: E2228-Series

Lot Number: L0030015

Expiration Date: 31 JUL 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.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality	<i>Karen Hamilla</i>	13 APR 2020