



## RFE MEMO

Page 1 of 1

**Product Description(s):** U-PLEX® Linker 9

**Catalog Number(s) / Size(s):** E2234-Series

**Lot Number:** L0090004

Based on additional testing, the original date of expiration for this kit linker has been extended by 21 months. The expiration date of L0090004 has been extended to 29 Jul 2019.

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar	<i>Dawn Stollar</i>	02 AUG 2018

This memo supersedes all other statements regarding date of expiration for this product.



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-05980

**Product Description:** U-PLEX<sup>®</sup> Linker 9

**Catalog Number:** E2234-3

**Lot Number:** L0090004

**Expiration Date:** 30 SEP 2017

**Storage Condition:** 2-8°C

**Statement:** The above lot was tested using MSD procedures and conforms to MSD's specifications. It is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Monika Jain	Quality Assurance	<i>Monika Jain</i>	14 SEP 2015