



RFE MEMO


Page 1 of 1

Product Description(s): MSD TAG™ NHS Ester 150 nMoles

Catalog Number(s) / Size(s): R91BN-1/150 nmol

Lot Number: T0030004

Based on additional testing, the original date of expiration for this reagent has been extended by 12 months. The expiration date of T0030004 has been extended to 31 MAR 2020.

Function	Name	Signature	Date
Quality	Dawn Stollar		22 MAY 2019

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



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD TAG™ NHS Ester 150 nmol
Catalog #/Size: R91BN-1
Lot Number: T0030004
Expiration Date: 31 MAR 2019

Purity specification (by HPLC): > 95%
Measured Average Purity: 100%

Quantity: 100% to 120% of nominal
Measured Average Quantity: 113%

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	09 OCT 2017