



CERTIFICATE OF ANALYSIS

Product Description: MSD TAG NHS Ester 500 nMoles
Catalog Number: R91BN-2
Lot Number: T0040010
Expiration Date 28 FEB 2029

Functional Results:

	Specification	Measured Average
Purity (by HPLC):	≥ 95%	100%
Quantity	100% to 120% of nominal	102%

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		15 MAY 2024