




CERTIFICATE OF ANALYSIS

Product Description: MSD TAG NHS Ester 500 nMoles
Catalog Number: R91BN-2
Lot Number: T0040011
Expiration Date 30 NOV 2030

Functional Results:

	Specification	Measured Average
Purity (by HPLC)	$\geq 95\%$	99%
Quantity	100% to 120% of nominal	107%

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		14 NOV 2025