



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: U-PLEX[®] 2-Assay, 96-Well SECTOR[®] Plate

Catalog Number: N05227A-1

Lot Number: Z00U0081A


Expiration Date: 31JUL2021

Storage Condition: 2-8°C

Acceptance Data:

Metric	Specifications	All Spots
Average Intraplate CV	$\leq 10.0\%$	Pass
CV of Intraplate Averages	$\leq 13.0\%$	Pass
Signal Range at Binding Capacity	$\leq 30.0\%$	Pass

Statement: The above plate lot is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance		22 MAR 2019