



## CERTIFICATE OF CONFORMANCE

Page 1 of 1

**Product Description:** U-PLEX<sup>®</sup> Human PD-L1 (Epitope 1) Antibody Set

**Catalog Number:** B22Z7-Series

**Lot Number:** SET01675A

**Expiration Date:** 31 DEC 2024

**Storage Temperature:** 2-8°C

Antibody Set Components		
Description	Lot Number	Expiration Date
Biotin Hu PD-L1 (Epitope 1) Antibody	B00U1675A	31 DEC 2024
SULFO-TAG <sup>™</sup> Human PD-L1 (Epitope 1) Antibody	D00U1675A	31 DEC 2024

**Statement:** This product was manufactured according to approved MSD documents. The above antibody set lot is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Xiomara Talavera	Quality		06 JAN 2022