




### CERTIFICATE OF CONFORMANCE

**Product Description:** 96-Well 1-Spot High Bind SECTOR Plates  
**Catalog Number:** L15XB-Series  
**Lot Number:** 22Q0IA  
**Expiration Date:** 30 APR 2028  
**Storage Temperature:** Room Temperature

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.

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Karen Hamilla	Quality		21 MAY 2026