



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: MULTI-ARRAY® 96-Well Small Spot GAM Plate

Catalog Number: L45MA-1

Lot Number: Z0070146

Expiration Date: 30 JUN 2022

Storage Temperature: 2-8°C

This product was manufactured according to MSD documents. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Xiomara Talavera	Quality		29 OCT 2021