




CERTIFICATE OF CONFORMANCE

Product Description: Human ACE2 Protein
Catalog Number: D21ADG-Series
Lot Number: D0082224
Expiration Date: 31 OCT 2027
Storage Temperature: 2-8°C

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		30 MAY 2024