



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: cAMP Lysis Buffer

Catalog Number: R60AE-1

Lot Number: Y0450033

Expiration Date: 30 NOV 2020

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	Karen Hamilla	Quality	<i>Karen Hamilla</i>	12 NOV 2019