



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: MSD cAMP Lysis Buffer


Catalog Number: R60AE-1

Lot Number: Y0450037

Expiration Date: 31 JUL 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 | Karen Hamilla | Quality |  | 08 NOV 2021 |