



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: cAMP Standard

Catalog Number: 3-3048-020001

Lot Number: A0010481

Expiration Date: 31 MAY 2022

Storage Temperature: $\leq -10^{\circ}\text{C}$

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	<i>Karen Hamilla</i>	21 NOV 2019