




RFE MEMO

Page 1 of 1

Product Description(s): cAMP Standard
Catalog Number(s) / Size(s): 3-3048-020001
Lot Number: A0010481

Based on additional testing, the original date of expiration for this Calibrator has been extended by 17 months. The expiration date of A0010481 has been extended to 31 OCT 2023.

Function	Name	Signature	Date
Quality	Karen Hamilla		31 OCT 2022

This memo supersedes all other statements regarding date of expiration for this product.



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