



CERTIFICATE OF ANALYSIS

COA - 06425

Product Description: cAMP Standard
Catalog #/Size: 3-3048-020001 / 50 μ L
Lot Number: A0010402
Expiration Date: 30 APR 2019
Statement: This product was manufactured according to approved MSD documents.

Concentration/Specific Activity: 1mM
Storage buffer: N/A
Storage conditions: $\leq -10^{\circ}$ C

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Xiomara Talavera	Quality Control	<i>Talavera</i>	04 JAN 2017
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	04 JAN 2017