

AlphaScreen®

cAMP Detection Kit**Product number:** 6760635M**Lot Number:** 3400495**Product Format:** 6760635D: 1000 assay points

6760635M: 10 000 assay points

6760635R: 50 000 assay points

The number of assay points is based on an assay volume of 25 µL in 384-well assay plates using a final bead concentration of 20 µg/mL.

Manufacturing date: April 15, 2025**Document version:** 1**Kit Components**

Component	6760635D	6760635M	6760635R
Anti-cAMP Acceptor Beads at 5 mg/mL in 1X PBS, 0.05% Kathon, pH 7.2	1 x 100 µL (6760100B)	1 x 1 mL (6760101)	1 x 5 mL (6760101B)
Streptavidin Donor Beads at 5 mg/mL in 1X PBS, 0.05% Kathon, pH 7.2	1 x 100 µL (6760003B)	1 x 1 mL (6760004)	1 x 5 mL (6760004B)
Biotinylated-cAMP Tracer Solid	1 x 10 nmol (6760301)	1 x 10 nmol (6760301)	1 x 65 nmol (6760301B)
cAMP standard at 50 µM	1 x 1 mL (FP1283)	1 x 1 mL (FP1283)	1 x 1 mL (FP1283)
10x Buffer : AlphaLISA Immunoassay Buffer**	1 x 2 mL (6760040G)	2 x 10 mL (6760040G)	1 x 100 mL (6760040G)

** Extra buffer can be ordered separately (cat # AL000C: 10 mL, cat # AL000F: 100 mL).

1X PBS pH 7.4 is not included. It can be ordered from Gibco (cat# 10010-023)

Product Information

Antibody: The cAMP antibody is a rabbit monoclonal antibody highly specific for cyclic AMP. Its cross-reactivity with cGMP is < 0.005%.

Stability: This kit is stable for at least 12 months from the date of manufacture when stored in its original packaging and the recommended storage conditions.

Storage: Store undiluted at 4°C protected from light. Freeze-thaw is not recommended and cause the beads to form aggregates.

Recommended Use: AlphaScreen® Donor beads are light-sensitive. All Alpha assays using the Donor beads should be performed under subdued laboratory lighting (< 100 lux). Green filters (LEE 090 filters) can be applied to light fixtures.

Quality Control

Alpha maximum signal, minimum signal, LDL and IC50 are determined by a cAMP competition assay performed on an EnVision® detection instrument. We certify that these results meet our requirements.

Maximum Signal:	125538 counts
Minimum Signal:	137 counts
LDL:	0.002 nM
IC50:	9.51 nM

Recommended Assay Conditions

Note: This protocol provides a method to verify kit performance and is not representative of an assay. Sufficient biotinylated cAMP-Tracer and 10x buffer is provided to perform 3 competition curves in triplicate as described.

1x Buffer:	Add 500 µL 10x buffer to 4.45 mL Milli-Q® H2O (or equivalent).
Acceptor Beads:	Add 5 µL Anti-cAMP Acceptor beads to 495 µL 1x buffer.
Biotinylated cAMP tracer:	Dissolve biotinylated cAMP-Tracer (10 nmol) in 1 mL 1X PBS (final 10 µM) Further dilute 10X into 1 X PBS, pH 7.4 to obtain a 1 µM solution **Solids may not be visible deposited along the walls of the tube**
Donor beads:	Add 6.25 µL Streptavidin Donor beads to 579.75 µL 1x buffer.
cAMP Standard:	From the 50 µM cAMP standard, prepare a ½ log dilution series (5 µM to 50 pM) in 1x buffer. Include a buffer only control.

Titration Protocol

To a white opaque OptiPlate-384:

- 1) Add 10 µL of Anti-cAMP Acceptor beads.
- 2) Add 5 µL cAMP standard dilutions (from lowest to highest concentration).
- 3) Pre-mix the biotinylated cAMP-Tracer and Streptavidin Donor beads from previously diluted solutions by adding 39 µL of biotinylated cAMP-Tracer to the Streptavidin Donor beads dilution.

Incubate the pre-mix and the plate in the dark at room temperature for 30 minutes

- 4) Add 10 µL of premix biotinylated cAMP-Tracer /Streptavidin Donor beads.

Incubate in the dark at room temperature for 60 minutes and analyze on your Alpha capable detection reader.

Typical Product Data

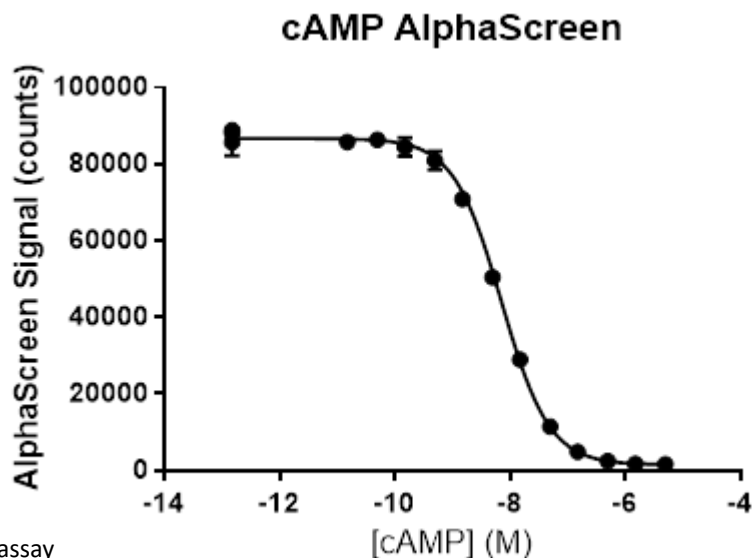


Figure 1: Competition assay
384-well cAMP competition curve (25 μ L final volume; Reader: Envision).

Note: Alpha signal will vary depending on instrument detection protocol, incubation temperature and incubation time.

Please visit our website for additional information on AlphaLISA technology at www.revvy.com

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

6760635-R Rev 01

revvity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA

(800) 762-4000 www.revvy.com

For a complete listing of our global offices, visit www.revvy.com
Copyright ©2023, Revvity, Inc. All rights reserved.