



6042 Cornerstone Ct. West, Ste. B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet **G9a Homogeneous Assay Kit** Catalog #52051

DESCRIPTION: The *G9a Homogeneous Assay Kit* is designed to measure G9a activity for screening and profiling applications. G9a is a histone methyltransferase that exhibits methylation activity toward H3-K9. The *G9a Homogeneous Assay Kit* comes in a convenient AlphaLISA® format, with biotinylated histone H3 peptide substrate, primary antibody, methylation assay buffer, and purified G9a for 384 enzyme reactions. The key to the *G9a Homogeneous Assay Kit* is a highly specific antibody that recognizes methylated substrate. With this kit, only three simple steps on a microtiter plate are required for methyltransferase detection. First, a sample containing G9a enzyme is incubated with the biotinylated substrate for two hours. Next, acceptor beads and primary antibody are added, then donor beads, followed by reading the Alpha-counts.

COMPONENTS:

Catalog #	Components	Amount	Storage	Storage
51001	G9a enzyme	8 µg	-80°C	(Avoid freeze/thaw cycles!)
52120	100 µM S-adenosylmethionine	2 x 250 µl	-80°C	
52140E	Primary antibody 5	2 x 12.5 µl	-80°C	
	Biotinylated histone H3 peptide substrate (G9α)	10 µl	-80°C	
	4x G9a assay buffer 1A (add DTT before experiment)	3 ml	-20°C	
	4x Detection buffer 1	2 ml	-20°C	

MATERIALS REQUIRED BUT NOT SUPPLIED:

DTT (Dithiothreitol), 0.5M (Sigma, Cat. # D0632)
AlphaLISA anti-mIgG acceptor beads, 5 mg
AlphaLISA anti-mIgG acceptor beads, 5 mg/ml (PerkinElmer #AL105C)
AlphaScreen Streptavidin-conjugated donor beads, 5 mg/ml (PerkinElmer #6760002)
Optiplate -384 (PerkinElmer #6007290)
AlphaScreen microplate reader

APPLICATIONS: Great for studying enzyme kinetics and HTS applications.

CONTRAINDICATIONS: Green and blue dyes that absorb light in the AlphaScreen signal emission range (520-620 nm), such as Trypan Blue. Avoid the use of the potent singlet oxygen quenchers such as sodium azide (NaN₃) or metal ions (Fe²⁺, Fe³⁺, Cu²⁺, Zn²⁺ and Ni²⁺). The presence of >1% RPMI 1640 culture medium leads to a signal reduction due to the presence of excess biotin and iron in this medium. MEM, which lacks these components, does not affect AlphaScreen assays.

STABILITY: At least one year from date of receipt when stored as directed.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Ct. West, Ste. B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

REFERENCES: Dillon SC, *et al.* 2005. *Genome Biology* 6:227.

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

Step 1:

- 1) Dilute **Biotinylated histone H3 peptide substrate** 40-fold with water. Dilute only the amount required for the assay. Discard any unused diluted **Biotinylated histone H3 peptide substrate** after use.
- 2) Add 125 μ l 0.5M DTT (not provided) to 3-ml tube with **4x G9a Assay Buffer**. Prepare **1x G9A buffer** by adding 1 part of **4x G9A buffer** to 3 parts water (v/v).
- 3) Prepare the master mixture: N wells x (2 μ l **4x G9A buffer** + 1 μ l **S-adenosylmethionine** (100 μ M) + 1 μ l **Biotinylated substrate**). Add 4 μ l to wells designated "Positive Control", "Test Sample", and "Blank". To wells labeled "Substrate Control", add 2 μ l 4x G9A buffer + 1 μ l Biotinylated substrate plus 1 μ l water.
- 4) Add 3 μ l of Inhibitor solution of each well labeled as "Test Inhibitor". For the "Positive Control", "Substrate Control" and "Blank", add 3 μ l of the same solution without inhibitor (Inhibitor buffer).
- 5) Thaw **G9a** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Aliquot **G9a** enzyme into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C. *Note: G9a is very sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme. Pre-incubation of enzyme with the inhibitor before starting the reaction may provide better results.*
- 6) Dilute **G9a** in **1x G9A assay buffer** at 0.3-0.9 ng/ μ l. Keep diluted enzyme on ice until use. Discard any unused diluted enzyme after use.

	Positive Control	Substrate Control	Test Sample	Blank
4x G9A assay buffer	2 μ l	2 μ l	2 μ l	2 μ l
100 μ M S-adenosylmethionine	1 μ l	-	1 μ l	1 μ l
Biotinylated substrate (diluted)	1 μ l	1 μ l	1 μ l	1 μ l
H ₂ O	1 μ l	2 μ l	1 μ l	1 μ l
Test Inhibitor/Activator	-	-	3 μ l	-
Inhibitor Buffer (no inhibitor)	3 μ l	3 μ l		3 μ l
G9a (0.3-0.9 ng/ μ l)	2 μ l	2 μ l	2 μ l	-
1x G9A assay buffer	-	-	-	2 μ l
Total	10 μl	10 μl	10 μl	10 μl

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com

- 7) To the wells designated as "Blank", add 2 μ l of **1x G9A buffer**.
- 8) Initiate reaction by adding 2 μ l of diluted **G9a** enzyme to the wells designated "Positive Control", "Substrate Control", and "Test Sample". Incubate for 2 hours at 30°C.

Protect your samples from direct exposure to light for steps 2 and 3!

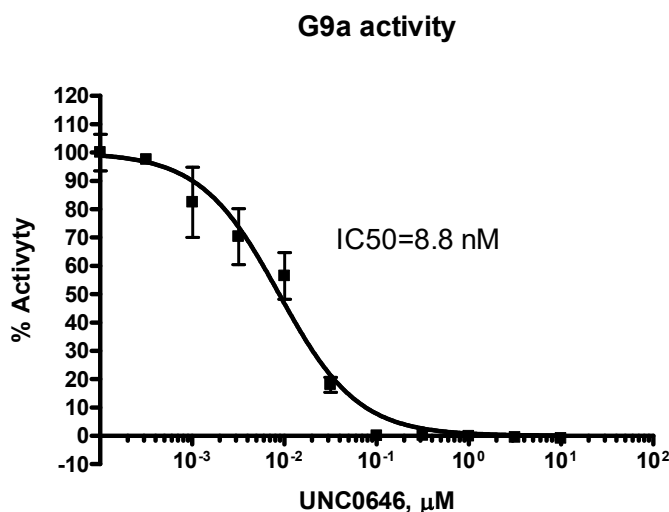
Step 2:

- 1) Dilute anti-Mouse Acceptor beads (PerkinElmer #AL105C) 1:250-fold with **1x Detection buffer 1** (made by diluting **4x Detection buffer 1** 1:4 in distilled water). Add 5 μ l per well. Shake plate briefly.
- 2) Dilute "**Primary antibody 5**" 100-fold with **1x Detection buffer 1**. Add 5 μ l per well. Shake plate. Incubate 30 min at room temperature.
(Alternatively, dilute anti-Mouse Acceptor beads (1:500) and Primary antibody 5 (1:200) with 1x Detection buffer in one step. Add 10 μ L of acceptor beads/antibody mixture per well.)

Step 3:

- 1) Dilute Streptavidin-conjugated donor beads (PE #6760002) 125-fold with **1x Detection buffer 1**. Add 10 μ l per well. Incubate for 10-15 min. at room temperature.
- 2) Read Alpha-counts. The "Blank" value is subtracted from all other values.

Example of Assay Results:



G9a inhibition by UNC0646, measured using the G9a Homogeneous Assay Kit, BPS Bioscience Cat. #52051. *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com*

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Ct. West, Ste. B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

RELATED PRODUCTS:

	Cat #:	Size:
G9a Chemiluminescent Assay Kit	52001L	96 reactions
G9a recombinant protein (insect)	51001	20 µg
G9a recombinant protein (<i>E. coli</i>)	51000	50 µg
SUV39H1 recombinant protein	51070	50 µg
SUV39H1, full length recombinant protein	51071	5 µg
SUV39H2 recombinant protein	51080	50 µg
SUV39H1 Chemiluminescent Assay Kit	52006L	96 reactions
SUV39H2 Chemiluminescent Assay Kit	52008	96 reactions
Histone H3(K9) Universal Methyltransferase Assay Kit	52072	96 reactions

Note: AlphaScreen® and AlphaLISA® are registered trademarks of PerkinElmer, Inc.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com