



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

CBX1 Inhibitor Screening Assay Kit

Catalog # 79052

DESCRIPTION: The *CBX1 Inhibitor Screening Assay Kit* is designed to measure the inhibition of CBX1 binding to its substrate. The kit comes in a convenient AlphaLISA[®] format, with enough biotinylated peptide substrate, assay buffer, detection buffer and purified GST-tagged CBX1 chromodomain to perform a total of 384 enzyme reactions. The key to the kit is the specific binding of the CBX1 chromodomain to the methylated-peptide substrate. With this kit, only three simple steps on a microtiter plate are required. First, a sample containing CBX1 and an inhibitor of choice is incubated with the biotinylated substrate for one hour. Next, acceptor beads are added, then donor beads, followed by reading the Alpha-counts.

COMPONENTS:

Catalog #	Component	Amount	Storage	
55009	CBX1, GST-tag (2-185)	1 µg	-80°C	(Avoid freeze/thaw cycles!)
	Methylated Ligand 1	400 µl	-80°C	
	Non-Methylated Ligand 1	200 µl	-80°C	
	3x CBX1 Assay Buffer	4 ml	-20°C	
	3x CBX1 Detection Buffer	3 ml	-20°C	

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

Glutathione AlphaLISA Acceptor Beads, 5 mg/ml (PerkinElmer #AL109C)
AlphaScreen Streptavidin-conjugated donor beads, 5 mg/ml (PerkinElmer #6760002S)
Optiplate-384 (PerkinElmer #6007290)
AlphaScreen microplate reader
Adjustable micropipettor and sterile tips

APPLICATIONS: Useful for the study of chromodomain binding assays, screening inhibitors, and selectivity profiling.

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 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 **1.858.829.3082**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

REFERENCE: Kaustov, L., *et al. J. Biol. Chem.* 2011; **286**(1):521-9.

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate. Use slow shaking for all incubations.

Step 1:

- 1) Prepare the master mixture: N wells × (2.5 µl **3x CBX1 Assay Buffer** + 1 µl **Methylated Ligand 1** + 1.5 µl distilled **H₂O**). Add 5 µl of master mixture to each well designated for the “Positive Control”, “Test Inhibitor”, and “Blank”. For the “Substrate Control”, add 2.5 µl **3x CBX1 Assay Buffer** + 1.5 µl distilled **H₂O** +1 µl of **Non-methylated Ligand 1**.
- 2) Prepare **1x CBX1 Assay Buffer** by diluting **3x CBX1 Assay Buffer** with water. Dilute only enough **3x CBX1 Assay Buffer** as required for the assay. Store remaining **3x CBX1 Assay Buffer** at -20°C in single-use aliquots.
- 3) Add 2.5 µl of **Inhibitor solution** to each well designated “Test Inhibitor”. For the “Positive Control”, “Substrate Control” and “Blank”, add 2.5 µl of the same **solution without inhibitor** (inhibitor buffer). *Note: Keep DMSO concentration below 1%.*

	Blank	Substrate Control	Positive Control	Test Inhibitor
3x CBX1 assay buffer	2.5 µl	2.5 µl	2.5 µl	2.5 µl
Methylated Ligand 1	1 µl	–	1 µl	1 µl
Non-Methylated Ligand 2	–	1 µl	–	–
H ₂ O	1.5 µl	1.5 µl	1.5 µl	1.5 µl
Test Inhibitor	–	–	–	2.5 µl
Inhibitor buffer (no inhibitor)	2.5 µl	2.5 µl	2.5 µl	–
1x CBX1 assay buffer	2.5 µl	–	–	–
CBX1 (0.8-1 ng/µl)	–	2.5 µl	2.5 µl	2.5 µl
Total	10 µl	10 µl	10 µl	10 µl

- 4) Thaw **CBX1** on ice. Upon first thaw, briefly spin tube containing protein to recover full content of the tube. Aliquot protein into single use aliquots. Store remaining

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

To place your order, please contact us by Phone **1.858.829.3082**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com

undiluted protein in aliquots at -80°C immediately. *Note: **CBX1** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted protein.*

- 5) Dilute **CBX1** in **1x CBX1 Assay Buffer** to 0.8-1 ng/ μl . Keep diluted protein on ice until use. Discard any unused diluted protein after use.
- 6) Add 2.5 μl of **1x CBX1 Assay Buffer** to the well designated "Blank".
- 7) Initiate reaction by adding 2.5 μl of **diluted CBX1** prepared as described above to the wells labeled "Positive Control", "Test Inhibitor", and "Substrate Control". Incubate at room temperature for 60 minutes.

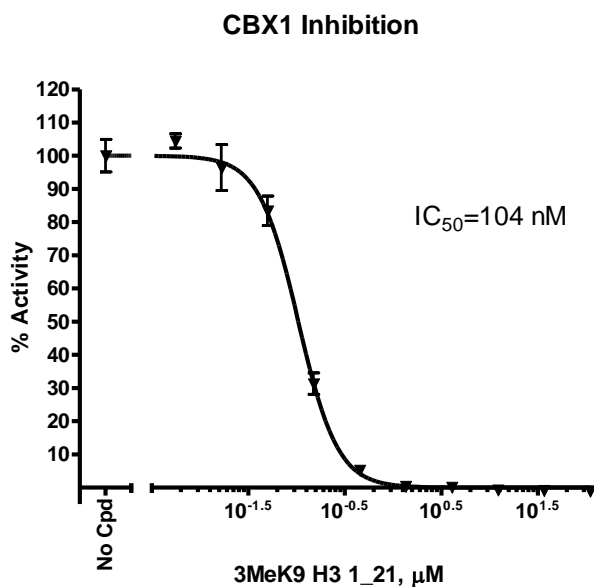
Step 2:

Note: Protect your samples from direct exposure to light!

- 1) Dilute Glutathione AlphaLISA Acceptor Beads (PerkinElmer #AL109C) 250-fold with **1x CBX1 Detection Buffer**. Add 10 μl per well. Shake plate briefly. Incubate at room temperature for 30 minutes.

Step 3:

- 1) Dilute Streptavidin-conjugated donor beads (PE #6760002S) 125-fold with **1x CBX1 Detection Buffer**. Add 10 μl per well. Incubate at room temperature for 5 - 10 minutes.
- 2) Read Alpha-counts.



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

To place your order, please contact us by Phone **1.858.829.3082**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Inhibition of CBX1 binding, measured using the CBX1 Inhibitor Screening Assay Kit, BPS Bioscience, Catalog #79052. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com.

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
CBX1, GST-tag	55009	100 µg
CBX2, GST-tag	55011	100 µg
CHD2, GST-tag	55005	100 µg
CDY1, GST-tag	55007	100 µg
L3MBTL1, GST-tag	55000	100 µg
L3MBTL1, His-tag	55002	100 µg
L3MBTL2, His-tag	55018	100 µg
L3MBTL3, GST-tag	55013	100 µg
L3MBTL3, His-tag	55014	100 µg
UHRF1 (2-793), His-Flag tag	55001	50 µg
UHRF1 (108-286), His-tag	55004	100 µg
UHRF1 (108-286), GST-tag	55003	100 µg
L3MBTL1 TR-FRET Assay Kit	55200	384 rxns.
L3MBTL1 Inhibitor Screening Assay Kit	55100	384 rxns.

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 **1.858.829.3082**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com