



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

Bromodomain Non-acetylated Ligand 4 Catalog #: 33011

DESCRIPTION: Bromodomains function as acetyl-lysine binding domains, and bromodomain/ acetyl-lysine recognition can regulate histone modification, DNA replications, chromatin remodeling, and gene transcription. The *Bromodomain Non-acetylated Ligand 4* is control peptide for the acetylated Bromodomain Ligand 4, Cat. #33010. It is derived from histone H3 and contains a C-terminal GG linker, followed by a biotinylated Lys. This peptide is suitable for use as a negative control for bromodomains including CREBBP.

APPLICATIONS: Study of bromodomain binding assays, screening inhibitors and selectivity profiling.

PURITY: ≥ 95%

SUPPLIED AS: Aqueous solution

QUANTITY: 0.5 ml @ 20 µM

STORAGE: Store in aliquots at -20°C. Avoid freeze/thaw cycles.

REFERENCE(S): Muller, S., Filippakopoulos, P., Knapp, S., *Expert Rev. Mol. Med.* 2011 Sep 13;13:e29.

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
CREBBP (1081-1197), GST-tag	31128	100 µg
CREBBP (1081-1197), His-tag	31119	100 µg
CREBBP Inhibitor Screening Assay	32607	384 reactions
Bromodomain Ligand 4	33010	0.5 ml

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