



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

## Data Sheet

### ***Bromodomain Ligand 4*** Catalog #: 33010

**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 Ligand 4* is a peptide derived from histone H3 that includes an acetylated lysine. It contains a C-terminal GG linker, followed by a biotinylated Lys. This peptide is suitable for use as a ligand for several 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 Non-acetylated Ligand 4	33011	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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)