



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

### ***BET Bromodomain Ligand*** Catalog #: 33000

**DESCRIPTION:** The *BET Bromodomain Ligand* is peptide histone H4 (1-21), acetylated at lysine 5, 8, 12 and 16. It contains a C-terminal GG linker, followed by a biotinylated Lys. Bromodomains function as acetyl-lysine binding domains, and bromodomain/acetyl-lysine recognition can regulate histone modification, DNA replications, chromatin remodeling, and gene transcription. This peptide is suitable for use as a ligand for several bromodomains including members of the BET bromodomain family: BRD2, BRD3, BRD4, and BRDT. It also binds to other bromodomains which are not included in the BET bromodomain family.

**APPLICATIONS:** Study of bromodomain binding assays, screening inhibitors and selectivity profiling. It is suitable use several BPS Bioscience BRD Assay Kits.

**PURITY:** > 95%

**MOLECULAR WEIGHT:** 2728.2

**SEQUENCE:** SGRG-K(Ac)-GG-K(Ac)-GLG-K(Ac)-GGA-K(Ac)-RHRKVGG-K(Biotin)

**SUPPLIED AS:** Aqueous solution

**QUANTITY:** 0.5 ml @ 20 µM

**STORAGE:** -20°C

#### **REFERENCES:**

1. Jasencakova, Z. *et al. Plant Cell* **12**, 2087 (2000).
2. Johnson, C. *et al. Nucl Acids Res* **26**, 994 (1998).

#### **RELATED PRODUCTS:**

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
ATAD2A (981 – 1108), GST-tag*	31121	100 µg
ATAD2B (953 – 1080), GST-tag*	31125	100 µg
BRD2 (339 – 459), GST-tag*	31023	100 µg
BRD3 (29 – 145), GST-tag*	31032	100 µg
BRD3 (306 – 417), GST-tag*	31033	100 µg
BRD4 (49 – 170), GST-tag*	31040	100 µg
BRD4 (342 – 460), GST-tag*	31041	100 µg
TAF1 (1400 – 1651), GST-tag*	31124	100 µg

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)



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)

**RELATED PRODUCTS continued:**

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
TAF1L (1398 – 1649), GST-tag*	31107	100 µg
TAF1L (1517 – 1649), GST-tag*	31106	100 µg
ATAD2A Inhibitor Screening Kit	32601	384 rxns
ATAD2B Inhibitor Screening Kit	32605	384 rxns
BRD2 (BD2) Inhibitor Screening Kit	32522	384 rxns
BRD3 (BD1) Inhibitor Screening Kit	32513	384 rxns
BRD3 (BD2) Inhibitor Screening Kit	32523	384 rxns
BRD4 (BD1) Inhibitor Screening Kit	32514	384 rxns
BRD4 (BD2) Inhibitor Screening Kit	32524	384 rxns
TAF1 (BD1+BD2) Screening Kit	32604	384 rxns
TAF1L (BD1+BD2) Screening Kit	32603	384 rxns
TAF1L (BD2) Inhibitor Screening Kit	32602	384 rxns

\*Also available with His-tag

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)