



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 1 Catalog #: 33005

DESCRIPTION: The *Bromodomain Non-acetylated Ligand 1* is peptide histone H4 (1-21). It contains an 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 negative control for several bromodomains including members of the BET bromodomain family: BRD2, BRD3, BRD4, and BRDT.

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

PURITY: ≥ 95%

MOLECULAR WEIGHT: 2560

SEQUENCE: SGRGKGGKGLGKGGAKRHRKVGG-K(Biotin)

SUPPLIED AS: Aqueous solution

QUANTITY: 0.5 ml @ 20 µM

STORAGE: -20°C. Do not freeze thaw.

REFERENCE(S): Park, P., *et al.*, *Am J Physiol Gastrointest Liver Physiol.* **289**, G1124 (2005).

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
BRD2 (339 – 459), His-tag	31020	100 µg
BRD3 (29 – 145), His-tag	31030	100 µg
BRD3 (306 – 417), His-tag	31031	100 µg
BRD4 (49 – 170), GST-tag	31040	100 µg
BRD4 (342 – 460), GST-tag	31041	100 µg
BRD4 (49 – 170), His-tag	31042	100 µg
BRD4 (342 – 460), His-tag	31043	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

Please visit our website at: www.bpsbioscience.com