

Data Sheet

GCN5 (727 - 837)

Human, recombinant, N-terminal HIS-tag

Catalog #: 31114

Lot #: 130411 Conc.: 2.13 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, <0.01 % Tween-
20, 20% glycerol.

Stability: >6 months at -80 °C

References:

1. Li, S., *et al.*, *J Biol Chem.* 2009 Apr 3;
284(14): 9411-7.
2. Santillan, D.A., *et al.*, *Cancer Res.* 2006
Oct 15; **66(20)**: 10032-9.

Description:

Human GCN5 or KAT2A containing bromodomain, amino-acids 727 – 837 (end), GenBank Accession No. NM_021078 with N-terminal HIS-tag, MW = 14.1 kDa, expressed in an *E. coli* expression system.

Application:

Useful for the study of bromodomain binding assays, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

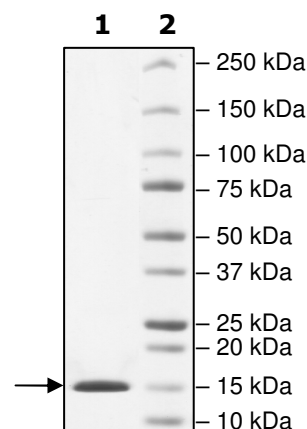
2 µg GCN5

Lane 2:

Protein Marker

MW: 14.1 kDa

Purity: ≥99%



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



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

Protein Sequence:

MHHHHHHLKDPDQLYTTLKNLLAQIKSHPSAWPFMEPVKKSEAPDYEEVIRFPIDLKTM
TERLRSRYVTRKLFVADLQRVIANCREYNPPDSEYCRCASALEKFFYFKLKEGGLIDK

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