

## Data Sheet

### ZMYND11

Human, recombinant, N-terminal GST-tag

**Catalog #:** 31144

Lot #: 130820-G2      Conc.: 0.16 mg/ml

**Formulated in:** 40 mM Tris-HCl pH 8, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, and 20% glycerol.

**Stability:** >6 months at  $-70^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

**References:**

1. Yang, H., *et al.*, *Ann Hematol.* 2010 Oct;**89(10)**:959-64.
2. Wei, G., *et al.*, *Anticancer Res.* 2003 May-Jun;**23(3A)**:2713-8.

**Description:**

Human bromodomain zinc finger, MYND-type containing 11 (ZMYND11), also known as BRAM1 and BS69, GenBank Accession No. NM\_006624, a.a. 150 - 270, with N-terminal GST-tag, MW = 41 kDa, expressed in an *E. coli* expression system.

**Application:**

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

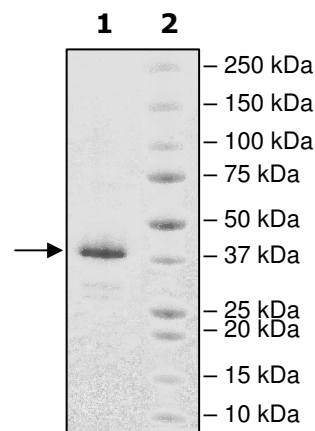
## Quality Assurance

**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
2  $\mu\text{g}$  ZMYND11

**Lane 2:**  
Protein Marker

**MW:** 41 kDa  
**Purity:**  $\geq 80\%$



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)

**Sequence:**

MSPILGYWKIKGLVQPTRLLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYID  
GDVKLTQSMARIIRYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRAYSKDFETLKVDF  
LSKLPPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLVCFKK  
RIEAIQIDKYLKSSKYIAWPLQGQWQATFGGGDHPPKSDLEVLFGGPLGSKKNTNKQEM  
GTYLRFIVSRMKERAIDLNKKGKDNKHPMYRRLVHSAVDVPTIQEKVNEGKYRSYEEFK  
ADAQLLLHNTVIFYGADSEQADIARMLYKDTCHELDELQLCKNCFYLSNARPD

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)