

## Data Sheet

### ZMYND8

Human, recombinant, N-terminal GST-tag

**Catalog #:** 31143

**Lot #:** 130828-G      **Conc.:** 0.28 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°C. *Avoid freeze/thaw cycles.*

#### References:

1. Zeng, W., *et al. Biochem Biophys Res Commun.* 2010 Apr 16;**394(4)**:1024-9.
2. Panagopoulos, I., *et al. PLoS One.* 2013 May 7;**8(5)**:e63663.

#### Description:

Human bromodomain zinc finger, MYND-type containing 8 (ZMYND8), also known as CTCL Tumor Antigen Se14-3, PRKCBP1, and RACK7, GenBank Accession No. NM\_183047, a.a. 160 - 280, 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.

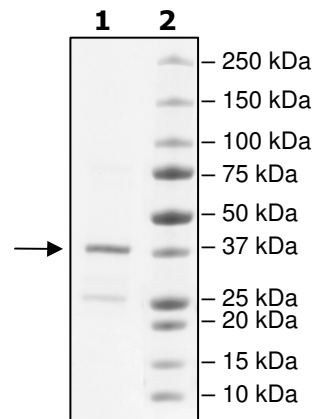
## Quality Assurance

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

**Lane 1:**  
2 µg ZMYND8

**Lane 2:**  
Protein Marker

**MW:** 41 kDa  
**Purity:** ≥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:**

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI  
DGDVKLTQSMARIYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRAYSKDFETLK  
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV  
CFKKRIEAIQIDKYLKSSKYIAWPLQGWQATFGGGDHPPKSDLEVLFGGPLGSETQSK  
AMTMLTIEQLSYLLKFAIQKMKQPGTDAFQKVPLEQHPDYAEYIFHPMDLCTLEKNA  
KKKMYGCTEAFLADAKWILHNCIIYNGGNHKLQIAKVVIKICEHEMNEIEVCPECYL

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)