

## Data Sheet

### TRIM24 (896-1014)

Human, recombinant, N-terminal His-tag

**Catalog #:** 31116

Lot #: 130403      Conc.: 3.7 mg/ml

**Formulated in:** 50 mM Tris-HCl, pH 8.0,  
138 mM NaCl, 2.7 mM KCl, and 10%  
glycerol.

**Stability:** >6 months at -80 °C

#### References:

1. Tsai, W.W., *et al.*, *Nature*. 2012 Dec 16;  
**468(7326):** 927-32.
2. Kikuchi, M., *et al.*, *Biochim Biophys Acta*.  
2009 Dec; **1793(12):** 1828-36.

#### Description:

Human bromodomain tripartite motif containing 24 (TRIM24), also known as TIF1, amino-acids 896 - 1014, (GenBank Accession No. NM\_015905) with N-terminal HIS-tag, MW = 15 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:

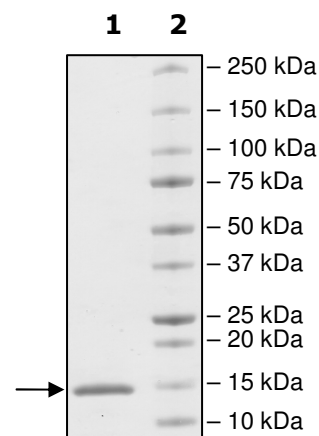
2 µg TRIM24

Lane 2:

Protein Marker

**MW:** 15 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](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)

**Protein Sequence:**

**MHHHHHGLVKLTPIDKRKRCERLLLFLYCHEMSLAFQDPVPLTVPDYKIIKNPMDLS  
TIKKRLQEDYSMYSKPEDFVADFRLIFQNCAEFNEPDSEVANAGIKLENYFEELLKNLY  
PEKRFPKPE**

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)