

Data Sheet

TRIM24 (896-1014)

Human, recombinant, N-terminal HIS-tag

Catalog #: 31116

Lot #: 120413-G Conc.: 1.7 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:

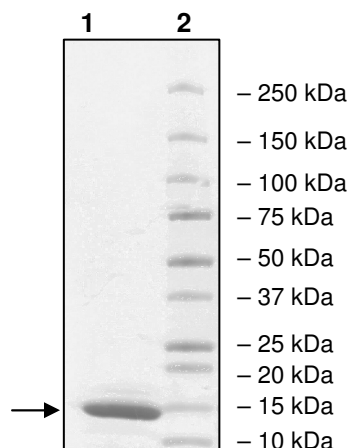
3 µg TRIM24

Lane 2:

Protein Marker
Biorad (#161-0373)

MW: 15 kDa

Purity: ≥93%



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:

MHHHHHGLVKLTPIDKRKRCERLLLFLYCHEMSLAFQDPVPLTVPDYYKIIKNPMDLS
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
Please visit our website at: www.bpsbioscience.com