

Data Sheet

TRIM24 (896-1014)

Human, recombinant, N-terminal GST-tag
Catalog #: 31127
 Lot #: 130207-D Conc.: 1.2 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, and 20%
 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.

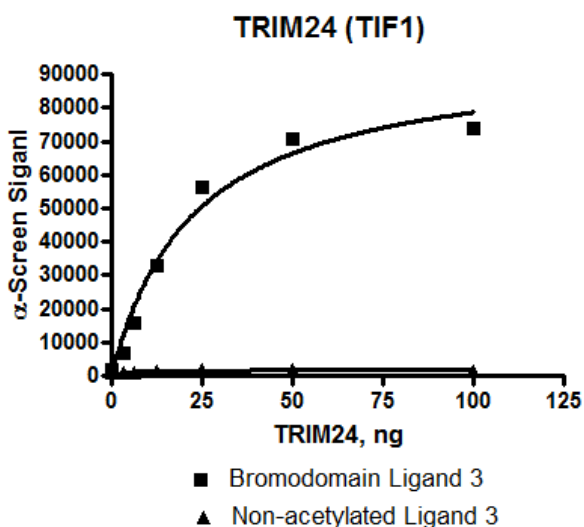
Description:

Human bromodomain tripartite motif containing 24 (TRIM24), also known as TIF1, amino-acids 896 - 1014, (GenBank Accession No. NM_015905) with N-terminal 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



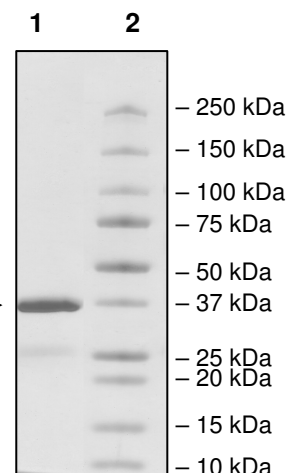
4-20% SDS-PAGE Coomassie staining

Lane 1:
3.6 µg TRIM24

Lane 2:
Protein Marker

MW: 41 kDa

Purity: ≥90%



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:

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMAIRYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRAYSKDFETLK
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIQIDKYLKSSKYIAWPLQGQWQATFGGGDHPPKSD**LEVLFGGPLGSGLVKL**
TPIDKRKCERLLLFLYCHEMSLAFQDPVPLTVPDYKIIKNPMDLSTIKKRLQEDYSMY
SKPEDFVADFRLLIFQNCAEFNEPDSEVANAGIKLENYFEELLKNLYPEKRFPKPE

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