

Data Sheet

TRIM28 (685-815)

Human, recombinant, N-terminal GST-tag

Catalog #: 31139

Lot #: 130828-G1 Conc.: 0.23 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. Herquel, B., *et al.*, *Transcription*. 2011 Sep-Oct;**2(5)**:231-236.
2. Zeng, L., *et al.*, *Nat. Struct. Biol.* 2008;**15(6)**:626-633.

Description:

Human bromodomain tripartite motif containing 28 (TRIM28), also known as KRIP-1, KAP-1, TIF1-Beta, and RNF96, GenBank Accession No. NM_005762, a.a. 685 - 815, 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.

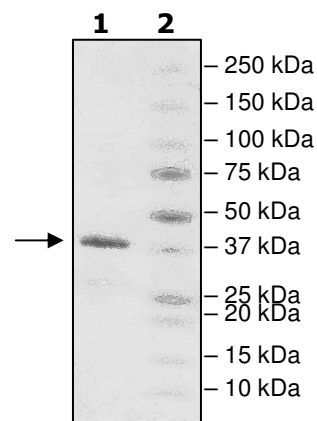
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg TRIM28

Lane 2:
Protein Marker

MW: 41 kDa
Purity: $\geq 90\%$



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

Sequence:

MSPILGYWKIKGLVQPTRLLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMARIYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRAYSDFETLK
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIPQIDKYLKSSKYIAWPLQGQATFGGGDHPPKSD**LEVLFQGPLGSDGAD**
STGVVAKLSPANQRKCERVLLALFCHEPCRPLHQLATDSTFSLDQPGGTLDLTLIRAR
LQEKLSPPYSSPQEFAQDVGRMFKQFNKLTEDKADVQSIIGLQRFFETRMNEAFGDT
KFSAVLVEPPPM

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