

Data Sheet

TRIM28 (685-815)

Human, recombinant, N-terminal His-tag

Catalog #: 31146

Lot #: 130717-G3 **Conc.:** 0.06 mg/ml

Formulated in: 40 mM Tris-HCl pH 8, 110 mM NaCl, 2.2 mM KCl, 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 His-tag, MW = 15.6 kDa, expressed in an *E. coli* expression system.

Application:

Useful for the study of bromodomain binding assays, screening inhibitors, and selectivity profiling.

Quality Assurance

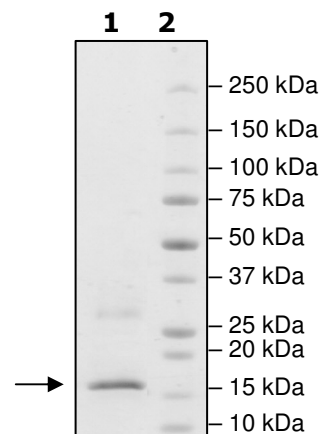
4-20% SDS-PAGE Coomassie staining

Lane 1:
1.8 μg TRIM28

Lane 2:
Protein Marker

MW: 15.6 kDa

Purity: $\geq 87\%$



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:

**MHHHHHHDGADSTGVVAKLSPANQRKCERVLLALFCHEPCRPLHQLATDSTFSLDQ
PGGTLDLTLIRARLQEKLSPPYSSPQEFAQDVGRMFKQFNKLTEDKADVQSIIGLQRF
FETRMNEAFGDTKFSAVLVEPPPM**

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