

## Data Sheet

### FALZ/BPTF

Human, recombinant, N-terminal His-tag  
**Catalog #:** 31131  
Lot #: 120412      Conc.: 1.58 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0,  
124 mM NaCl, 2.4 mM KCl, 225 mM  
imidazole, 3 mM DTT, and 10% glycerol.

**Stability:** >6 months at -80°C

**References:**

1. Buganim, Y., *et al.*, *Structure. PLoS One*. 2010;11;5(3):e9657.
2. Ruthenburg, A.J., *et al.*, *Cell*. 2011;145(5):692-706.

**Description:**

Human FALZ, also known as BPTF, GenBank Accession No. NM\_182641, a.a. 2791-2911, with N-terminal His-tag, MW=15.2 kDa, expressed in an *E. coli* expression system.

**Application:**

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

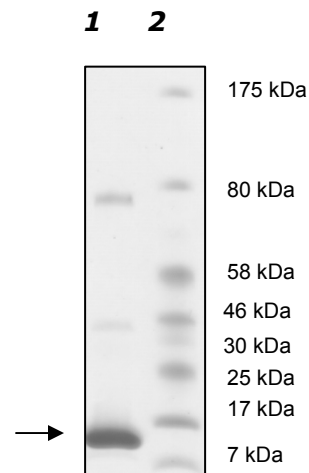
### Quality Assurance

**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
8 µg FALZ/BPTF  
**Lane 2:**  
Protein Marker

**MW:** 15.2 kDa

**Purity:** ≥88%



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:**

MHHHHHHSTEDAMTVLTPLTEKDYEGLKRVLRSLQAHKMAWPFLEPVDPNDA PDYYG  
VIKEPMDLATMEERVQRRYYEKLTEFVADMTKIFDNCRYYNPSDSPFYQCAEVLESFFV  
QKLKGFKASRSH

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)