

Data Sheet

FALZ (BPTF)

Human, recombinant, N-terminal GST-tag
Catalog #: 31134
Lot #: 130718-D **Conc.:** 4.1 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. 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 Fetal Alzheimer Antigen (FALZ), also known as Fetal Alz-50 Clone 1 Protein or bromodomain PHD finger transcription factor (BPTF), GenBank Accession No. NM_182641, a.a. 2791-2911, with N-terminal GST-tag, MW=41 kDa, expressed in an *E. coli* expression system.

Application:

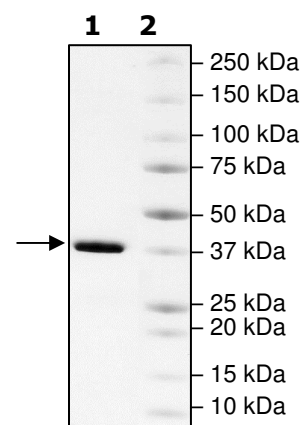
Useful for the study of bromodomain binding assays, screening inhibitors and selectivity profiling.

Quality Assurance

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

Lane 1:
3 µg FALZ
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

Sequence:

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMARIYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRAYSKDFETLK
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIQIDKYLKSSKYIAWPLQGQATFGGGDHPPKSDLEVLFGGPLGSSTEDA
MTVLTPLTEKDYEGLKRVLRSLQAHKMAWPFLEPVDPNDA PDYYGVIKEPMDLATME
ERVQRRYYEKLTEFVADMTKIFDNCRYYNPSDSPFYQCAEVLESFFVQKLKGFKASR
SH

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