

Data Sheet

BAZ1B (1335-1450)

Human, recombinant, N-terminal GST-tag
Catalog #: 31137
Lot #: 130731-G1 Conc.: 0.61 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
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. Pascual, J., *et al.*, *J Mol Biol.* 2000
Dec 15;**304(5)**:723-9.
2. Lu, X., *et al.*, *Genomics.* 1998 Dec
1;**54(2)**:241-9.

Description:

Human Bromodomain adjacent to zinc
finger domain 1B (BAZ1B), also known
as WBSR9, WBSR10, HWALp, and
WSTF, GenBank Accession No.
NM_032408, a.a. 1335-1450, with N-
terminal GST-tag, MW = 40.5 kDa,
expressed in an *E. coli* expression
system.

Application:

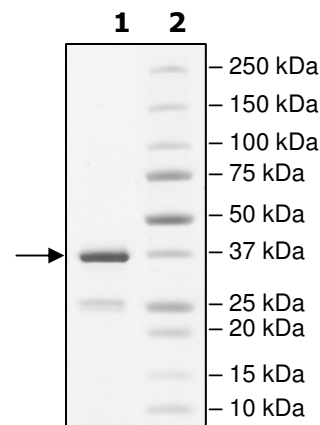
Useful for the study of bromodomain
binding assays, screening inhibitors, and
selectivity profiling.

Quality Assurance

4-20% SDS-PAGE
Coomassie staining

Lane 1:
2 μg BAZ1B
Lane 2:
Protein Marker

MW: 40.5 kDa
Purity: $\geq 80\%$



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
DGDVKLTQSMARIYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRAYSDFETLK
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIQIDKYLKSSKYIAWPLQGWQATFGGGDHPPKSDLEVLFGGPLGSKRSSR
RQSLELQKCEEILHKIVKYRFSWPFREPVTREDAEDYYDVITHPMDFQTVQNKCSGGS
YRSVQEFLTDMKQVFTNAEVYNCRGSHVLSCMVKTEQCLVALLHKHLPGHPYV

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