

Data Sheet

PRDM12

Human, recombinant, N-terminal GST-tag
Catalog #: 50320
Lot #: 130924-G1 Conc.: 0.09 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 -80°C. Avoid freeze/thaw cycles.

References:

1. Reid, A.G., Nacheva E.P., *Leukemia*. 2004 Jan;**18(1)**:178-80.
2. Xiao, B., *et al.*, *Curr Opin Struct Biol*. 2013 Dec;**13(6)**:699-705.

Description:

Human PR domain containing 12 (PRDM12), also known as PFM9 and PR Domain Zinc Finger Protein 12 (GenBank Accession No. NM_021619), a.a. 2-243, with N-terminal GST-tag, MW=54 kDa, expressed in an *E. coli* cell expression system.

Application:

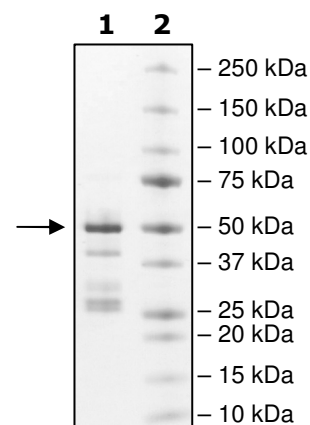
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2 µg PRDM12
Lane 2:
Protein Marker

MW: 54 kDa
Purity: ≥50%





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
DGDVKLTQSMAIRYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRIAYSKDFETLK
VDFLSKLPPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIPOIDKYLKSSKYIAWPLQGQATFGGGDHPPKSDLEVLFQ^GPLGSMGSV
LPAEALVLKTGLKAPGLALAEVITSDILHSFLYGRWRNVLGEQLFEDKSHHASPKTAF
TAEVLAQSFSGEVQKLSSLVLPAEVIAQSSIPGEGLGIFSKTWIKAGTEMGPFTGRVIA
PEHVDICKNNNLMWEVFNEDGTVRYFIDASQEDHRSWMTYIKCARNEQEQNLEVVQI
GTSIFYKAIEMIPPDQELLVWYGNSHNTFLGIPGVPGLEEDQKKNKHEDFHPADSAAG
PAGRM

[LEVLFQ^GP](#): PreScission Protease cleavage site

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