

## Data Sheet

### PRDM14

Human, recombinant, N-terminal GST-tag  
Catalog #: 50321  
Lot #: 130925-G1      Conc.: 0.13 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. Nishikawa, N., *et al.*, *Cancer Res.* 2007 Oct 15; **67(20)**:9649-57.
2. Tsuneyoshi, N., *et al.*, *Biochem Biophys Res Commun.* 2008 Mar 21; **367(4)**:899-905.

**Description:**

Human PR domain containing 14 (PRDM14), also known as PFM11 and PR Domain Zinc Finger Protein 14 (GenBank Accession No. NM\_024504), a.a. 2-412, with N-terminal GST-tag, MW=72 kDa, expressed in an *E. coli* cell expression system.

**Application:**

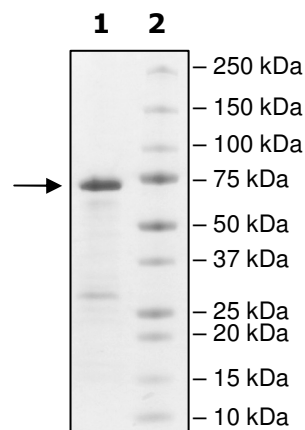
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### **Quality Assurance**

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

**Lane 1:**  
2 µg PRDM14  
**Lane 2:**  
Protein Marker

**MW:** 72 kDa  
**Purity:** ≥70%





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)

**Sequence:**

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI  
DGDVKLTQSMAIRYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRIAYSKDFETLK  
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV  
CFKKRIEAIQIDKYLKSSKYIAWPLQGQWQATFGGGDH**PPKSDLEVL**FQ^G**PLGS**ALPR  
PSEAVPQDKVCYPPESSQNLAAYYTPFPSYGHYRNSLATVEEDFQPFRQLEAAASA  
APAMPPFFFRMAPLLSPGLGLQREPLYDLPWYSKLPWYPIPHVPREVPPFLSSSH  
EYAGASSEDLGHQIIGGDNESGPCCGPDTLIPPPADASLLPEGLRRTSQLLPCSPSKQ  
SEDGPKPSNQEGKSPARFQFTEEDLHFVLYGVTPSLEHPASLHHAISGLLVPPDSSGS  
DSLQTLDKDSLQLPEGLCLMQTVFGEVPHFGVFCSSFIKGVRFQPFQKVVNASE  
VKTYGDNMWEIFEDGHLSHFIDGKGGTGNWMSYVNCARFPKEQNLVAVQCQGHI  
FYESCKEIHQNQELLVWYGDCYEKFLDIPVSLQVTEPGKQPSGPSEESAEGYRCERC  
GKVFTYK

[LEVL\*\*FQ^G\*\*P](#): 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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)