

Data Sheet

PRDM9, GST-Tag

Human, Recombinant, N-terminal GST-Tag

Catalog #: 100078

Lot #: 180522-2 **Conc.:** 0.55 mg/ml

Formulated in: 40 mM Tris-HCl pH8.0, 110 mM NaCl, 2.2 mM KCl, 20mM glutathione, 3mM DTT and 20% glycerol

Stability: At least 6 months at -80°C. *Avoid freeze/thaw cycles.*

References:

1. Baudat, F., *et al.*, *Science*. 2010 Feb; **327(5967)**: 836-40.
2. Parvanov, E., *et al.*, *Science*. 2010 Feb; **317(5967)**: 835.

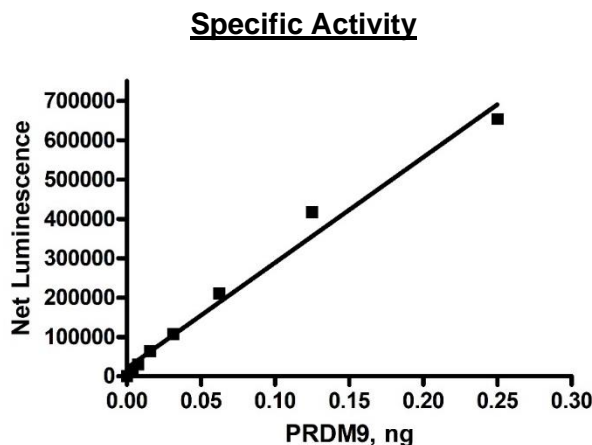
Description: Human PRDM9, also known as PR domain zinc finger protein 9, MEISETZ, and PR/SET domain 9, GenBank Accession No. NM_020227, a.a. 2-400, with an N-terminal GST-tag, expressed in an *E. coli* expression system. MW = 73 kDa.

Specific Activity: 4.86 pmol/min/μg

Assay Conditions: For standard, H3, 1-21 K4 (2Me)-Biotin was coated onto a neutravidin plate at varying concentrations. For test lots, H3,1-21-Biotin was coated onto a neutravidin plate at 0.4 μg/ml. Reaction mixture contained 50 mM Tris buffer pH 8.8, 5 mM MgCl₂, 4 mM DTT, 0.05% Tween-20, 40 μM S-adenosylmethionine, and various amounts of PRDM9—incubated for 1 hour at room temperature. Followed by 1 hour incubation with primary antibody against H3 K4 (2Me), and 30 minute incubation with secondary antibody. Luminescence was developed with HRP substrate A and B (Pierce).

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

Quality Assurance



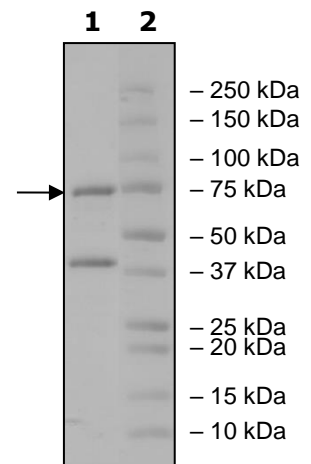
**4-20% SDS-PAGE
Coomassie staining**

Lane 1: PRDM9

Lane 2: Protein Marker

MW: 73 kDa

Purity: 46%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com