



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

PRMT4 peptide substrate, biotinylated

Catalog # 50352

DESCRIPTION: Histone peptide substrate, biotin conjugate.

SUPPLIED AS: Lyophilized Powder

QUANTITY: 200 μ l @ 400 μ M

STORAGE: -20°C

PURITY: >90% by HPLC

APPLICATIONS: Study enzyme kinetics, and screen small molecular inhibitors of PRMT4 histone methyltransferase for drug discovery and HTS applications.

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