



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

200 μ M H3K4me1 Substrate **Catalog #: 50101**

DESCRIPTION: Histone H3 peptide methylated at Lysine 4, substrate for LSD1

PURITY: >90% by HPLC

MOLECULAR WEIGHT: 2283

APPLICATIONS: Study enzyme kinetics, and screen small molecular inhibitors of LSD1 for drug discovery and HTS applications

SUPPLIED AS: water solution

QUANTITY: 500 μ l

STORAGE: -20°C or -70°C

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