



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

Data Sheet

H3K4me2 Substrate

Catalog #: 50158

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

Purity: >90% by HPLC

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

Supplied as: Water solution

Concentration: 200 μ M

Size: 500 μ l

Storage: -80°C

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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com