



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

Biotinylated Histone H3 Peptide Substrate

Catalog #: 79656

Description: Biotinylated histone H3 peptide substrate optimized for use in the PHF8 (KDM7B) Homogeneous Assay Kit (BPS Bioscience Catalog #50515).

Purity: >95%

Applications: Study enzyme kinetics, and screen small molecular inhibitors of PHF8 for drug discovery and HTS applications

Supplied as: Powder

Concentration: 5 μ M after reconstitution in 500 μ l of water

Size: 500 reactions

Storage: -20°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