



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

4x Detection Buffer

Catalog #: 79389

DESCRIPTION: The 4x Detection Buffer is optimized for use with the PRMT5 Homogeneous Assay Kit (BPS Bioscience #52052).

APPLICATIONS: Studying enzyme kinetics, and screening small molecular inhibitors of S-adenosyl homocysteine hydrolase and histone methyltransferases for drug discovery and HTS applications.

SUPPLIED AS: Tris-buffered solution

QUANTITY: 10 ml

STORAGE: -20°C. Avoid freeze/thaw cycles.

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
PRMT5 Homogeneous Assay Kit	52052	384 Reactions

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