



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

4x HDM Assay Buffer 2 **Catalog #52407**

DESCRIPTION: The *4x HDM Assay Buffer 2* is optimized for use with the following BPS Bioscience histone demethylase assay kits:

- JARID1A (KDM5A) Homogeneous Assay Kit, BPS Cat. #50510
- JARID1B (KDM5B) Homogeneous Assay Kit, BPS Cat. #50512
- JMJD1A (KDM3A) Homogeneous Assay Kit, BPS Cat. #50412
- JMJD2B (KDM4B) Homogeneous Assay Kit, BPS Cat. #50414

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

SUPPLIED AS: HEPES-based buffer containing reducing agent

QUANTITY: 10 ml

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

STABILITY: Stable for at least 1 year when stored as directed.

INSTRUCTIONS FOR USE: Please see appropriate assay kit protocol for detailed instructions.

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
4x HDM Incomplete Assay Buffer	52410	5 ml
JARID1A (KDM5A), FLAG-tag	50110	20 µg
JARID1B (KDM5B), His-FLAG-tag	50121	20 µg
JMJD1A (KDM3A), FLAG-tag	50130	20 µg
JMJD1B (KDM3B), FLAG-tag	50421	20 µg
JMJD2B (KDM4B), GST-tag	50111	100 µg
JARID1A (KDM5A) Homogeneous Assay Kit	50510	384 reactions
JARID1B (KDM5B) Homogeneous Assay Kit	50512	384 reactions
JMJD1A (KDM3A) Homogeneous Assay Kit	50412	384 reactions
JMJD2B (KDM4B) Homogeneous Assay Kit	50414	384 reactions

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

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

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

Please visit our website at: www.bpsbioscience.com