



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

Data Sheet

Blocking Buffer 1

Catalog #: 79556

Description: This Blocking Buffer 1 is optimized for minimizing background in multiple BPS Bioscience Epigenetics Assay Kits.

The BPS Bioscience Blocking Buffer improves S/B ratio of HMTs NSD2 and SetD2 ELISAs with respect to competitor blocking buffer by both decreasing background (0% Activity control, no enzyme) and increasing assay signal (100% Activity control, enzyme without inhibitor).

Applications: Suitable for use with enzymatic assays to study enzyme kinetics or screen for small molecular inhibitors of methyltransferases, demethylases, or TETs.

Supplied As: Tris-buffered solution

Format: Aqueous buffer solution

Quantity: 50 ml

Storage/Stability: Store at 4°C. Avoid freeze/thaw cycles. At least 12 months from date of receipt, when stored as directed.

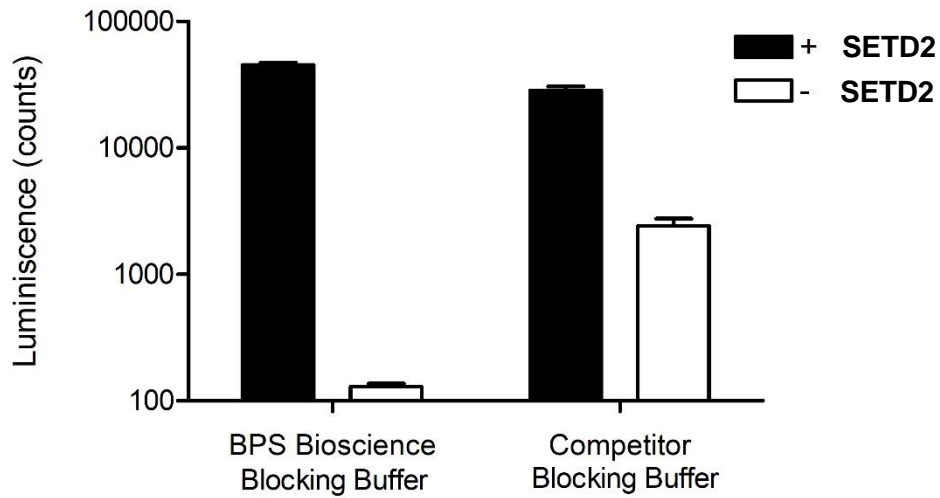
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

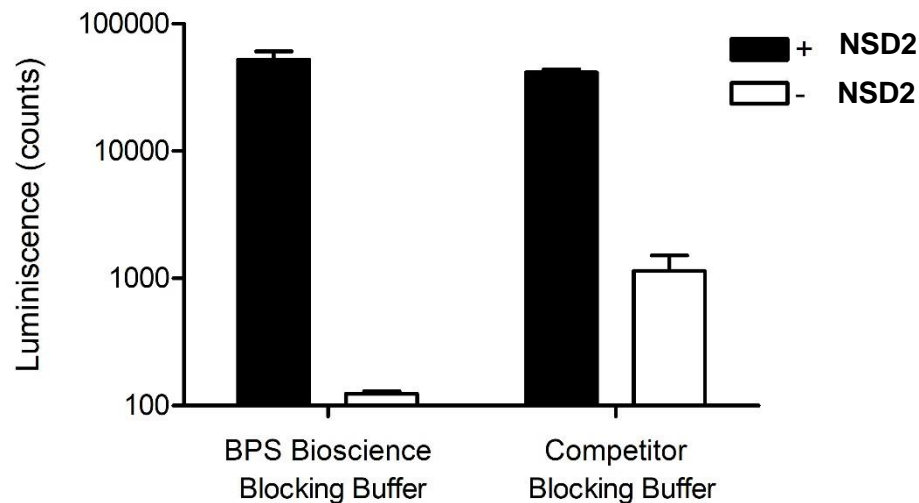
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

SetD2 Assay



NSD2 Assay



Product	Cat. #	Size
SETD2 Chemiluminescent Assay Kit	52060	96 rxns
NSD2 Chemiluminescent Assay Kit	53009	96 rxns
FTO Chemiluminescent Assay Kit	79344	96 rxns
PRMT5 Chemiluminescent Assay Kit	52002L	96 rxns

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