



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 4

Catalog #: 52100

Description: This Blocking Buffer 4 is optimized for minimizing background in various chemiluminescent methyltransferase, demethylase, and TET assays.

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**](mailto:info@bpsbioscience.com)

Please visit our website at: [**www.bpsbioscience.com**](http://www.bpsbioscience.com)