



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Data Sheet

Blocking Buffer 3

Catalog #: 79743

Description: This Blocking Buffer 3 is optimized for minimizing background in multiple PARP BPS Bioscience Assay Kits.

Applications: Suitable for use with enzymatic assays to study enzyme kinetics or screen for small molecular inhibitors in select BPS Bioscience PARP assay kits.

Supplied As: Phosphate-Buffered solution

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

Optimized for use with:

<u>Product Name</u>	<u>Cat #</u>	<u>Size</u>
PARP1 Chemiluminescent Assay Kit	80551	96 rxns
PARP14 Chemiluminescent Assay Kit	80568	32 rxns
PARP6 Chemiluminescent Assay Kit	80556	32 rxns
TNKS2 Histone Ribosylation Assay Kit (Antibody Detection)	80576	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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com