



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

10x PARP Assay Buffer Catalog #80602

DESCRIPTION: The *10x PARP Assay Buffer* is optimized for use with BPS Bioscience's poly(ADP) ribose polymerase (PARP) assay kits:

- PARP1 Chemiluminescent Assay Kit #80551 96 reactions
- PARP2 Chemiluminescent Assay Kit #80552 96 reactions
- PARP3 Chemiluminescent Assay Kit #80553 96 reactions
- PARP6 Chemiluminescent Assay Kit #80556 96 reactions
- PARP7 Chemiluminescent Assay Kit #80567 96 reactions
- PARP10 Chemiluminescent Assay Kit #80560 96 reactions
- PARP11 Chemiluminescent Assay Kit #80561 96 reactions

APPLICATIONS: Study enzyme kinetics, and screen small molecular inhibitors of poly(ADP) ribose polymerase for HTS applications in drug discovery.

SUPPLIED AS: Sodium phosphate-buffered solution

QUANTITY: 1 ml

STORAGE: -20 °C

RELATED PRODUCTS:

5x Histone Mixture	#52029	1 ml
Activated DNA	#80605	500 µl
10 x PARP Assay Mixture	#80601	300 µl
PARP1 Enzyme	#80501	10 µg
PARP2 Enzyme	#80502	10 µg
PARP3 Enzyme	#80503	10 µg
PARP6 Enzyme	#80506	10 µg
PARP7 Enzyme	#80507	10 µg
PARP10 Enzyme	#80510	10 µg
PARP11 Enzyme	#80511	10 µg

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