



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

PARP Substrate Mixture 2

Catalog #: 78371

SIZE: 250 μ l

DESCRIPTION: Aqueous buffer containing biotinylated substrate optimized for use with BPS Bioscience's poly(ADP) ribose polymerase (PARP) assay kits:

- | | | |
|-------------------------------------|--------|--------------|
| • PARP6 Chemiluminescent Assay Kit | #80556 | 96 reactions |
| • PARP7 Chemiluminescent Assay Kit | #79729 | 96 reactions |
| • PARP11 Chemiluminescent Assay Kit | #80561 | 96 reactions |
| • PARP14 Chemiluminescent Assay Kit | #80568 | 96 reactions |

APPLICATIONS: Suitable for use with cell extracts and purified PARP enzymes.

SUPPLIED AS: Tris-buffered solution

PURITY: \geq 98% purity by HPLC.

STORAGE: Store at -80°C. Avoid freeze/thaw cycles.

STABILITY: Stable for at least 6 months when stored as recommended.

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