



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 1

Catalog #: 78366

SIZE: 250 µl

DESCRIPTION: Aqueous buffer containing biotinylated substrate 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
- PARP10 Chemiluminescent Assay Kit #80560 96 reactions

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

SUPPLIED AS: Tris-buffered solution

PURITY: ≥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