



6405 Mira Mesa Blvd Ste 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

## Data Sheet

### ***Opti-PARP Assay Mixture***

**Catalog #: 79814**

**SIZE:** 300 µl

**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 | 384 reactions |
| • PARP11 Chemiluminescent Assay Kit | #80561 | 96 reactions  |
| • PARP14 Chemiluminescent Assay Kit | #80568 | 32 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.

#### **RELATED PRODUCTS:**

<u>Name</u>	<u>Catalog #</u>	<u>Amount</u>
5x Histone Mixture	52029	1 ml
Activated DNA	80605	500 µl
10 x PARP Assay Buffer	80602	1 ml
PARP6 Enzyme	80506	10 µg
PARP7 Enzyme	80527	10 µg
PARP11 Enzyme	80511	10 µg
PARP14 Enzyme	80514	10 µg

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

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