



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## **Data Sheet**

### ***CRBN Assay Buffer***

**Catalog #: 79972**

**Description:** The *CRBN Assay Buffer* is optimized for use with the Cereblon Binding Assay Kit (BPS Bioscience #79899).

**Applications:** Studying enzyme kinetics and screening small molecular inhibitors for drug discovery and HTS applications.

**Supplied As:** Liquid

**Format:** Aqueous buffer solution

**Storage/Stability:** Store at -20°C. Avoid freeze/thaw cycles. Stable at least 6 months from date of receipt, when stored as directed.

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: [\*\*info@bpsbioscience.com\*\*](mailto:info@bpsbioscience.com)

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