



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

4X GLS Assay Buffer

Catalog #: 79948

DESCRIPTION: The 4X GLS Assay Buffer is optimized for use with GLS1 Inhibitor Screening Assay Kit (BPS Bioscience #79596), or GLS2 Inhibitor Screening Assay Kit (BPS Bioscience #79925).

APPLICATIONS: Ideal for detection of Kidney-type glutaminase activity, and screening of small molecule inhibitors for drug discovery and HTS applications.

SUPPLIED AS: Aqueous solution

STORAGE: Store in aliquots at -20°C. Avoid freeze/thaw cycles. Stable at least 12 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
Please visit our website at: www.bpsbioscience.com