



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

Western ECL Substrate

Catalog #: 79572

Description: This substrate is ideal for quick, sensitive chemiluminescent detection of horseradish peroxidase (HRP). It's particularly useful for detecting HRP-labeled secondary antibodies or streptavidin-HRP in immunoblotting applications.

Applications: Western blots

Storage: Room temperature

Western Blot Assay Protocol:

1. After final wash of your membrane, prepare the fresh ECL substrate solution by mixing the equal volume of Solution A and Solution B. *Note: It is highly recommended to prepare the substrate immediately before use. Prepare only the amount required for the assay.*
2. Carefully apply the substrate solution onto the membrane.
3. Drain excess substrate solution and expose the membrane to the imaging system.

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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

Quality Assurance:

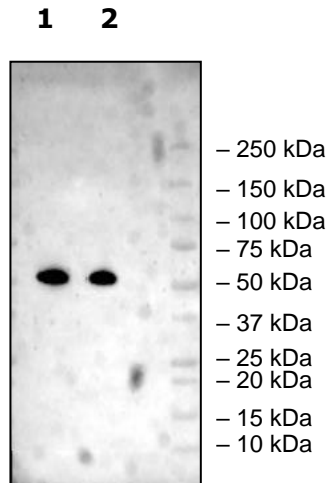


Figure 1: Acetylated Tubulin from Tissue

7.5 µg/well of total proteins from mouse spleen tissue samples were analyzed using an anti-acetyl tubulin antibody and an HRP conjugated secondary antibody (exposure time: 1 minute)

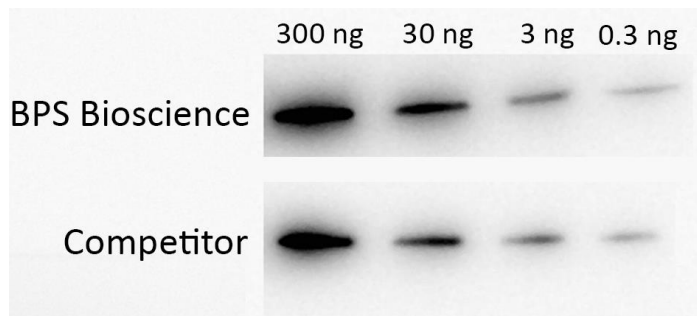


Figure 2: Purified recombinant mouse GST-PTP1B was analyzed by an anti-GST antibody and an HRP-conjugated secondary antibody (exposure time: 1 minute)

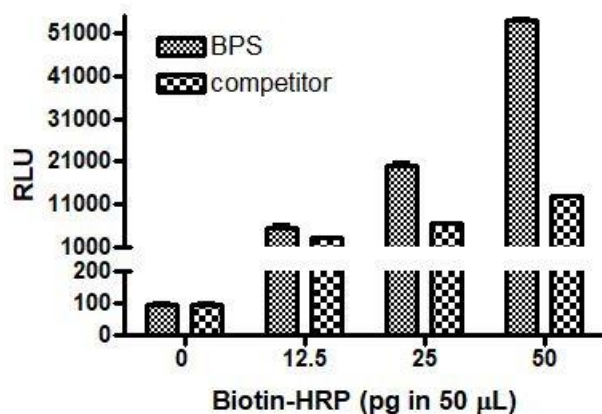


Figure 3: Acetylated Tubulin from Tissue

Biotin-HRP (Pierce) was diluted in PBS and coated onto the BPS Neutravidin plate for 30 min at room temperature. After washing 5 times with PBST, 100 µl of ECL substrate mixtures were added and read luminescence using a microplate reader.