



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

ELISA ECL Substrate

Catalog #: 79670

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: ELISA

Storage: Room temperature

ELISA Assay Protocol:

1. Just before use, mix on ice 50 μ l ELISA ECL Substrate A and 50 μ l ELISA ECL Substrate B per well. Note: It is highly recommended to prepare the substrate immediately before use. Prepare only the amount required for the assay.
2. Add 100 μ l of the substrate solution mixture into each well.
3. Immediately read sample in a luminometer or microtiter-plate capable of reading chemiluminescence. "Blank" value is subtracted from all readings.

Reading Chemiluminescence:

Chemiluminescence is the emission of light (luminescence) which results from a chemical reaction. The detection of chemiluminescence requires no wavelength selection because the method used is emission photometry and is not emission spectrophotometry. To properly read chemiluminescence, make sure the plate reader is set for LUMINESCENCE mode. Typical integration time is 1 second, delay after plate movement is 100 msec. Do not use a filter when measuring light emission. Typical settings for the Synergy 2 BioTek plate reader are: use the "hole" position on the filter wheel; Optics position: Top; Read type: endpoint. Sensitivity may be adjusted based on the luminescence of a control assay without enzyme (typically we set this value as 100).

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

Quality Assurance:

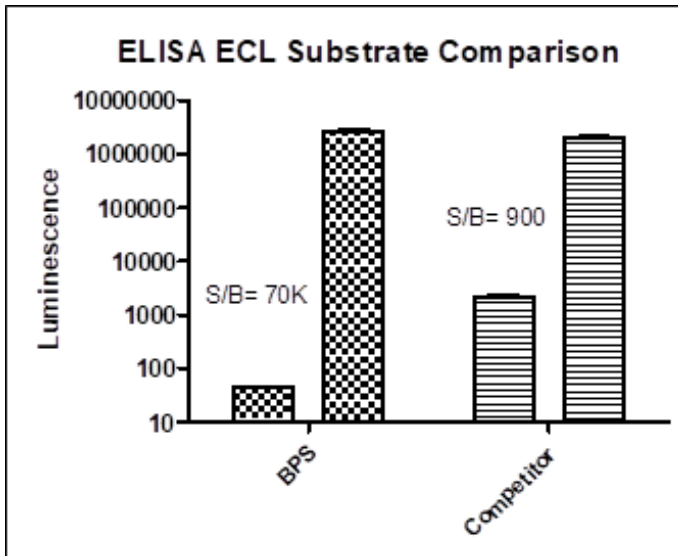


Figure 1: ELISA ECL Substrate Comparison

Signal/background comparison with BPS Bioscience's ELISA ECL Substrate versus the leading competitor.

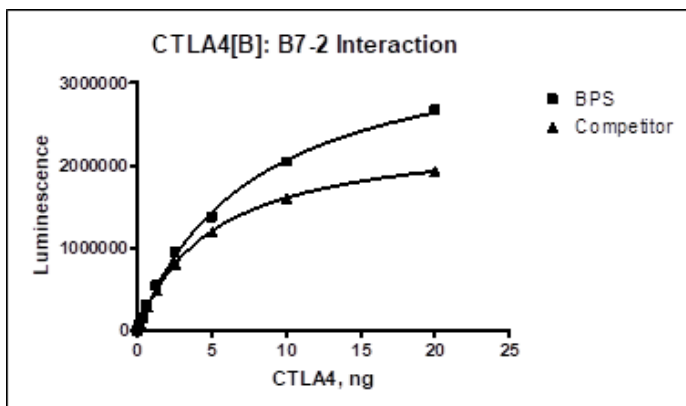


Figure 2: CTLA4 [Biotin] : B7-2 Interaction

Luminescence comparison with BPS Bioscience's ELISA ECL Substrate versus the leading competitor through using the CTLA4 [Biotinylated]: B7-2 Inhibitor Screening Assay Kit (BPS #72024).