



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 **HRP Colorimetric Substrate** Catalog #: 79651

Description: TMB (3,3', 5, 5' – tetramethylbenzidine) Substrate develops a deep blue color in the presence of horseradish peroxidase (HRP)- labeled conjugates measurable at 650 nm and turns yellow when stopped by acidification and read at 450 nm.

Supplied as: Liquid

Applications: The HRP Colorimetric substrate is used in a variety of BPS Assay Kits related to coronavirus, immunotherapy, PARP, and cytokine related research.

Storage/Stability: At least 6 months at +4°C. TMB substrate is light-sensitive and should be protected from direct sunlight or UV sources.

Optimized for Use With:

<u>Product</u>	<u>Catalog #</u>	<u>Size</u>
Spike S1 (SARS-CoV-2): ACE2 Inhibitor Screening Colorimetric Assay Kit	79994	96 reactions
SARS-CoV-2 IgG Detection Kit (Colorimetric Trimer Anti-Spike IgG detection)	79975	96 reactions
SARS-CoV-2 Spike Trimer (S1+S2):ACE2 Inhibitor Screening Colorimetric Assay Kit	79999	96 reactions
SARS-CoV-1 Spike Trimer (S1+S2):ACE2 Inhibitor Screening Colorimetric Assay Kit	78012	96 reactions

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