

TREVIGEN[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures.

Anti-Glyceraldehyde-3-Phosphate Dehydrogenase Rabbit Polyclonal Antibody

Catalog #: 2275-PC-1

Size: 100 μ l

Description: The abundance of glyceraldehyde-3-phosphate dehydrogenase (G3PDH) in eukaryotic cells is relatively unaffected by external factors. In Western blot analysis the level of G3PDH, a ~38 kDa protein, can be used as a reference value for comparisons between different cell lysates or gel loadings.

Physical State: Purified IgG is provided in phosphate buffered saline without preservative.

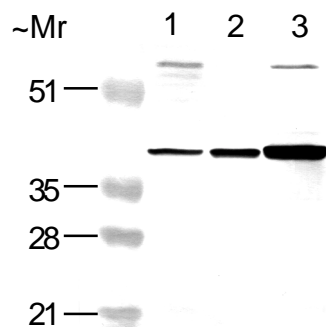
Immunogen: A synthetic peptide corresponding to a portion of the human G3PDH sequence.

Ig Class: G3PDH specific rabbit IgG.

Specificity: The antibody detects human and mouse G3PDH (other species not tested).

Storage: The antibody can be stored frozen at -20°C to -80°C for at least one year. Avoid repeated freezing and thawing by aliquoting smaller portions of the antibody into microtubes and storing at -20°C or -80°C.

Applications: For western blotting and immunoprecipitation, an antibody dilution of between 1:1000 and 1:5000 is recommended. Empirical testing may be required.



Cell lysates from L292 cells (lane 1), WEHI 7.1 cells (lane 2), and INT 407 cells (lane 3) were analyzed by SDS-PAGE and Western blot using a 1:1000 dilution of G3PDH in 5% nonfat milk in PBS, 0.05% Tween[®]20. The blot was developed using Trevigen's Blue Membrane Solution.

Cell Lysates for Western Blotting:

To prepare total cell lysates, cells are solubilized in 1X SDS gel sample buffer (20 mM dithiothreitol, 6% SDS, 250 mM Tris (pH 6.8), 10% glycerol, and 0.05% bromophenol blue) at 5×10^5 - 1×10^6 cells per ml. The extracts are heated in a boiling water bath for 5 minutes prior to electrophoresis on 12% Tris-Glycine SDS-PAGE gel.

Procedure for Immunoblotting using Peroxidase Detection:

Blocking solution: 5% (w/v) nonfat dry milk in PBS.

Antibody diluent: 5% (w/v) nonfat dry milk, 0.05% Tween[®] 20 in PBS.

Transfer the electrophoresed proteins to nitrocellulose membrane by Western transfer. Incubate the membrane for 1/2 hour at room temperature in blocking solution.

Incubate the membrane for 1 hour at room temperature (or overnight at 4°C) in 1:1000 to 1:5000 dilution of antibody in antibody diluent. Empirical determination of primary antibody concentration will be required for optimal results.

Wash the membrane at room temperature for 15 minutes with 3 changes of PBS, 0.05% Tween 20.

Incubate the membrane at room temperature for 1 hour in antibody diluent containing a dilution of anti-rabbit HRP. Empirical determination of secondary antibody concentration will be required for optimal results.

Wash the membrane for 15 minutes with 3 changes of 0.05% Tween 20 in PBS, then rinse in water.

Develop peroxidase reaction Trevigen's Blue Membrane Solution (Cat# 4857-20-13).

Tween 20 is a registered trademark of ICI Americas, Inc., Wilmington, DE

Anti-human G3PDH Polyclonal Antibody

Catalog #: 2275-PC-1

Storage: -20°C to -80°C

TREVIGEN[®]

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-216-2801 • e-mail: info@trevigen.com • www.trevigen.com

TREVIGEN[™]

1-800-873-8443