

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Anti-Glyceraldehyde-3-Phosphate Dehydrogenase Rabbit Polyclonal Antibody

Catalog #: 2275-PC-020

Volume: 20 µl

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 (figure 1).

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 Conditions: This antibody can be stored at -20°C or -80°C. Avoid repeated freeze-thawing by aliquoting into smaller portions.

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

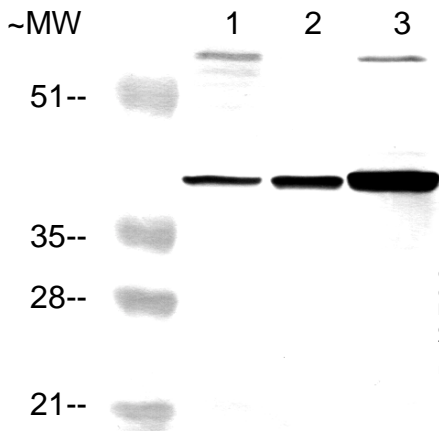


Figure 1. Cell lines vary regarding their G3PDH content. Cell lysates from L292 cells (lane 1), WEHI 7.1 cells (lane 2), and INT 407 cells (lane 3) were analyzed by SDS-PAGE and Western blot using a 1:1000 dilution of G3PDH in 5% nonfat milk in PBS, 0.05% Tween®20. The blot was developed using HRP membrane 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 with HRP membrane solution.

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

DNA Damage Antibodies:

Catalog #	Description	Size
4411-PC-100	γ -H2AX polyclonal	100 μ l
4410-PC-100	Fen-1 polyclonal	100 μ l
4350-MC-100	UVssDNA mAb (clone C3B6)	100 μ g
4354-MC-50	Anti-8-oxo-dG mAb (clone 2E2)	50 μ l
4335-MC-100	Anti-PAR polymer mAb (clone 10HA)	100 μ l
4336-BPC-100	Anti- PAR polymer polyclonal	100 μ l
4338-MC-50	Anti-human/murine-PARP mAb (clone C2-10)	50 μ g

Anti-human G3PDH Polyclonal Antibody

Catalog#: 2275-PC-020

Storage: -20 to -80°C

TREVIGEN®
1-800-873-8443