

Product Information

Anti-DHFR (N-ter)

Developed in Rabbit, Affinity Isolated Antibody

Product Number: **D1067**

Product Description

Anti-DHFR (N-ter) is developed in rabbit using a synthetic peptide corresponding to amino acids 27-40 of mouse DHFR, conjugated to KLH via a C-terminal added cysteine residue as immunogen. The immunizing sequence is conserved in mouse, human and rat. The antibody is affinity purified on the immunizing peptide immobilized on agarose.

Anti-DHFR (N-ter) antibody reacts specifically with DHFR. Staining of the DHFR band by immunoblotting is inhibited by the immunizing peptide.

DHFR (Dihydrofolate reductase) is a 187 amino acids enzyme that catalyzes the reduction of folic acid in two NADPH-dependent steps, which first yield 7,8-dihydrofolate (DHF) and then 5,6,7,8-tetrahydrofolate (THF).^{1,2} THF is the coenzyme for thymidilate synthetase in the biosynthesis of thymidine, participating in the biosynthesis of amino acids and purines as well. Since folic acid reduction is essential to generate active coenzyme, DHFR becomes critical for DNA replication.^{2,3} For this reason, much research has been done to determine the structure and mechanism of action of the enzyme.^{3,4} Inhibitors of DHFR, such as methotrexate, are folate analogs which can bind to the active site and deactivate the enzyme. Methotrexate has important implications for cancer therapy.⁵ From another angle, structural and enzymatic properties of DHFR led to development of a variety of screenings, in which DHFR can function in a fashion that resembles reporter genes. The screenings are based on the fact that DHFR can be dissected into two halves that can reassemble to form an active enzyme.⁶ Thus, each of the two halves of DHFR can be expressed as two fusion proteins that when interacting with each other can restore the DHFR enzymatic activity. The readout of this protein-protein interaction and consequent enzymatic activity, can be either restoration of growth in bacterial, yeast and plant, or receptor activation.⁷⁻¹¹ The approach was proved to be useful for receptor-ligand, antigen-antibody and other interactions.^{8,10,11} Antibodies specific for DHFR are a useful tool for following and identifying such protein interactions.

Reagents

The product is provided as a solution in 0.01M phosphate buffered saline pH 7.4 containing 1% bovine serum albumin and 15 mM sodium azide as a preservative.

Antibody Concentration: approx. 1.0 mg/ml

Precautions and Disclaimer

Due to the sodium azide content a material safety sheet (MSDS) for this product has been sent to the attention of the safety officer of your institution. Consult the MSDS for information regarding hazards and safe handling practices

Storage/Stability

For continuous use store at 2-8 °C for up to one month. For extended storage, freeze in working aliquots. Repeated freezing and thawing is not recommended. Storage in "frost-free" freezers is not recommended. If slight turbidity occurs upon prolonged storage, clarify the solution by centrifugation before use. Working dilution samples should be discarded if not used within 12 hours.

Procedure

Procedure for Immunoblotting

Note: The whole procedure is performed at room temperature.

1. Separate extracts containing DHFR or DHFR fusion proteins from sample lysates using a standard sodium dodecyl sulfate polyacrylamide gel electrophoresis (SDS-PAGE) protocol.
Note: The amount of extract depends on the level of expression of the fusion protein and the specific application.
2. Transfer proteins from the gel to a nitrocellulose membrane.
3. Block the membrane using a solution of PBS containing 5% non-fat dry milk (PBS, Product No. D8537; non-fat dry milk, Product No. M7409) for at least 60 minutes.

4. Wash the membrane three times for 5 minutes each in PBS containing 0.05% Tween 20 (Product No. P3563).
5. Incubate the membrane with anti-DHFR antibody as the primary antibody in PBS containing 0.05% Tween 20, with agitation for 120 minutes
6. Wash the membrane three times for 5 minutes each in PBS containing 0.05% Tween 20.
7. Incubate the membrane with anti-rabbit IgG, peroxidase conjugate (Product No. A0545) as the secondary antibody at the recommended concentration in PBS, containing 0.05% Tween 20. Incubate with agitation for 60 minutes. Adjust the product concentration to maximize detection sensitivity and to minimize background.
8. Wash the membrane three times for 5 minutes each in PBS containing 0.05% Tween 20.
9. Treat the membrane with a peroxidase substrate.

Procedure for Immunoprecipitation

1. Centrifuge 40 μ l of a 1:1 suspension of protein A-agarose beads (Product No. P3476) for 1 min. at 12000g, and then wash twice with 1 ml RIPA buffer (50 mM Tris Base, 0.25% w/v Deoxycholate, 1% NP40, 150 mM NaCl, 1mM EDTA, pH 7.4) at 4°C.
2. Add anti-DHFR antibody diluted in PBS, and incubate by swinging head-over-tail for 1 hour at room temperature.
3. Centrifuge 1 min at 12,000g, wash twice with 1 ml RIPA at 4°C.
4. Add 0.1-1.0 ml of cell extract containing DHFR tagged protein to the beads, or purified DHFR (see note), and incubate from 2 hours to overnight at 4°C, while swinging head-over tail.
Note: The amount of cell extract depends on the level of expression of the tagged protein and the specific application.
5. Spin down beads; remove supernatant.
6. Wash beads four times with 1ml RIPA buffer and once with PBS by vortex and short spin.

7. Resuspend pellet in 25 μ l 2XSDS-PAGE sample buffer. Boil sample for 5 min and spin down. The sample is ready to be loaded on an SDS-PAGE gel.

Product Profile

0.5-1.0 μ g/ml of the antibody detects, by immunoblotting, about 100 ng of purified recombinant DHFR, using a chemiluminiscent substrate.

0.5-1.0 μ g of the antibody can precipitate between 100-200 ng of purified DHFR.

Note: In order to obtain best results in different techniques and preparations we recommend determining optimal working dilutions by titration test.

References

1. Oefner, C., et. al., *Eur. J. Biochem.*, 174, 77-385 (1988).
2. Chen, M. J., et. al., *J. Biol. Chem.*, 259, 3933-3943 (1984).
3. Mariani, B.D., et. al., *Proc. Natl. Acad. Sci. USA*, 78, 4985-4989 (1981).
4. Davies, J.F., et. al., *Biochemistry*, 29, 9467-9479 (1990).
5. Banerjee, D., et. al., *Biochem. Biophys. Acta*, 1587, 164-173 (2002).
6. Michnick, S.W., et. al., *Curr. Opin. Struct. Biol.*, 11, 472-477 (2001).
7. Pelletier, J. N., et. al., *Nat. Biotechnol.*, 17, 639-640 (1999).
8. Mossner, E., et. al., *J. Mol. Biol.*, 308, 115-122 (2001).
9. Subramaniam, R., *Nat. Biotechnol.*, 19, 769-722 (2001).
10. Tucker, C.L., *Nature*, 19, 1042-1046 (2001).
11. Remy, I., et. al., *Science*, 283, 990-993 (1999).

NV/JK 9/2003

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.