



Mouse anti-6xHistidine IgG conjugated to Biotin

Product Number B1-1711
Amount 100 μ g
Store at 4°C

Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at 4°C in the dark. Phycobiliproteins are sensitive to freeze-thaw cycles; after reconstitution, store at 2-8°C in the dark, do not freeze.

Handling

Prior to use, reconstitute to 1 ml with 50% glycerol, vortex gently and allow material to reconstitute for 20 minutes. Avoid exposure to heat.

Buffer

Upon rehydration with 1 ml of 50% glycerol, the product is in 25mM Tris (pH 7.2), 150mM sodium chloride, and 2mM sodium azide as a preservative.

Stability

Lyophilized material is stable for up to one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 6 months. If further dilution of the conjugate is required, use diluted material within one week.

Antibody/Antigen Information

Mouse IgG₁ clone AD1.1.10

Note

For research use only, not for diagnostic or therapeutic use.

References

Coward, P., Conn, M., Tang, J., Xiong, F., Menjares, A., & Reagan, J. D. (2009). Application of an allosteric model to describe the interactions among retinol binding protein 4, transthyretin, and small molecule retinol binding protein 4 ligands. *Anal Biochem* (Vol. 384).

Day, J. E. H., Sharp, S. Y., Rowlands, M. G., Aherne, W., Workman, P., & Moody, C. J. (2010). Targeting the Hsp90 chaperone: synthesis of novel resorcylic acid macrolactone inhibitors of Hsp90. *Chemistry* (Vol. 16).

Vassilev, L. T., Tovar, C., Chen, S., Knezevic, D., Zhao, X., Sun, H., et al. (2006). Selective small-molecule inhibitor reveals critical mitotic functions of human CDK1. *PNAS* (Vol. 103).

Zhou, V., Han, S., Brinker, A., Klock, H., Caldwell, J., & Gu, X. -ju. (2004). A time-resolved fluorescence resonance energy transfer-based HTS assay and a surface plasmon resonance-based binding assay for heat shock protein 90 inhibitors. *Anal Biochem* (Vol. 331).

