



Rabbit anti-T7 tag IgG conjugated to Biotin

Product Number B1-1832
Amount 100 µg
Store at 4°C

Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at -20° Celsius.

Handling

Prior to use, reconstitute to 1 ml with distilled water, vortex gently and allow material to reconstitute for 20 minutes. After reconstitution store the product at 4°C.

Buffer

Upon rehydration with 1 ml distilled water the product is in PBS (pH 7.4) and 50 mM sucrose.

Stability

Lyophilized material is stable for up to one year.

Western blot detection

Recommended dilution of 1:500 - 1:10000 depending on target protein concentration.

Antibody/Antigen Information

Rabbit polyclonal versus sequence MASMTGGQQMG.

Note

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

References

Itoh T, Fujita N, Kanno E, Yamamoto A, Yoshimori T, Fukuda M. Golgi-resident Small GTPase Rab33B Interacts with Atg16L and Modulates Autophagosome Formation. *Mol Biol Cell*. 2008 Jul; 19(7): 2916-2925.

Day B, Dahlbeck D, Staskawicz BJ. NDR1 Interaction with RIN4 Mediates the Differential Activation of Multiple Disease Resistance Pathways in Arabidopsis. *Plant Cell*. 2006 Oct; 18(10): 2782-2791.

Rayala SK, Martin E, Sharina IG, Molli PR, Wang X, Jacobson R, Murad F, Kumar R. Dynamic interplay between nitration and phosphorylation of tubulin cofactor B in the control of microtubule dynamics. *PNAS*. 2007 Dec 4; 104(49): 19470-19475.



COLUMBIA BIOSCIENCES