



Rabbit anti-S tag IgG conjugated to Biotin

Product Number B1-1831
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 KETAAAKFERQHMS.

Note

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

References

Plagmann I, Chalaris A, Kruglov AA, Nedospasov S, Rosenstiel P, Rose-John S, Scheller J. Transglutaminase-catalyzed covalent multimerization of Camelidae anti-human TNF single domain antibodies improves neutralizing activity. *J Biotechnol.* 2009 Jun 15;142(2):170-8.

Carson SD, Kim KS, Pirruccello SJ, Tracy S, Chapman NM. Endogenous low-level expression of the coxsackievirus and adenovirus receptor enables coxsackievirus B3 infection of RD cells. *J Gen Virol.* 2007 Nov;88(Pt 11):3031-8.

Cao L, Shen G, Zhu Y, Wang W, Zhao X, Ye Q, Zhu H, Lu Z, Si J. Characterization of a single-chain variable fragment (scFv) antibody directed against the human asialoglycoprotein receptor. *Biotechnol Appl Biochem.* 2006 May;44(Pt 2):65-72.



COLUMBIA BIOSCIENCES