

Rabbit anti-S tag IgG conjugated to Alkaline Phosphatase

Product Number	AP-1831
Lot Number	AP014-21-002
Amount	100 µg
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

Rabbit polyclonal versus sequence KETAAAKFERQHMDS.

<u>Note</u>

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.

4985 Winchester Blvd. Frederick, MD 21703 301.732.5415 www.columbiabiosciences.com For technical inquiries: info@columbiabiosciences.com For sales inquiries: sales@columbiabiosciences.com