



Rabbit anti-S tag IgG conjugated to Dylight® 650

Product Number D10-1831
Lot Number XXXX
Approximately 5-7
Amount 100 µg
Store at 2-8°C

Immunogen

KETAAAKFERQHMS

Form/Shipping & Storage

Supplied as a liquid. Upon receipt, store at 2-8°C in the dark, do not freeze. Phycobiliproteins are sensitive to freeze-thaw cycles.

Handling

We recommend that the investigator determine the appropriate working concentration for their specific application. Avoid exposure to heat and light.

Buffer

100mM NaPO₄ (pH 7.4), 100mM NaCl, 2mM Sodium Azide

Stability

Product should be stored at 2-8°C in the dark and be used within 1 year. If further dilution of the conjugate is required, use diluted material within one month.

Antigen Info

Antibody was raised against S-tag with amino acid sequence KETAAAKFERQHMS.

Note

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

Spectral Characteristics

Visible absorption maxima 652
Emission maximum 672

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