



Rabbit anti-S tag conjugated to Dylight® 550

Product Number D9-1831
Lot Number XXXX
Approximately 5-7
Amount 100 µg
Concentration 1 mg/mL
Antigen Amino acid sequence KETAAKFERQHMS
Store at 2-8°C

Form/Shipping & Storage

Upon receipt, store at 2-8 °C. Fluorescent dyes are sensitive to light. Please store in the dark and avoid exposure to sunlight.

Handling

We recommend that the investigator determine the appropriate working concentration for their specific application.

Buffer

0.1M sodium phosphate pH7.4, 0.1M NaCl, 2mM NaN₃

Stability

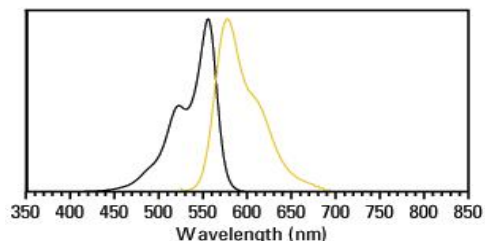
Product should be stored at 2-8°C in the dark and be used within 1 year. If further dilution of the conjugate is required, use diluted material within one month.

Note

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

Spectral Characteristic

Visible absorption maxima 562 nm
Emission maximum 576 nm



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