



Rabbit Anti-EP4 Receptor (C-term.) conjugated to R-Phycoerythrin

Product Number D5-1866
Lot# RPE102-17-003
Amount 50 µg total protein

Form/ Storage

Conjugate supplied as a lyophilized powder. Upon receipt, store at 2-8°C in the dark.

Handling

Avoid exposure to heat and light. Prior to use reconstitute to 0.5 ml with distilled deionized water, vortex and allow it to sit on ice for 20 minutes.

Buffer

Upon reconstitution, the product is in 100 mM sodium phosphate (pH 7.4), 50 mM sucrose, 100 mM sodium chloride, 0.1% BSA, and 2 mM sodium azide. The concentration of the conjugate is 100 µg/ml.

Stability

Lyophilized material is stable for one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 3 months.

Note

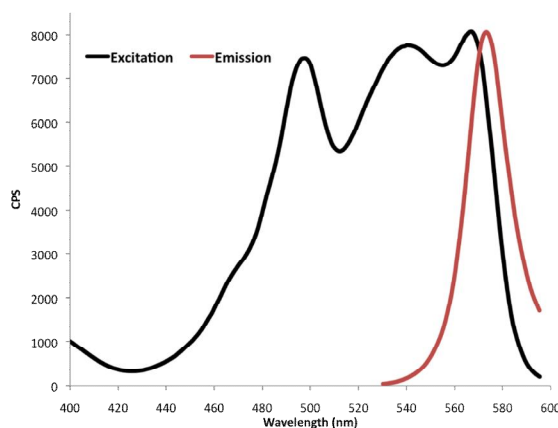
For research use only, not for diagnostic or therapeutic use.

Antigen Info

EP4 receptor C-terminal amino acids 459-488 (GSGRAGPAPKGSSLQVTFPSETLNLSEKCI)

Reactivity

Human, murine, rat, and ovine EP4 receptor; non-reactive with EP1, EP2, and EP3 receptors; other species not tested.



Fluorescence excitation and emission spectra of R-phycoerythrin in 100 mM sodium phosphate (pH 7.2) + 1 mM EDTA and 1 mM sodium azide. Emission scan was taken with excitation at 498 nm. Excitation scan was taken with emission at 575 nm.

Spectral Characteristics

Visible absorption maxima 565>540>498
Emission maximum 578 nm

Fluor:Protein -0.91:1

References

- An, S., Yang, J., Xia, M., et al. Cloning and expression of the EP2 subtype of human receptors for prostaglandin E2. *Biochem. Biophys. Res. Commun.* 197, 263-270 (1993).
- Bastien, L., Sawyer, N., Grygorczyk, R., et al. Cloning, functional expression, and characterization of the human prostaglandin E2 receptor EP2 subtype. *J. Biol. Chem.* 269, 11873-11877 (1994).
- Coleman, R.A., Eglén, R.M., Jones, R.L., et al. Classification of prostanoid receptors IUPHAR receptor compendium. *IUPHAR Compendium 1-12* (1997).
- Narumiya, S., Sugimoto, Y., and Ushikubi, F. Prostanoid receptors: structures, properties, and functions. *Physiol. Rev.* 79, 1193-1226 (1999).

4985 Winchester Blvd. Frederick, MD 21703
301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com