



Rabbit anti-15-Hydroxy PGDH IgG conjugated to R-Phycoerythrin

Product Number D5-1864
Amount 100 µg total protein
Store at 4°C

Form/ Storage

Supplied as a lyophilized powder. Upon receipt, store at 2-8°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

Avoid exposure to heat and light. Prior to use reconstitute to 1 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, 150 mM sodium chloride, 0.1% BSA as a stabilizer, and 2 mM sodium azide as a preservative.

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.

Antigen Info

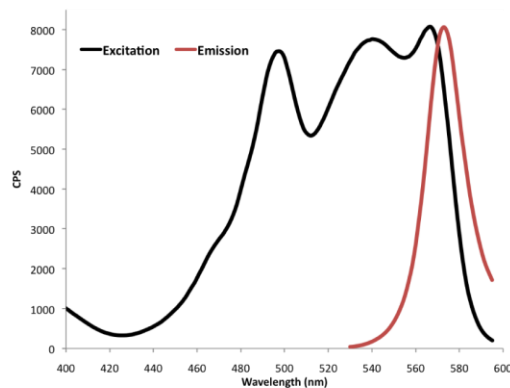
NAD⁺-dependent 15-hydroxy PGDH amino acids 92-105 (AGVNNEKNWEKTLQ)

Reactivity

Human, bovine, guinea pig, and baboon 15-hydroxy PGDH; other species not tested.

Note

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



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

Concentration

After reconstitution to 1.0 ml
0.1 mg/mL
Fluor:Protein = ~2:1

References:

Ensor, C.M., Yang, J.-Y., Okita, R.T., et al. J. Biol. Chem. 265, 14888-14891 (1990).
Duffy, D.M., Dozier, B.L., and Seachord, C.L. J. Clin. Endocrinol. Metab. (2004).
Hansen, H.S. Prostaglandins 12, 647-679 (1976).
Jörnvall, H., Persson, B., Krook, M., et al. Biochemistry 34, 6003-6013 (1995).
Krook, M., Marekov, L., and Jörnvall, H. Biochemistry 29, 738-743 (1990).
Wermuth, B. NADP-dependent Prostaglandins 44, 5-9 (1992).

